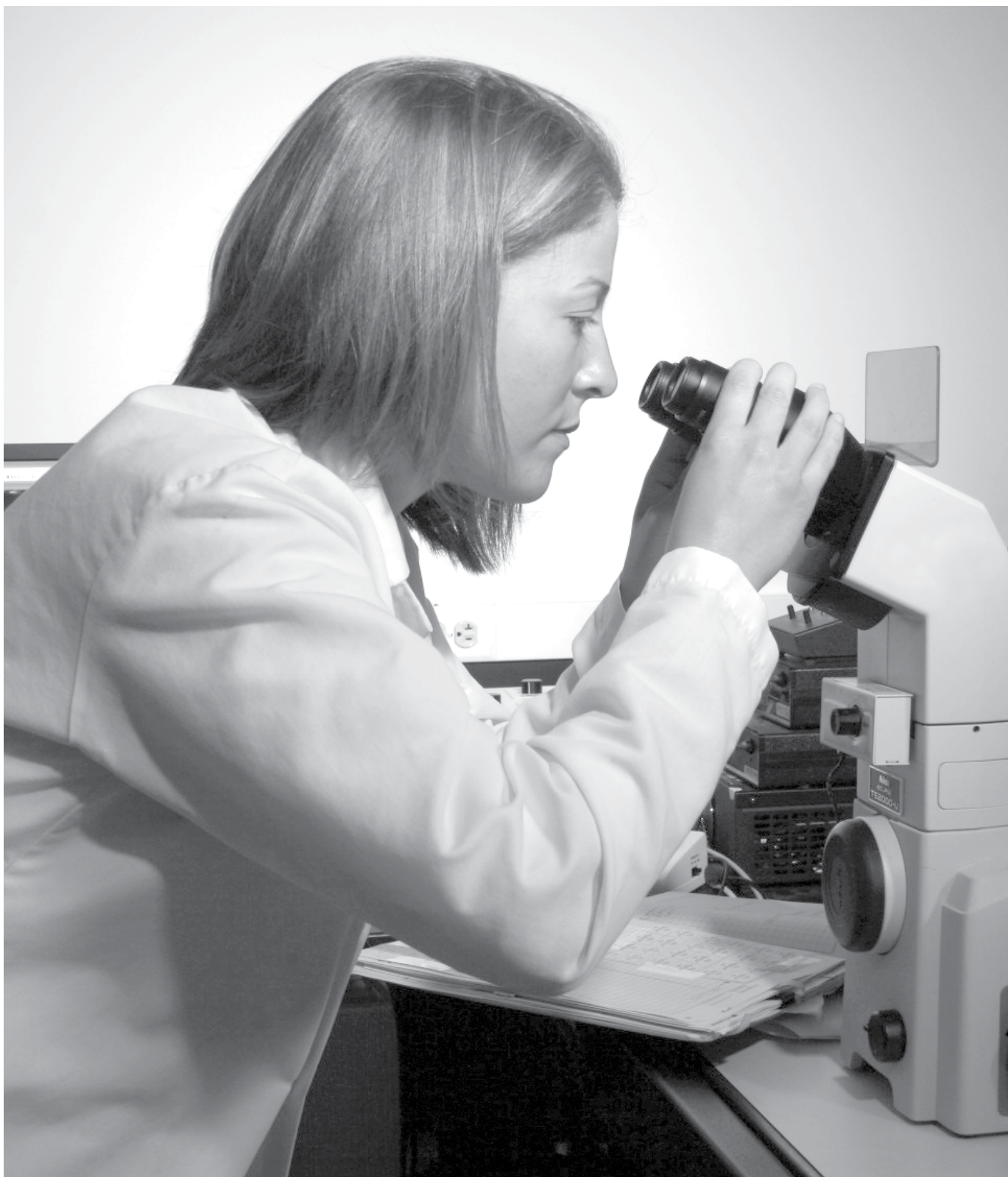




# A GUIDE TO SUCCESSFUL IMMUNOFLUORESCENCE

FROM CELL SIGNALING TECHNOLOGY | [www.cellsignal.com](http://www.cellsignal.com)

# IF



Immunofluorescence is a powerful tool for elucidating the complex signaling events that underlie biological processes and disease. This guide highlights critical steps in the immunofluorescence protocol and demonstrates how protocol changes can affect the final outcome of your experiment.



**Cell Signaling**  
TECHNOLOGY®



# Introduction

Immunofluorescence (IF) combines the use of antibodies with fluorescence imaging techniques to visualize target proteins and other biomolecules within fixed cell or tissue samples. This process can reveal the localization, relative expression, and even activation states of target proteins. The power of IF is that it provides data that is both graphical and quantifiable.

When performing IF experiments, proteins of interest can be detected using either primary antibodies covalently conjugated to fluorophores (direct detection) or a two-step approach with unlabeled primary antibody followed by fluorophore-conjugated secondary antibody (indirect detection). Either method allows the user to combine multiple fluorophores (multiplex analysis), making IF ideal for investigating protein co-localization, changes in subcellular localization, differential activation of proteins within a cell, identification of different cell subsets, and other analyses.

At Cell Signaling Technology (CST), our goal is to provide highly specific antibodies that yield strong, specific signal with minimal background. Our scientists screen a large number of antibodies and recommend only those best suited for the application. Our validation efforts include extensive protocol optimization and antibody titration to determine the best working conditions for each antibody. In addition, our scientists validate supporting reagents, such as fluorophore-conjugated secondary antibodies, to enhance antigen detection and improve the efficiency of IF protocols.

In this application guide, we will highlight the critical steps in our protocol for IF, introduce important concepts about antibody performance and design of controls, and provide supporting data to explain our recommendations.

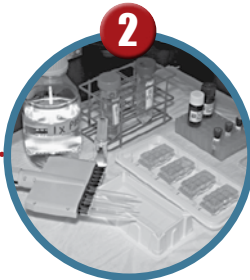
Immunofluorescence techniques referenced in this guide		
Technique	Starting material	CST protocols available
<b>IF-IC</b> (immunocytochemistry)	Cultured cell lines or primary cells	IF Standard, IF Methanol-fixed, IF Methanol-perm
<b>IF-F</b> (frozen)	Frozen tissue	IF Standard, IF Methanol-perm
<b>mIHC</b> (multiplexed immunohistochemistry)	Formalin-fixed, paraffin-embedded (FFPE) tissue	mIHC/Paraffin

# 9 Step Protocol for a Successful Immunofluorescence Experiment



## Design Your Experiment

- Experimental controls
- Selecting antibodies and fluorophores:  
direct vs. indirect detection methods
- Multiplexing



## Prepare Your Cells or Tissue

- Tissue
- Plating conditions
- Optimizing cell density and cell health



3

## Fixation – Stabilize and Preserve Sample

- Choice of fixative
- Fixation incubation time

4

## Permeabilization

- Choice of detergents or alcohols
- For paraffin-embedded tissue: deparaffinization/rehydration  
and antigen retrieval steps

5

## Blocking

- Choice of blocking reagent

6

## Immunostaining

- Primary antibody dilution
- Washing and secondary incubation



7

## Counterstain (Optional)

8

## Mount Sample



9

## Data Collection and Imaging

For more data on experimental controls visit

[www.cellsignal.com/ifvalidation](http://www.cellsignal.com/ifvalidation)

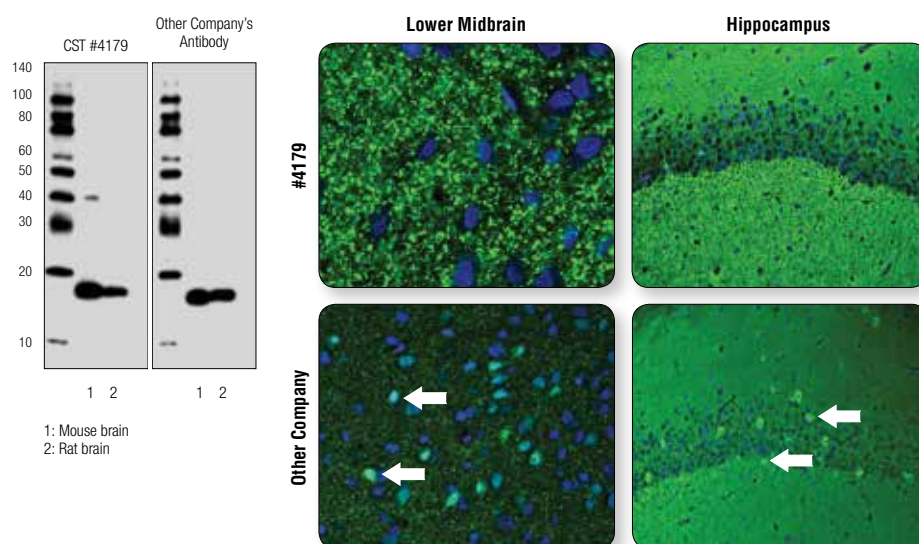
# The Importance of a Well-Validated Antibody

The primary antibody is a critical component of an IF experiment, and its performance directly affects data quality. Detection of a specific band on a western blot (WB) is not sufficient to guarantee a chosen antibody will perform in IF (see figure on this page). WB analysis subjects proteins to harsh reducing/denaturing conditions that alter protein structure, so an epitope recognized by an antibody approved for WB may be buried and/or inaccessible for IF, where proteins remain in their native state.

