

ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer

Catalog number: 45091
Unit size: 100 mL

Component	Storage	Amount (Cat No. 45091)
Component A: ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer 01	Refrigerated (2-8 °C)	50 mL
Component B: ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer 02	Refrigerated (2-8 °C), Minimize light exposure	50 mL

OVERVIEW

ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer is formulated to effectively remove fluorescence signals generated from tyramide or styramide amplification steps in immunofluorescence workflows. This reagent enables removal of both the fluorophore and bound antibody complex, allowing reuse of tissue sections for additional rounds of labeling without requiring heat-based antigen retrieval. The stripping process is performed at room temperature, offering a gentle and efficient alternative to traditional thermal methods. ReadiUse™ Stripping Buffer streamlines multiplexed TSA and PSA protocols by preserving tissue morphology and reducing assay complexity between cycles.

AT A GLANCE

1. Prepare the ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer working solution
2. Perform immunofluorescence with tyramide or styramide according to the manufacturer's instructions and image the tissue
3. Wash with PBS, incubate with ReadiUse™ TSA/PSA Stripping Buffer for 30–60 minutes at room temperature, then wash again
4. Proceed to the next immunofluorescence cycle starting from the blocking step

PREPARATION OF WORKING SOLUTION

ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer working solution:

Mix 1:1 ratio of ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer 01 (Component A) and ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer 02 (Component B) to make final working solution.

Note: For best results, mix both components prior to use.

Note: Mixing Component A and Component B may result in effervescence; exercise caution during mixing.

Note: Discard the unused working solution after each experiment. Do not store it.

SAMPLE EXPERIMENTAL PROTOCOL

1. Perform Immunofluorescence with tyramide or styramide on tissue samples as per manufacturer's recommendations.
2. Observe and record the image of the tissue samples.
3. Briefly wash the tissue samples with PBS buffer.
4. Add 100 µL of ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer directly to cover the sample and incubate at room temperature for 30 to 60 minutes.
5. Remove the ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer and wash the tissue samples with PBS.
6. Proceed to the next cycle of immunofluorescence from blocking step.

IMPORTANT NOTES

1. ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer completely removes the fluorescence from the previous cycle of tyramide or styramide reaction.
2. ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping may also remove residual primary and secondary antibodies, potentially reducing the need for heat-induced antigen retrieval in subsequent staining rounds. Users should optimize experimental conditions as needed for their specific samples and protocols.

EXAMPLE DATA ANALYSIS AND FIGURES

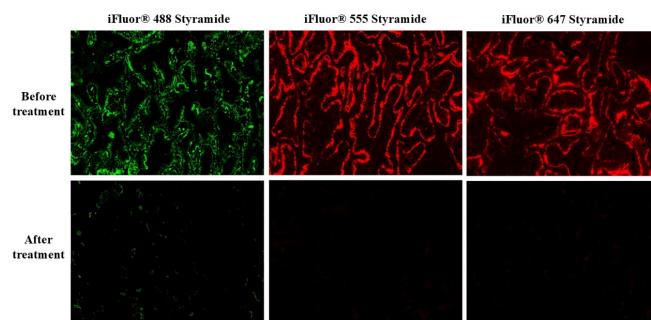


Figure 1. Fluorescence IHC of formaldehyde-fixed, paraffin-embedded human lung adenocarcinoma tissue using PSA™ and TSA amplification methods. Tissue sections were stained with rabbit anti-EpCAM antibody, followed by polyHRP-conjugated goat anti-rabbit IgG secondary antibody and detection with iFluor® 488 Styramide™ (Cat# 45020), iFluor® 555 Styramide™ (Cat# 45027), or iFluor® 647 Styramide™ (Cat# 45045). Fluorescence signal was evaluated before and after treatment with ReadiUse™ Tyramide (TSA)/Styramide (PSA) Stripping Buffer (Cat# 45091).

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 if you have any questions.