

**XFD532 tyramide**

 Catalog number: 11072  
 Unit size: 200 Slides

Component	Storage	Amount (Cat No. 11072)
XFD 532 tyramide	Freeze (< -15 °C), Minimize light exposure	200 Slides

**OVERVIEW**

Tyramide signal amplification (TSA) has proven to be a particularly versatile and powerful enzyme amplification technique with improved assay sensitivity. TSA is based on the ability of HRP, in the presence of low concentrations of hydrogen peroxide, to convert labeled tyramine-containing substrate into an oxidized, highly reactive free radical that can covalently bind to tyrosine residues at or near the HRP. The signal amplification conferred by the turnover of multiple tyramide substrates per peroxidase label translates ultrasensitive detection of low-abundance targets and the use of smaller amounts of antibodies and hybridization probes. In immunohistochemical applications, sensitivity enhancements derived from TSA method allow primary antibody dilutions to be increased to reduce nonspecific background signals, and can overcome weak immunolabeling caused by suboptimal fixation procedures or low levels of target expression. XFD 532 tyramide contains the Alexa Fluor® 532 fluorophore that can be readily detected with the Alexa Fluor® 532/ATTO 532 filter set (Alexa Fluor® is the trademark of ThermoFisher). XFD 532 tyramide has intense yellow-orange fluorescence color.

**AT A GLANCE**
**Protocol Summary**

1. Fix/permeabilize/block cells or tissue
2. Add primary antibody in blocking buffer
3. Add HRP-conjugated secondary antibody
4. Prepare tyramide working solution and apply in cells or tissue for 5-10 minutes at room temperature

**KEY PARAMETERS**
**Fluorescence microscope**

Emission	Cy3/TRITC filter set
Excitation	Cy3/TRITC filter set
Recommended plate	Black wall/clear bottom

**PREPARATION OF STOCK SOLUTIONS**

*Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles*

**Tyramide stock solution (200X)**

Add 100 µL of DMSO to the vial of XFD532 tyramide and mix well.

**Note:** Make single-use aliquots and store unused 200X stock solution at 2-8 °C, protected from light. Avoid repeat freeze-thaw cycles.

**PREPARATION OF WORKING SOLUTION**
**Tyramide working solution (1X)**

Add 100 µL of the tyramide stock solution into 20 mL of a buffer of

your choice containing 0.003% H<sub>2</sub>O<sub>2</sub>.

**Note:** For optimal performance, use Tris Buffer, pH=7.4.

**Note:** A 20 mL solution is good for 200 tests. The tyramide working solution should be used immediately and made fresh on the day of use. Avoid direct exposure to light.

**Secondary antibody-HRP working solution**

Make an appropriate concentration of secondary antibody-HRP working solution per the manufacturer's recommendations.

**SAMPLE EXPERIMENTAL PROTOCOL**

This protocol is applicable for both cells and tissues staining.

**Cell fixation and permeabilization**

1. Fix the cells or tissue with 3.7% formaldehyde or paraformaldehyde, in PBS at room temperature for 20 minutes.
2. Rinse the cells or tissue with PBS twice.
3. Permeabilize the cells with 0.1% Triton X-100 solution for 1-5 minutes at room temperature.
4. Rinse the cells or tissue with PBS twice.

**Tissue fixation, deparaffinization and rehydration**

Deparaffinize and dehydrate the tissue according to the standard IHC protocols. Perform antigen retrieval with the preferred specific solution/protocol as needed. A protocol can be found at:

<https://www.aatbio.com/resources/guides/paraffin-embedded-tissue-immunohistochemistry-protocol.html>

**Peroxidase labeling**

1. Optional: Quench endogenous peroxidase activity by incubating cell or tissue sample in peroxidase quenching solution (such as 3% hydrogen peroxide) for 10 minutes. Rinse with PBS twice at room temperature.
2. Optional: If using HRP-conjugated streptavidin, it is advisable to block endogenous biotins by biotin blocking buffer.
3. Block with preferred blocking solution (such as PBS with 1% BSA) for 30 minutes at 4 °C.
4. Remove blocking solution and add primary antibody diluted in recommended antibody diluent for 60 minutes at room temperature or overnight at 4 °C.
5. Wash with PBS three times for 5 minutes each.
6. Apply 100 µL of secondary antibody-HRP working solution to each sample and incubate for 60 minutes at room temperature.

**Note:** Incubation time and concentration can be varied depending on the signal intensity.

7. Wash with PBS three times for 5 minutes each.

**DISCLAIMER**

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.

**Tyramide labeling**

1. Prepare and apply 100 µL of Tyramide working solution to each sample and incubate for 5-10 minutes at room temperature.

**Note:** If you observe a non-specific signal, you can shorten the incubation time with Tyramide. You should optimize the incubation period using positive and negative control samples at various incubation time points. Or you can use a lower concentration of the tyramide reagent in the working solution.

2. Rinse with PBS three times.

**Counterstain and fluorescence imaging**

1. Counterstain the cell or tissue samples as needed. AAT provides a series of nucleus counterstain reagents as listed in Table 1. Follow the instruction provided with the reagents.

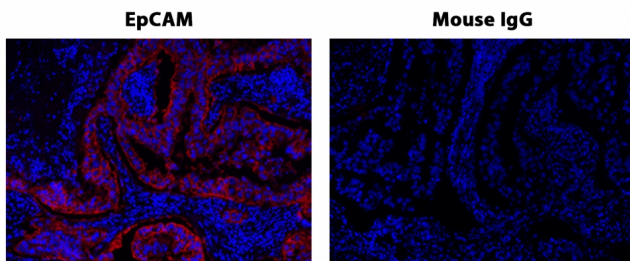
2. Mount the coverslip using a mounting medium with anti-fading properties.

**Note:** To ensure optimal results, it is recommended to use either ReadiUse™ microscope mounting solution (Cat. 20009) or FluoroQuest™ TSA/PSA Antifade Mounting Medium \*Optimized for Tyramide and Styramide Imaging\* (Cat. 44890) instead of Vectashield® mounting media. There are instances where Vectashield® mounting media may not be suitable for certain TSA/PSA conjugates.

3. Use the appropriate filter set to visualize the signal from the Tyramide labeling.

**Table 1.** Products recommended for nucleus counterstain

Cat#	Product Name	Ex/Em (nm)
17548	Nuclear Blue™ DCS1	350/461
17550	Nuclear Green™ DCS1	503/526
17551	Nuclear Orange™ DCS1	528/576
17552	Nuclear Red™ DCS1	642/660

**EXAMPLE DATA ANALYSIS AND FIGURES**

**Figure 1.** Fluorescence IHC of formaldehyde-fixed, paraffin-embedded human lung adenocarcinoma positive tissue using PSA™ amplified methods. Human lung adenocarcinoma positive tissue sections were stained with Mouse anti-EpCAM or Control Mouse IgG antibody and then incubated with polyHRP-labeled Goat anti-Mouse IgG secondary antibody followed by XFD 532 tyramide (Cat#11072).