

Tyramide signal amplification (TSA) typically provides 2-3 logs of sensitivity enhancement over standard detection methods (for example, fluorophore-labeled secondary antibodies or enzymatic detection with a chromogenic substrate) while maintaining resolution and allowing reduced consumption of primary antibodies. Protocol optimization for TSA detection is critical for success with the technique.

## Important information

- Horseradish peroxidase (HRP) is necessary to drive the TSA labeling reaction.
- If present, endogenous peroxidases should be quenched to prevent non-specific detection.
- If HRP concentration is too high, formation of TSA dimers will be favored over the labeling reaction, leading to reduced signal and higher background. For this reason, it is critical to optimize dilution of primary antibody and HRP reagent.
- Use of a blocking reagent is highly recommended. PerkinElmer offers a blocking reagent that is optimal for use with TSA (catalog number FP1012).

- TSA stock solution should be diluted 1:50 in amplification diluent. Problems with excess signal can be resolved by using more dilute antibody solutions for detection. This approach ensures the highest specificity in detection.
- First time TSA users should apply this to a proven IHC system, and include an unamplified control slide.

## Suggested optimization procedure for first time use of TSA

TSA is ideal for detection of targets that have weak signals using standard detection methods. It enables use of much more dilute antibody solutions for improved specificity. If the signal is strong with standard detection methods, dilute the primary antibody until the signal begins to disappear and then proceed with TSA optimization in the same manner.

1. Primary antibody optimization	2. Introduction of HRP: Select an option	3. Detection
<p><b>Slide 1:</b> Dilution specified by current method or recommended by primary Ab manufacturer</p> <p><b>Slide 2:</b> 5-fold dilution from slide 1</p> <p><b>Slide 3:</b> 5-fold dilution from slide 2</p> <p><b>Slide 4:</b> 5-fold dilution from slide 3 (further dilution may be needed)</p> <p><b>Slide 5:</b> Negative control (no primary antibody)</p>	<p>Secondary antibody HRP conjugates</p> <p>Anti-mouse-HRP, (catalog number NEF822001EA)</p> <p>Anti-rabbit-HRP, (catalog number NEF812001EA)</p> <p>Recommended starting range of 1:500 to 1:2000.</p> <p>Further dilution may be needed</p> <p>SA-HRP reagent (catalog number NEL750001EA)</p> <p>1:1250-1:2500 dilution</p> <p>Anti-fluorescein-HRP (catalog number NEF710001EA)</p> <p>1:100-1:500 dilution</p> <p>HRP reagents from other manufacturers</p> <p>Begin with recommended range for slide applications.</p> <p>Further dilution may be needed.</p>	<p>TSA 1:50 in amplification diluent</p>

## TSA Stock Solutions

Kit	Catalog Numbers	Solvent	Volume per tube
TSA Plus Biotin	NEL749A, NEL749B	DMSO	300 µL
TSA Plus DNP, 25-75 slides	NEL746B, NEL747B	DMSO	150 µL
TSA Plus DNP, 50-150 slides	NEL746A, NEL747A	DMSO	300 µL
TSA Plus Fluorescence Kits	Kits beginning with NEL741-745	DMSO	150 µL
TSA Plus Multi-Fluor Kits	Kits beginning with NEL752-756	DMSO	150 µL
TSA Plus Palette Kit	NEL760	DMSO	60 µL
TSA Biotin, 200-600 slides	NEL700	DMSO	1200 µL
TSA Biotin, 50-150 slides	NEL700A	DMSO	300 µL
TSA Fluorescence Kits, 100-300 slides	NEL701, NEL702, NEL703	DMSO	600 µL
TSA Fluorescein, 50-150 slides	NEL701A	DMSO	300 µL
TSA Cyanine 3 or 5, 50-150 slides	NEL704A, NEL705A	H <sub>2</sub> O	300 µL

Dilute stock solutions 1:50 in amplification diluent to make working solution. DMSO should be molecular biology or HPLC grade.

## Protocol for TSA IHC/ICC/IF

<b>Standard IHC</b>	Quench endogenous peroxidase activity (if needed) Block slides for 30 min. in TNB buffer at room temperature Incubate slides in primary antibody for 30-60 minutes at room temperature Wash slides 3X for 5 min. in TNT buffer at room temperature		
<b>Incorporation of HRP</b>	Incubate slides in HRP labeled secondary antibody for 30 min. at room temperature	Incubate slides in biotinylated secondary antibody 30-60 min. at room temperature	
		Wash slides 3X for 5 min. in TNT buffer Incubate slides in SA-HRP for 30 min.	
	Wash slides 3X for 5 min. in TNT buffer at room temperature		
<b>TSA Plus Amplification</b>	Incubate in Amplification Reagent working solution for 3 to 10 min. at room temperature		
	Wash slides 3X for 5 min. in TNT buffer at room temperature		
<b>Visualization</b>	<b>Direct Fluorescence</b>	<b>Indirect Fluorescence</b>	<b>Chromogenic</b>
		Incubate slides in fluorophore conjugate for 30 min.	Incubate slides in HRP or AP conjugate for 30 min.
	Wash slides 3X for 5 min. in TNT buffer at room temperature	Wash slides 3X for 5 min. in TNT buffer at room temperature	Wash slides 3X for 5 min. in TNT buffer at room temperature
	Counterstain and mount for fluorescence microscopy	Counterstain and mount for fluorescence microscopy	Add appropriate chromogen, counterstain and mount for microscopy

PerkinElmer, Inc.  
 940 Winter Street  
 Waltham, MA 02451 USA  
 P: (800) 762-4000 or  
 (+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2011, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

009671\_02

May, 2011