## Validate Specificity to Avoid Misleading IF Results

The  $\alpha$ -Synuclein protein is highly expressed in the brain and its dysfunction plays a role in neurodegenerative diseases, such as Parkinson's disease. In healthy tissue,  $\alpha$ -Synuclein is expected to be localized to presynaptic terminals where it associates with synaptic vesicles. CST scientists compared  $\alpha$ -Synuclein (D37A6) XP<sup>®</sup> Rabbit mAb #4179 side-by-side with another company's  $\alpha$ -Synuclein antibody, using each at the manufacturer's recommended dilution. Both antibodies performed as expected in WB. In IF, punctate staining consistent with presynaptic localization of  $\alpha$ -Synuclein was observed in midbrain sections for #4179, while the other company's antibody showed a less punctate pattern, but critically mis-reported localization in nuclei or soma (arrows). This experiment illustrates the importance of application-specific validation. **(A)**

**(A) A "clean" WB is not sufficient to ensure performance and reliability of an antibody in IF analysis.** WB analysis of extracts from mouse and rat brain using  $\alpha$ -Synuclein (D37A6) XP<sup>®</sup> Rabbit mAb #4179 or the other company's antibody to  $\alpha$ -Synuclein (left). Confocal IF analysis (right) of [mouse] lower midbrain and hippocampus sections using #4179 (upper row) or the other company's antibody (lower row). Neuronal soma/nuclei that are mis-labeled for  $\alpha$ -Synuclein with the other company's antibody are noted with white arrows.



## TROUBLESHOOTING TIP

**Setting up a small-scale pilot experiment is a good way to check how an antibody will perform in a protocol.** This can be used to optimize parameters and troubleshoot critical steps in the protocol before moving on to larger-scale experiments, and can save time and reagents.



# Critical Steps in the Recommended Protocol

An optimized protocol is necessary to achieve consistent, reliable IF results. We have tested common variations in fixation, permeabilization, and antibody concentration for many of our IF-approved antibodies. This section of the guide will highlight some of the data we use to support our IF protocol recommendations.

## Design Your Experiment

### Experimental Controls

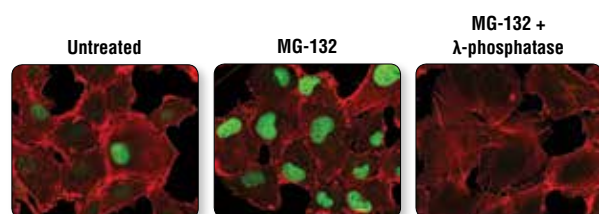
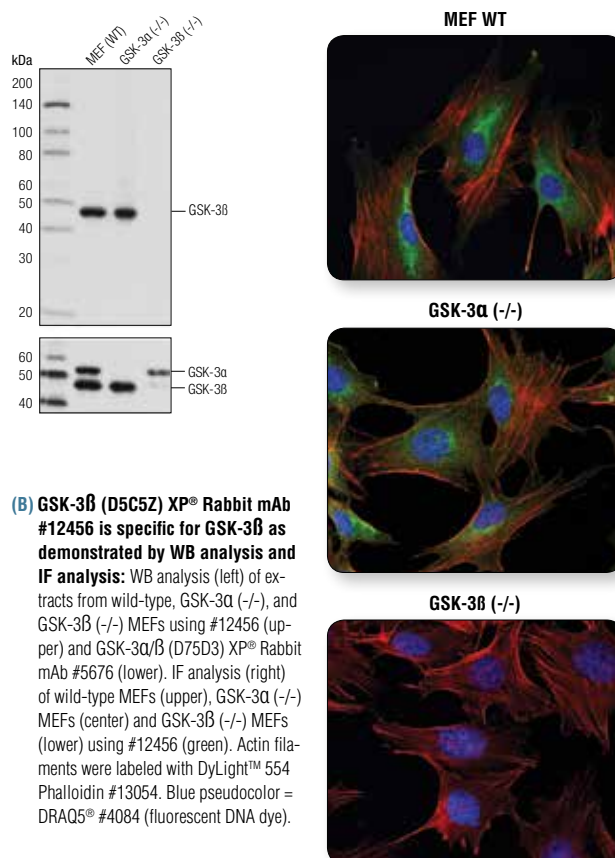
Incorporation of appropriate controls is important to confirm that the only changes between samples are in the experimental variable(s), and that the reagents – including antibodies – are performing as expected. Such controls could involve pharmacological treatments, addition of extracellular ligands to modulate signaling pathways, or comparison of cells with differential gene expression (knockout, siRNA, etc.). Typically, variables and controls are performed in parallel such that the fixation step and subsequent processing can be performed at the same time. The type of control used is dependent on the type of experiment. In the following sections we provide examples of controls our scientists use to evaluate antibody performance in IF-IC, some of which you may want to incorporate to confirm specificity in your cell type.

#### Knockout Cell Lines to Verify Target Specificity

If you are investigating a target with multiple isoforms, it may be useful to perform controls using knockout cell lines. For example, two genes encode glycogen synthase kinase (GSK-3), designated GSK-3 $\alpha$  and GSK-3 $\beta$ , which are regulated at different serine residues. When performing an IF experiment, you would want to know if the antibody you're using recognizes one isoform, both, or neither. This can be confirmed using cells that are known to be positive or negative for expression of the target of interest. Using wild-type (positive for GSK-3 $\beta$ ), GSK-3 $\alpha$  knockout (positive for GSK-3 $\beta$ ), and GSK-3 $\beta$  knockout (negative for GSK-3 $\beta$ ) mouse embryonic fibroblasts (MEFs) demonstrates that GSK-3 $\beta$  (D5C5Z) XP<sup>®</sup> Rabbit mAb #12456 specifically recognizes the correct isoform in IF. **(B)**

#### Confirming Phospho-specificity by Modulating Target Phosphorylation State

Antibodies specific for post-translational modifications (PTMs), including phosphorylation, acetylation, ubiquitination, cleavage, and others, can reveal important information about the biological function of the target protein. In some cases, changes in PTM may coincide with changes in expression and/or localization. For this reason, confirmation of PTM changes in cells can be done using PTM-specific antibodies in conjunction with enzymatic or chemical agonists and/or inhibitors to modulate the activation state of the target protein. **(C)**



**(C) Experimental perturbations confirm PTM-specificity of an antibody.** Confocal IF analysis of HT-1080 cells untreated (left) or treated with Protease MG-132 alone (center) or with MG-132 followed by  $\lambda$ -phosphatase (right), using Phospho-Cyclin D1 (Thr286) (D29B3) XP<sup>®</sup> Rabbit mAb #3300 (green). Actin filaments were labeled with DyLight<sup>™</sup> 554 Phalloidin #13054.

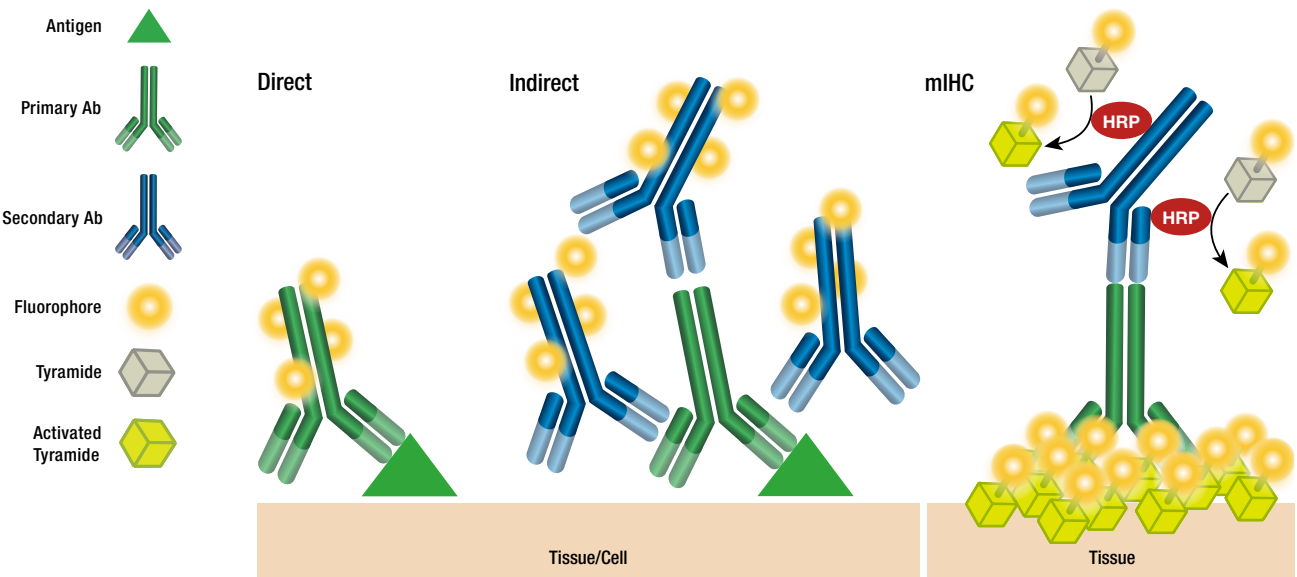
# Selecting Antibodies and Fluorophores

Detection of antigens can be performed by either applying fluorophore-conjugated primary antibodies (direct detection) or by applying primary antibodies first, followed by washing and applying secondary antibodies (indirect detection). Direct detection methods save time and allow multiplexing, using antibodies raised in the same host, while indirect detection traditionally offers higher sensitivity due to signal amplification (multiple secondary antibodies recognizing one primary). However, with the advent of more sensitive cameras used in fluorescent microscopy, the difference in signal quality has been shrinking.

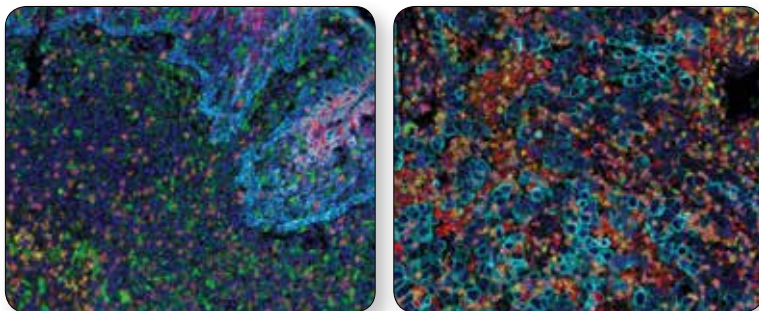
Selecting appropriate primary-secondary pairs and fluorophore conjugates that will work with your imaging system is a critical part of experimental design. For multiplexed IF-IC, you will need to consider the wavelengths of excitation sources and spectral characteristics of filter sets available to you when selecting fluorophores for your experiments. It is prudent to minimize spectral overlap in order to avoid bleed-through fluorescence from the shorter wavelength channel into the longer wavelength channel. Careful examination of the spectral characteristics of a dye's excitation and emission curve can go a long way to minimize mixed signals that include the target antibody and an antibody in a shorter wavelength channel. A good pairing is Alexa Fluor 488® with Alexa Fluor® 647. Keep in mind that if you use an organelle counterstain (see Counterstaining on pg. 10), it will mean one less channel available for multiplexing.

For tissue sections embedded in paraffin, Fluorescent Multiplex Immunohistochemistry (mIHC) is an alternative protocol to chromogenic IHC that employs tyramide signal amplification (TSA®) (serial deposition of tyramide-fluorophore conjugates) at the site of the target. **(D)** Sequential rounds of primary and HRP-conjugated secondary antibody incubation, tyramide-fluorophore deposition, and antibody stripping enables multiple antibodies raised in the same host (e.g., rabbit monoclonals) to be used together. CST antibodies that have been approved for use in IHC-P are compatible with mIHC. CST's protocol for mIHC is available at [www.cellsignal.com/CSTProtocols](http://www.cellsignal.com/CSTProtocols).

There are advantages and disadvantages to both direct and indirect immunofluorescence methods. The table below outlines some of these. The specific experiment and the reagents available will determine which one you choose.



<b>Antigen Detection</b>	Direct immunofluorescence with primary antibody conjugated to a fluorophore	Indirect immunofluorescence with secondary antibody conjugated to a fluorophore	Indirect immunofluorescence HRP-conjugated secondary antibody with tyramide amplification
<b>Multiplex Protocol</b>	Parallel staining	Parallel staining	Serial staining
<b>Primary Antibody</b>	Same host species can be used for multiple targets	Different host species or isotope for each target	Same host species can be used
<b>Secondary Antibody</b>	No	Yes	Yes
<b>Signal Amplification</b>	None	Moderate	High
<b>Incubation Steps</b>	1	2	Dependent on the number of targets
<b>Background</b>	Lower	Potentially higher for samples containing endogenous IgG's	Potentially higher
<b>Notes</b>	Dependent on availability of conjugated antibodies; limited sensitivity	Offers flexibility but still dependent on availability of validated antibodies from different species/isotypes	Requires antigen retrieval and antibody removal steps; spectral detection required for >3 targets



**(D) Examples of mIHC performed on paraffin-embedded tissue.**

Fluorescent mIHC analysis of paraffin-embedded human tonsil (left) using PD-L2 #82723 (orange), PD-L1 #13684 (red), CD8 #70306 (magenta), PD-1 #43248 (yellow), CK #4545 (cyan), CD68 #76437 (green); and paraffin-embedded human lung cancer (right) using TIM-3 #45208 (yellow), CD8 #70306 (magenta), CD68 #76437 (red), Lag-3 #15372 (orange), PD-1 #43248 (green), CK #4545 (cyan).

## Prepare Your Cells or Tissue

### Tissue

Tissue sections can be prepared either by quick freezing in optimal cutting temperature (OCT) medium (IF-F) or by embedding in paraffin (IHC-P). Frozen samples should be sectioned using a cryostat and allowed to air dry on the slide 10–15 minutes prior to fixation in the next step. Formalin-fixed, paraffin-embedded (FFPE) tissue blocks are another common preparation in the IHC-P protocol, which require deparaffination and antigen retrieval steps before incubation with antibodies.

### Cells: Plating Conditions (IF-IC)

Because the IF-IC protocol ends with imaging fixed and stained cells on a fluorescence microscope, it must begin with seeding of cells (either passaged from immortalized cell lines, or isolated primary cells) on a support material compatible with fluorescence microscopy. Typical support formats include glass-bottom cell culture dishes, glass coverslips (kept in plastic culture dishes) prepared with polylysine and/or extracellular matrix components to support adherent cell culture, and commercially available multiwell chambers mounted on glass slides that are compatible with microscopes.

### Cells: Optimizing Cell Health and Density (IF-IC)

Culture conditions may affect cell health, morphology, expression/location of target antigen, and ultimately dictate the quality of your IF data. Ensure your cells remain healthy by regularly checking the culture media for pH changes and inspecting for signs of cell stress (such as multinucleated cells) at low magnification on a microscope. Also check that the confluence of the cells is appropriate for the cell type and target. Some proteins exhibit confluence-dependent changes in localization, such as YAP. **(E)**

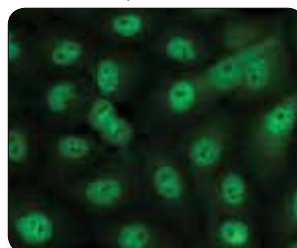
## Fixation and Permeabilization

**At CST, we test different fixation and permeabilization conditions to identify the combination that gives the optimal signal for each antibody.** The fixation and permeabilization of your samples are key steps that can determine your experiment's failure or success. The ideal fixative preserves a "life-like" snapshot while quickly stopping the degradative process of autolysis by crosslinking and inhibiting endogenous enzymes. Consult the datasheet for CST's recommended protocol, which varies by product.

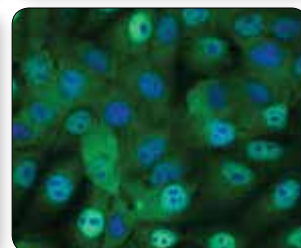
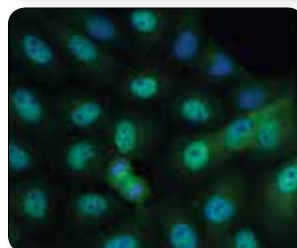
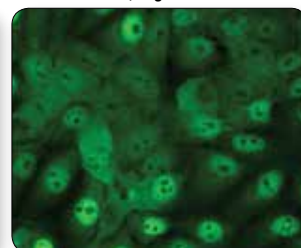
### TROUBLESHOOTING TIP

Cell lines that are grown in suspension can be ported into the IF protocol by low-speed centrifugation onto coverslips after performing the fixation step in suspension, or by performing the fixing and staining steps in suspension and washing by centrifugation before pipetting into chamber slides or multiwell dishes for imaging.

**MCF 10A, low confluence**



**MCF 10A, high confluence**



**(E) Confluence affects cytoplasmic vs. nuclear localization of YAP protein.**

Confocal IF analysis of low confluence MCF 10A cells (left) and high confluence MCF 10A (right) using YAP (D8H1X) XP® Rabbit mAb #14074 (green). Blue pseudocolor in lower images = DRAQ5® #4084 (fluorescent DNA dye). Increased nuclear and decreased cytoplasmic localization of YAP protein is seen in low confluence (proliferating) cells.

## Fixation

### Tissue

IF-F samples that have been fresh frozen and cryostat sectioned should now be treated with fixative. Alternatively, tissue samples may be preserved by fixing first with transcardial perfusion, post-fixation, and cryo-preservation steps, followed by freezing and sectioning.

For fluorescent IHC, tissue is preserved via perfusion as indicated above for IF-F, followed by paraffin embedding and sectioning. Prior to incubation with antibody, the sections must be deparaffinized with xylene and treated by heat or microwave in an antigen/epitope retrieval step.

### Cells

In IF-IC, the samples should be rapidly fixed such that cellular structures and target localization are faithfully preserved, while also allowing antibodies to recognize and report their target. This is accomplished by quickly exchanging media for fixative solution. Different approaches to fixation in IF-IC are available (see below), and the optimal fixation method depends on the cell type, target protein, and the antibody being used.

### Choice of Fixative

Aldehyde-based fixatives such as formaldehyde, formalin (a mixture of dissolved formaldehyde with a lower percentage of methanol), and glutaraldehyde are most commonly used. For most antibodies, CST recommends fixation with 4% formaldehyde. Aldehydes react with and crosslink cellular proteins, stabilizing and hardening the sample. The degree of crosslinking is dependent upon conditions (see below). Aldehydes cross the plasma membrane and fix soluble proteins better than alcohols, but some targets can lose their antigenicity with aldehyde crosslinking.

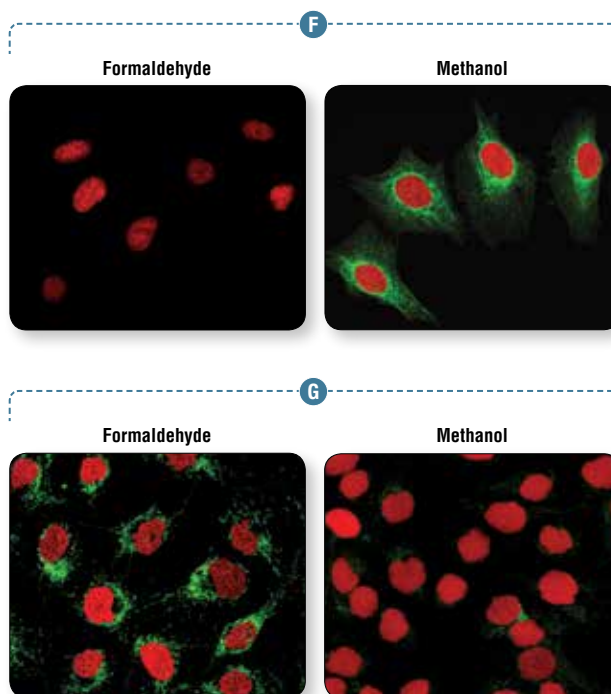
Dehydrating/denaturing fixatives displace water around cellular macromolecules, resulting in their denaturation and precipitation in situ. Denaturation of the target protein may expose normally buried epitopes, making this approach advantageous for some antibodies. **(F)** However, dehydrating fixatives are less suited for soluble targets and modification state-specific antibodies such as phospho-antibodies. **(G)** Check the product datasheet for optimal fixation method.

### Fixation Incubation Time

The degree of cross-linking, and hence the preservation of IF signal, is dependent upon time of fixation. In the following experiment, the concentration of formaldehyde was held at 4% and the duration of fixation was varied. Using fluorescence intensity as a readout, fixation for 15 min provides optimal signal for S6 Ribosomal Protein #2217. Incomplete fixation (5 min) results in delocalization of signal, and loss of signal intensity at longer fixation times. **(H)**

### What fix/perm conditions do I use when multiplexing?

If you are multiplexing with antibodies that call for different CST protocols (IF Standard vs. IF Methanol or IF Methanol-perm), you may need to prioritize which antibody to use at its optimal conditions. Performing a small-scale test run comparing different protocols may be informative before scaling up your experiments.



**Optimal fixative may vary depending on antibody/target.** Keratin 8/18 (C51) Mouse mAb #4546 works best with methanol fixation **(F)**. Confocal IF analysis of HeLa cells fixed with formaldehyde (left) or methanol (right) using #4546 (green). Red = Propidium Iodide (PI)/RNase Staining Solution #4087. AIF (D39D2) XP® Rabbit mAb #5318 works best with formaldehyde fixation **(G)**. Confocal IF analysis of HeLa cells fixed with formaldehyde (left) or methanol (right) using #5318 (green). Red = Propidium Iodide (PI)/RNase Staining Solution #4087.

### TROUBLESHOOTING TIP

Regularly replace old stocks of aldehyde fixative and prepare fresh dilutions; old solutions can contribute to elevated autofluorescence.



## Permeabilization

### Choice of Detergents or Alcohols

If a crosslinking fixative is used, the plasma membrane will still be intact, making intracellular targets inaccessible to antibodies. Therefore, permeabilization should be performed after crosslinking fixation for both IF-IC and IF-F unless your antibodies recognize extracellular epitopes. CST's IF Standard protocol (pg. 12) incorporates Triton® X-100 permeabilization after fixation with the blocking step (see next section). Triton and other detergents such as NP-40, TWEEN®, Saponin, Digitonin and DOTMAC remove different molecules from cellular membranes and create variable "pore" sizes to allow antibody access. Alternatively, alcohol permeabilization with ethanol or methanol may be performed after the fixation step. This method combines the rapid fixation of crosslinking fixatives with intermediate denaturation. This can improve signal for certain targets, particularly those associated with organelles or the cytoskeleton. As with fixation, the optimal permeabilization method varies depending on the antibody. Methanol permeabilization improves performance of some antibodies. **(I)** Consult product datasheet for recommended method.

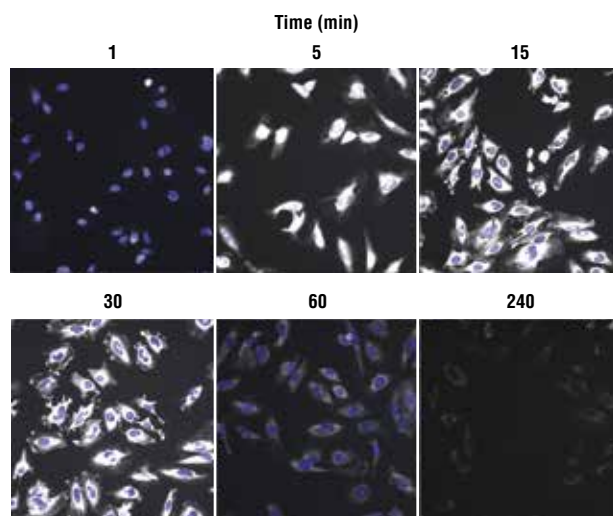
## Blocking

### Choice of Blocking Reagent

The blocking step reduces background signal caused by non-specific binding of primary and secondary antibodies to sites other than their intended target. Ideally, the blocking reagent will occupy "sticky" sites in the sample, such as exposed charged or hydrophobic protein surfaces, without interfering with primary/secondary antibody recognition of target epitopes, maximizing signal-to-noise ratio (S/N). The best blocking agent for the job varies by antibody; check product datasheet for recommendations. Our most common recommendation is 5% normal goat serum (or serum from the same species as the secondary antibody) in PBS + 0.3% Triton® X-100. CST also offers Image-IT® FX Signal Enhancer #11932, an alternate blocking reagent that reduces nonspecific charge-based interactions between the fluorophore and fixed sample. Ultimately, however, choosing high-quality antibodies that have been tested and approved for the application of choice will go a long way in reducing background, in concert with blocking.

### TROUBLESHOOTING TIP

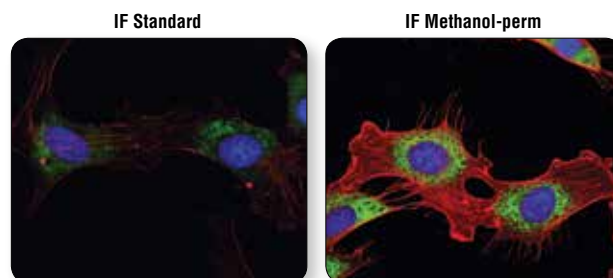
- Take care that the solution does not spill off the edge of the glass or dry out.
- Perform IF incubations in an airtight, light-blocking container lined with parafilm, and with a damp piece of paper towel to maintain a humid environment and protect fluorophore from photodamage.



**(H) Confocal IF analysis of HeLa cells fixed with formaldehyde for varying times as indicated, using S6 Ribosomal Protein (5G10) Rabbit mAb #2217 (white). Blue = Hoechst 33342 #4082 (fluorescent DNA dye).**

### Accounting for Autofluorescence

*Autofluorescence is a phenomenon where materials including tissues and cells emit background fluorescence in the absence of any experimental fluorophore, leading to degradation of signal-to-noise ratio. Sources of autofluorescence include NADH/NADPH, flavins and flavoproteins, the extracellular matrix proteins elastin and collagen, and lipofuscin, an insoluble lysosomal polymer that accumulates in aging tissues. The level of autofluorescence can vary by cell/tissue type and with choice of fixative. Autofluorescence tends to be strongest in shorter wavelength channels (green/yellow/orange emission). Although there are reported approaches to reduce autofluorescence in tissue such as short wavelength irradiation or quenching with sodium borohydride or pigments, these will also affect experimental fluorophores and target antigens in the sample as well. If you are concerned your sample may be autofluorescent, particularly for a dim signal, you can perform a control to compare intensities with and without antibody/fluorophore.*



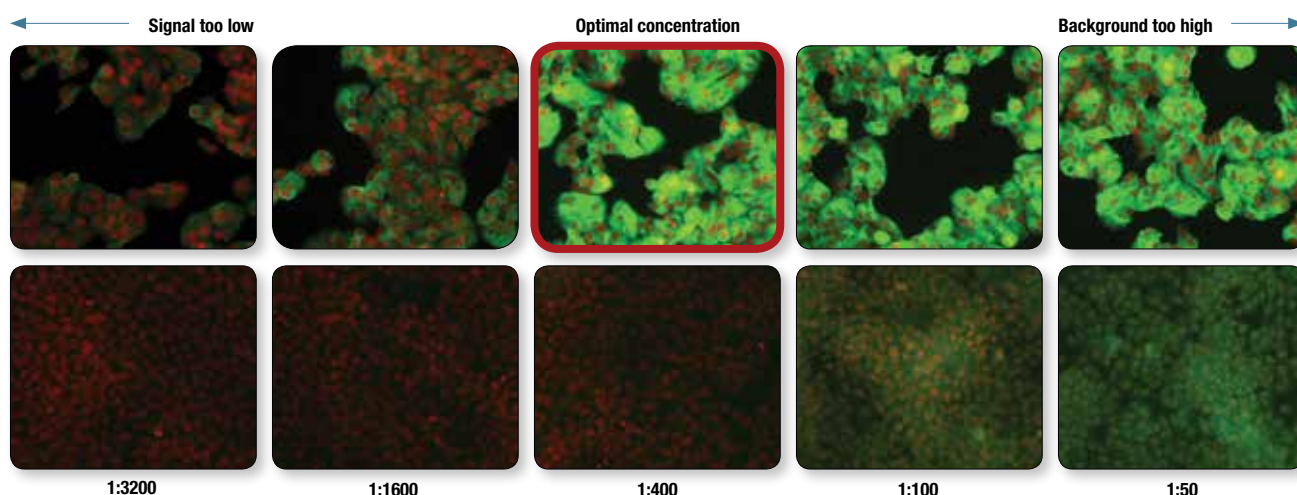
**(I) PDI (C81H6) Rabbit mAb #3501 and B-Actin (8H10D10) Mouse mAb #3700 work best with methanol permeabilization:** Confocal IF analysis of NIH/3T3 cells, permeabilized with 0.3% Triton™ X-100 (left) or methanol (right), using #3501 (green) and #3700 (red). Blue pseudocolor = DRAQ5® #4084 (fluorescent DNA dye).

# Immunostaining

## Primary Antibody Dilution

It is important to use an antibody at its recommended dilution in order to obtain a high signal-to-noise ratio (S/N). If the antibody is applied at too low of a concentration, the fluorescence signal will be too dim to distinguish from background noise. On the other hand, an excessively high concentration will contribute to background staining, decreasing S/N. CST scientists routinely perform titrations using positive and negative cell lines to provide you with the recommended dilution that gives optimal signal with minimal background staining. An example of an antibody titration is shown below, with the optimal concentration/recommended dilution outlined in red. The recommended antibody dilution for each IF-validated CST antibody is provided on its datasheet. Even though the sample has been blocked in the previous step, antibodies are always diluted and applied to samples in a dilution buffer containing a protein carrier (BSA). **(J)**

All of our recommended primary antibody dilutions are based upon overnight incubation at 4°C. This does not mean that CST antibodies will not work with a shorter incubation period, which is often used for automated platforms. The concentration of primary antibody may be increased to compensate for shorter incubation times; however, this will increase costs. The following experiment illustrates the dependence of signal intensity for anti-Vimentin #5741 on both incubation time and temperature. Optimal signal levels can be seen at the recommended incubation conditions of 4°C/overnight, while shorter incubation times at higher temperatures yield signal that is significantly lower. **(K)**



**(J) Optimization of MUC1 (D908K) XP® Rabbit mAb #14161 concentration for IF analysis:** IF analysis of ZR-75-1 cells (MUC1-positive, upper) and HCT 116 cells (MUC1-negative, lower) using #14161 (green) at the indicated dilutions. Red = Propidium Iodide (PI)/RNase Staining Solution #4087.

## Washing and Secondary Incubation

Thorough washing after fixation, in between primary and secondary incubations and/or counterstaining steps is critical for reducing background and increasing S/N. The antigen-binding region of an IF-approved antibody will have high affinity for its target epitope, but other regions of the antibody may still have lower affinity, non-specific interactions that could contribute to increased background. Washing the sample (3X for 5 min in PBS) helps to remove unbound antibody, prevent formation of primary-secondary immune complexes, and reduce unwanted background. Diluted secondary is applied for 1–2 hr at room temperature, protected from light.

## Counterstaining (Optional)

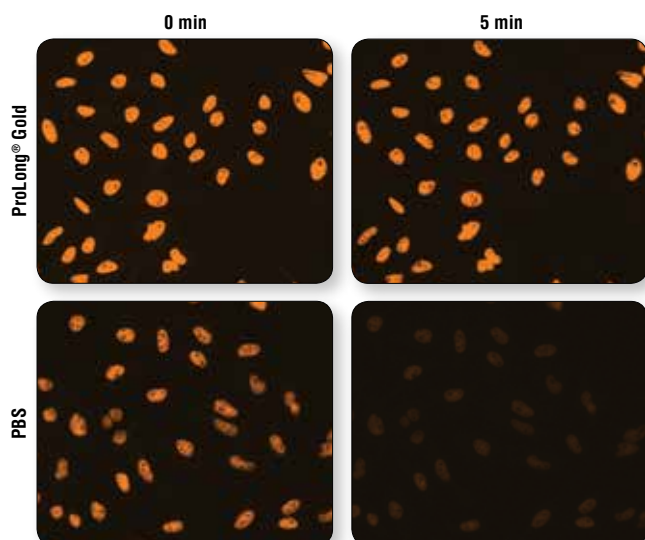
IF-IC detection is ideal for investigating changes in subcellular localization because protein location can be analyzed in relation to labeled organelle markers, other proteins of interest, or other cellular structures. Common counterstains used at CST are phalloidins (i.e., Alexa Fluor® 488 Phalloidin #8878), to label actin filaments, DRAQ5® #4084 to label DNA/nuclei, or MitoTracker® Red CMXRos #9082 to label mitochondria. **(L)** A list of Cellular Dyes/kits available from CST is on page 15. When using a counterstain, apply the same considerations of emission profile as you would with multiplexed antibodies to avoid bleed-through.

## Mount Sample

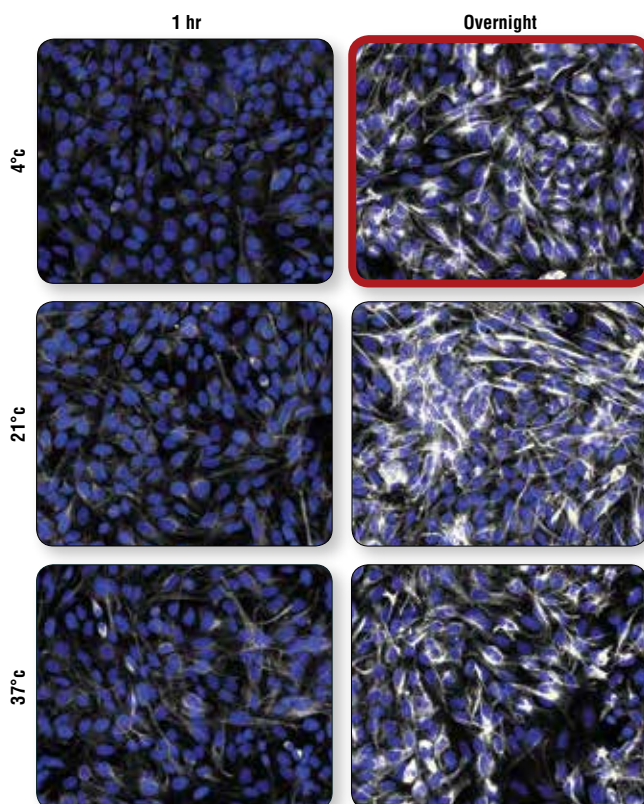
Exposure to high-energy light during image acquisition leads to oxidative damage and photobleaching of fluorophores. Moreover, crosslinking fixation is slowly reversible in aqueous media, potentially leading to delocalization of signal. Replacing the final PBS wash with a mounting agent such as ProLong® Gold Antifade Reagent #9071 provides several benefits: 1) displacing water to prevent crosslinker depolymerization, 2) curing (hardening) to stabilize the sample, 3) more closely matching the refractive index of the optics (for oil immersion objective lenses), and 4) reduction of photobleaching by free radical scavenging. The sample may be affixed to a glass slide by the mounting media and after it is cured, sealed with nail polish to prevent drying. The example below shows how the photobleaching effects of illumination can be reduced by ProLong® Gold. **(M)**

## Data Collection and Imaging

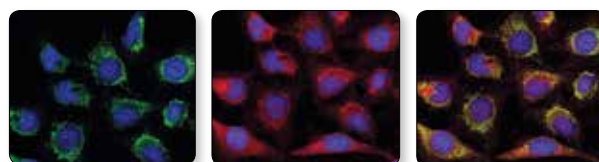
When performing IF analysis, it is prudent to consider imaging parameters early in the process, ideally with pilot experiments. Typically, these imaging parameters include exposure time, excitation intensity (e.g., by using neutral density filters), and camera gain. Parameters should be adjusted so that the signal is in the dynamic range of the camera and pixels are not saturated, while ensuring S/N is reasonably high. If you will be incorporating positive and/or negative controls, the same imaging protocol should be used for both sets of data, and images of the same antibody should always be shown using the same post-acquisition processing and analysis. Many journals have detailed guidelines for image acquisition and processing.



**(M) ProLong® Gold Antifade Reagent #9071 reduces photobleaching.** HeLa cells stained with ASH2L (D93F6) XP® Rabbit mAb #5019 and Anti-rabbit IgG (Alexa Fluor® 594) Conjugate and were mounted with #9071 (upper) or PBS (bottom) and imaged at the beginning (left) or end of 5 min illumination with epifluorescent excitation light.



**(K) Confocal IF analysis of vimentin-positive SNB-19 using Vimentin (D21H3) XP® Rabbit mAb #5741 (white).** Incubation of primary antibody at its recommended dilution was performed at 4°C, 21°C, or 37°C for 1 hr or overnight. Conditions recommended by CST for primary incubation (4°C overnight) yield maximum signal with little background (outlined in red). Blue = Hoechst 33342 #4082 (fluorescent DNA dye)



**(L) AIF colocalizes with mitochondria:** Confocal IF analysis of HeLa cells using AIF (D39D2) XP® Rabbit mAb #5318 (green), showing colocalization (yellow) with mitochondria that were labeled with MitoTracker® Red CMXRos #9082 (red). Blue pseudocolor = DRAQ5® #4084 (fluorescent DNA dye).

# Protocols

**IMPORTANT:** Please refer to the APPLICATIONS section on the front page of product datasheet or product webpage to determine if this product is validated and approved for use on cultured cell lines (IF-IC) or frozen tissue sections (IF-F). Alternatively, antibodies that have been approved for IHC can be used in paraffin-embedded sections (see [www.cellsignal.com/CSTprotocols](http://www.cellsignal.com/CSTprotocols) for our mIHC/Paraffin protocol). Please see product datasheet or product webpage for appropriate antibody dilution. The following protocols accommodate both indirect and direct IF.

**NOTE:** Some CST antibodies work optimally using an alternate protocol. Please see product datasheet for product-specific recommendations.

## A. Solutions and Reagents

**NOTE:** Prepare solutions with reverse osmosis deionized (RODI) or equivalent grade water.

- **1X Phosphate Buffered Saline (PBS):** To prepare 1 L 1X PBS pH 8.0, add 100 mL 10X PBS pH 7.4 (1.37M NaCl, 27mM KCl, 101 mM  $\text{Na}_2\text{HPO}_4$ , 18 mM  $\text{KH}_2\text{PO}_4$ ) to 900 mL dH<sub>2</sub>O, mix. NOTE: adjust pH to 8.0.
- **Formaldehyde: 16%, methanol free, Polysciences, Inc. (cat# 18814),** use fresh and store opened vials at 4°C in dark. Dilute 1 in 4 in 1X PBS to make a 4% formaldehyde solution.
- **Blocking Buffer:** (1X PBS/5% normal serum/0.3% Triton® X-100): To prepare 10 ml, add 0.5 ml normal serum from the same species as the secondary antibody (e.g., Normal Goat Serum (#5425) to 9.5 ml 1X PBS) and mix well. While stirring, add 30 µl Triton® X-100.
- **Antibody Dilution Buffer:** (1X PBS/1% BSA/0.3% Triton® X-100): To prepare 10 ml, add 30 µl Triton® X-100 to 10 ml 1X PBS. Mix well then add 0.1g BSA (#9998), mix.
- **Fluorophore-conjugated Secondary Antibodies:** (Anti-mouse #4408, #4409, #8890, #4410) (Anti-rabbit #4412, #4413, #8889, #4414) (Anti-rat #4416, #4417, #4418).
- **ProLong® Gold Antifade Reagent (#9071),** ProLong® Gold Antifade Reagent with DAPI (#8961).
- **Methanol:** 100%, ice-cold (for IF-Methanol-fixed and IF-Methanol-perm protocols)

**Proceed to one of the following protocols.**

## Immunofluorescence General Protocol (IF Standard)

### B. Specimen Preparation

#### I. Cultured Cell Lines (IF-IC)

**NOTE:** Cells should be grown, treated, fixed and stained directly in multi-well plates, chamber slides or on coverslips.

1. Aspirate liquid, then cover cells to a depth of 2–3 mm with 4% formaldehyde diluted in warm PBS.

**NOTE:** Formaldehyde is toxic, use only in a fume hood.

2. Allow cells to fix for 15 min at room temperature.
3. Aspirate fixative, rinse three times in 1X PBS for 5 min each.
4. Proceed with Immunostaining (Section C).

#### II. Frozen/Cryostat Sections (IF-F)

1. For fixed frozen tissue proceed with Immunostaining (Section C).
2. For fresh, unfixed frozen tissue, fix immediately, as follows:  
Cover sections with 4% formaldehyde diluted in warm 1X PBS.  
Allow sections to fix for 15 min at room temperature.  
Rinse slides three times in PBS for 5 min each.  
Proceed with Immunostaining (Section C).

## C. Immunostaining

**NOTE:** All subsequent incubations should be carried out at room temperature unless otherwise noted in a humid light-tight box or covered dish/plate to prevent drying and fluorochrome fading.

1. Block specimen in blocking buffer for 60 min.
2. While blocking, prepare primary antibody by diluting as indicated on datasheet in antibody dilution buffer.
3. Aspirate blocking solution, apply diluted primary antibody.
4. Incubate overnight at 4°C.
5. Rinse three times in 1X PBS for 5 min each.

**NOTE:** If using a fluorophore-conjugated primary antibody, then skip to Section C, Step 8.

6. Incubate specimen in fluorochrome-conjugated secondary antibody diluted in antibody dilution buffer for 1–2 hr at room temperature in the dark.
7. Rinse three times in 1X PBS for 5 min each.
8. Coverslip slides with ProLong® Gold Antifade Reagent (#9071) or ProLong® Gold Antifade Reagent with DAPI (#8961).
9. For best results, allow the mounting reagent to cure overnight at room temperature. For long-term storage, store slides flat at 4°C protected from light.



## Immunofluorescence Protocol with Methanol Fixation (IF Methanol-fixed)

**IMPORTANT:** Please refer to the APPLICATIONS section on the front page of the datasheet to determine if this product is validated and approved for use with this protocol.

### B. Specimen Preparation

#### Cultured Cell Lines (IF-IC)

**NOTE:** Cells should be grown, treated, fixed and stained directly in multi-well plates, chamber slides or on coverslips.

1. Aspirate liquid, then cover cells to a depth of 2–3 mm with ice-cold 100% methanol.
2. Allow cells to fix for 15 min at -20°C.
3. Aspirate fixative, rinse three times in PBS for 5 min each.
4. Proceed with Immunostaining (Section C).

### C. Immunostaining

**NOTE:** All subsequent incubations should be carried out at room temperature unless otherwise noted in a humid light-tight box or covered dish/plate to prevent drying and fluorochrome fading.

1. Block specimen in Blocking Buffer for 60 min.
2. While blocking, prepare primary antibody by diluting as indicated on datasheet in Antibody Dilution Buffer.
3. Aspirate blocking solution, apply diluted primary antibody.
4. Incubate overnight at 4°C.
5. Rinse three times in PBS for 5 min each.

**NOTE:** If using a fluorochrome-conjugated primary antibody, then skip to Section C, Step 8.

6. Incubate specimen in fluorochrome-conjugated secondary antibody diluted in Antibody Dilution Buffer for 1–2 hours at room temperature in dark.
7. Rinse in PBS as in step 5.
8. Coverslip slides with ProLong® Gold Antifade Reagent (#9071) or ProLong® Gold Antifade Reagent with DAPI (#8961).
9. For best results, examine specimens immediately using appropriate excitation wavelength. For long-term storage, store slides flat at 4°C protected from light.

## Immunofluorescence Protocol with Methanol Permeabilization (IF Methanol-perm)

### B. Specimen Preparation

#### I. Cultured Cell Lines (IF-IC)

**NOTE:** Cells should be grown, treated, fixed and stained directly in multiwell plates, chamber slides or on coverslips.

1. Aspirate liquid, then cover cells to a depth of 2–3 mm with 4% formaldehyde in PBS.

**NOTE:** Formaldehyde is toxic, use only in fume hood.

2. Allow cells to fix for 15 min at room temperature.
3. Aspirate fixative, rinse three times in PBS for 5 min each.
4. Proceed with Immunostaining (Section C).

#### II. Frozen/Cryostat Sections (IF-F)

1. For fixed frozen tissue proceed with Immunostaining (Section C).
2. For fresh, unfixed frozen tissue, please fix immediately, as follows:
  1. Cover sections with 4% formaldehyde in PBS.

**NOTE:** Formaldehyde is toxic, use only in fume hood.

2. Allow sections to fix for 15 min at room temperature.
3. Rinse slides three times in PBS for 5 min each.
4. Proceed with Immunostaining (Section C).

### C. Immunostaining

**NOTE:** All subsequent incubations should be carried out at room temperature unless otherwise noted in a humid light-tight box or covered dish/plate to prevent drying and fluorochrome fading.

1. Methanol Permeabilization Step: Cover cells or tissue sections with ice-cold 100% methanol (use enough to cover completely to a depth of 3–5 mm, DO NOT LET DRY), incubate in methanol for 10 min at -20°C, rinse in PBS for 5 min.
2. Block specimen in Blocking Buffer for 60 min.
3. While blocking, prepare primary antibody by diluting as indicated on datasheet in Antibody Dilution Buffer.
4. Aspirate blocking solution, apply diluted primary antibody.
5. Incubate overnight at 4°C.
6. Rinse three times in PBS for 5 min each.

**NOTE:** If using a fluorochrome-conjugated primary antibody, then skip to Section C, Step 9.

7. Incubate specimen in fluorochrome-conjugated secondary antibody diluted in Antibody Dilution Buffer for 1–2 hours at room temperature in dark.
8. Rinse in PBS as in step 6.
9. Coverslip slides with ProLong® Gold Antifade Reagent (#9071) or ProLong® Gold Antifade Reagent with DAPI (#8961).
10. For best results examine specimens immediately using appropriate excitation wavelength. For long term storage, store slides flat at 4°C protected from light.

# IF Troubleshooting Guide

WEAK OR NO SIGNAL	
Possible cause	Recommendations
Inappropriate storage of samples. Signal may fade if fluorophores are exposed to light for extended periods of time.	Perform incubations and store samples in the dark. Samples may be mounted in an anti-fade solution, such as ProLong® Gold Antifade Reagent #9071. Samples should be imaged immediately following mounting for the best results.
Cell/tissue samples stored for too long	Use freshly prepared slides/plates to avoid loss of antigenicity.
Inadequate fixation	Consult the CST product datasheet for the recommended protocol; remove media and quickly/thoroughly wash in fixative immediately after treatment. For phospho-specific antibodies, use at least 4% formaldehyde to inhibit endogenous phosphatases.
Incorrect antibody dilution (antibody too dilute)	Consult the CST product datasheet or cellsignal.com for the recommended antibody dilution.
Not using the recommended incubation time	Primary antibody incubation according to a rigorously tested protocol provides consistent, reliable results. CST antibodies have been developed and validated for optimal results when incubated at 4°C overnight.
Inappropriate testing model	If possible, protein expression should be confirmed by western blot or other means.
Target protein not induced properly	Optimal treatment conditions and controls should be determined for each target.
Low expression of protein of interest	Modify detection approach; consider signal amplification, or pair with a brighter fluorophore.
Wrong permeabilization method	Consult datasheet for recommended protocol.
Incorrect use of secondary antibody	Use recommended concentration and check secondary antibody is matched to host species of the primary antibody.
Wrong excitation wavelength	Ensure illumination and detection (laser/excitation/emission filter) matches excitation wavelength of fluorophore(s).
Low signal in multiplexed IHC	Optimization of deparaffination, antigen retrieval and signal amplification methods.

HIGH BACKGROUND	
Cause	Solution
Sample autofluorescence	Use unstained samples as controls to check autofluorescence levels. Check datasheet for correct fixation reagent. Old fixative may autofluoresce; replace old formaldehyde stocks and prepare fresh dilutions. Use EM-grade glutaraldehyde freshly diluted from ampules. Choose longer wavelength channels for low-abundance targets.
Insufficient blocking	Use normal serum from the same species as the secondary antibody used. Consider a charge-based blocker such as Image-iT® FX Signal Enhancer #11932, depending on the source of background.
Incorrect antibody dilution (primary or secondary antibody too concentrated)	Consult the CST product datasheet or cellsignal.com for the recommended antibody dilution.
Samples dried out	It is vital that sample remains covered in liquid throughout the staining procedure.
Insufficient washing	Wash to remove excess fixative, excess secondary antibody, and loosely bound, non-specific antibody interactions.
Secondary cross-reactivity	Use isotype control antibodies to determine whether your secondary antibody is cross-reacting.
Non-specific antibody binding	If available, compare to knockdown (siRNA) or knockout cells, or compare to cells known to express higher or lower levels of the target antigen.

# Recommended Application Solutions

Cell Signaling Technology offers reagents and kits to support your immunofluorescence experiments. These products are used by CST to validate primary antibodies for immunofluorescence.

## Blocking Reagents

BSA #9998  
Normal Goat Serum #5425  
Image-iT® FX Signal Enhancer #11932

## Primary Antibody Conjugates

Visit [cellsignal.com/IF](http://cellsignal.com/IF)

## Secondary Antibodies

Anti-mouse IgG (H+L), F(ab')<sub>2</sub> Fragment (Alexa Fluor® 488 Conjugate) #4408  
Anti-mouse IgG (H+L), F(ab')<sub>2</sub> Fragment (Alexa Fluor® 555 Conjugate) #4409  
Anti-mouse IgG (H+L), F(ab')<sub>2</sub> Fragment (Alexa Fluor® 594 Conjugate) #8890  
Anti-mouse IgG (H+L), F(ab')<sub>2</sub> Fragment (Alexa Fluor® 647 Conjugate) #4410  
Anti-rabbit IgG (H+L), F(ab')<sub>2</sub> Fragment (Alexa Fluor® 488 Conjugate) #4412  
Anti-rabbit IgG (H+L), F(ab')<sub>2</sub> Fragment (Alexa Fluor® 555 Conjugate) #4413  
Anti-rabbit IgG (H+L), F(ab')<sub>2</sub> Fragment (Alexa Fluor® 647 Conjugate) #4414  
Anti-rabbit IgG (H+L), F(ab')<sub>2</sub> Fragment (Alexa Fluor® 594 Conjugate) #8889  
Anti-rat IgG (H+L), (Alexa Fluor® 488 Conjugate) #4416  
Anti-rat IgG (H+L), (Alexa Fluor® 555 Conjugate) #4417  
Anti-rat IgG (H+L), (Alexa Fluor® 647 Conjugate) #4418

## Cellular Dyes

Alexa Fluor® 488 Phalloidin #8878  
Alexa Fluor® 555 Phalloidin #8953  
Alexa Fluor® 647 Phalloidin #8940  
DyLight™ 554 Phalloidin #13054  
DyLight™ 488 Phalloidin #12935

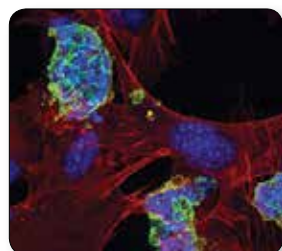
DyLight™ 650 Phalloidin #12956  
DyLight™ 350 Phalloidin #12848  
DyLight™ 594 Phalloidin #12877  
MitoTracker® Deep Red FM #8778  
MitoTracker® Red CMXRos #9082  
Hoechst 33342 #4082  
DAPI #4083  
DRAQ5® #4084  
Propidium Iodide (PI)/RNase Staining Solution #4087

## IF Kits

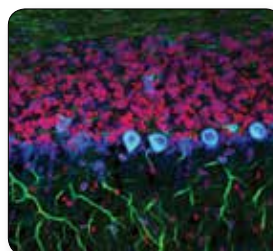
Immunofluorescence Application Solutions Kit #12727  
Cellular Localization IF Antibody Sampler Kit #4753  
Organelle Localization IF Antibody Sampler Kit #8653  
ER and Golgi-Associated Marker Proteins Antibody Sampler Kit #12718  
PathScan® Apoptosis and Proliferation Multiplex IF Kit #7851  
PathScan® EMT Duplex IF Kit #7771  
Neuronal Marker IF Antibody Sampler Kit #8572  
StemLight™ iPS Cell Reprogramming Antibody Kit #9092

## Mounting Reagents

ProLong® Gold Antifade Reagent #9071  
ProLong® Gold Antifade Reagent with DAPI #8961



Confocal IF analysis of mouse embryonic stem cells growing on mouse embryonic fibroblast (MEF) feeder cells using SSEA1 (MC480) Mouse mAb #4744 (green). Actin filaments have been labeled with DyLight™ 554 Phalloidin #13054 (red). Blue pseudocolor = DRAQ5® (fluorescent DNA dye).



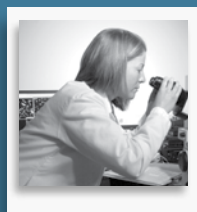
Confocal IF analysis of rat brain using S6 Ribosomal Protein (54D2) (Alexa Fluor® 647) Mouse mAb #5548 (blue pseudocolor) and β3-Tubulin (D71G9) XP® Rabbit mAb #5568 (green). Red = Propidium Iodide (fluorescent DNA dye).

For more data on experimental controls visit

[www.cellsignal.com/ifvalidation](http://www.cellsignal.com/ifvalidation)

# Technical Support

We hope this guide is a helpful resource for performing immunofluorescence in your own lab. Cell Signaling Technology prides itself in providing you with exceptional customer service and support, and we are happy to share our experience with you. Since all of our antibodies are produced in-house, the same scientists who develop and assay these reagents are available as technical resources for our customers. These scientists can be contacted directly and will personally provide technical assistance to you, our customer.



**FRONT COVER PHOTO:**  
Tara, Imaging Specialist, has  
been with CST since 2008.

## By Region

### UNITED STATES

Orders: 877-616-2355 | [orders@cellsignal.com](mailto:orders@cellsignal.com)  
Support: 877-678-8324 | [support@cellsignal.com](mailto:support@cellsignal.com)  
[www.cellsignal.com](http://www.cellsignal.com)

### CHINA

Tel: +86-21-58356288  
Support (China): 4006-473287/GreatQ | [tech@cst-c.com.cn](mailto:tech@cst-c.com.cn)  
Support (Asia Pacific): [csttechasia@cst-c.com.cn](mailto:csttechasia@cst-c.com.cn)  
[www.cst-c.com.cn](http://www.cst-c.com.cn)

### EUROPE, MIDDLE EAST & AFRICA

Tel: +31 (0)71 720 0200  
Support: [eusupport@cellsignal.eu](mailto:eusupport@cellsignal.eu)  
[www.cellsignal.com](http://www.cellsignal.com)

### JAPAN

Tel: 03-3295-1630 | Support: [info@cstj.co.jp](mailto:info@cstj.co.jp)  
[www.cstj.co.jp](http://www.cstj.co.jp)

### ORDER INFORMATION

Find order information online at  
[www.cellsignal.com/orderinfo](http://www.cellsignal.com/orderinfo)

♻️ Printed on recycled paper (25% post-consumer waste fiber) using vegetable inks and processed chlorine free.

© 2017 Cell Signaling Technology, Inc. All rights reserved.

Cell Signaling Technology, CST, PathScan, SignalSlide, SignalStain, StemLight, and XP are trademarks of Cell Signaling Technology, Inc. Alexa Fluor is a registered trademark of Life Technologies Corporation. DRAQ5 is a registered trademark of Biostatus Limited. DyLight is a trademark of Thermo Fisher Scientific, Inc. and its subsidiaries. Image-iT is a registered trademark of Life Technologies Corporation. MitoTracker is a registered trademark of Molecular Probes, Inc. Triton is a trademark of The Dow Chemical Company. TSA is a registered trademark of PerkinElmer.

Alexa Conjugates: This product is provided under an intellectual property license from Life Technologies Corporation. The transfer of this product is contingent on the buyer using the purchased product solely in research, including use with HCS or other automated imaging applications but excluding use in combination with DNA microarrays. The buyer must not sell or otherwise transfer this product or its components for (a) diagnostic, therapeutic or prophylactic purposes; (b) testing, analysis or screening services, or information in return for compensation on a per-test basis; (c) manufacturing or quality assurance or quality control, or (d) resale, whether or not resold for use in research. For information on purchasing a license to this product for purposes other than as described above, contact Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, CA 92008 USA or [outlicensing@lifetech.com](mailto:outlicensing@lifetech.com).

**For Research Use Only. Not For Use in Diagnostic Procedures.**

17-APS-011-BR01-E1