

ab326080 – Tyramide (TSA)/Styramide (PSA) Optimized Reaction Buffer

For research use only - not intended for diagnostic use.

Materials Supplied

Item	Quantity	Storage Condition
TSA/PSA Buffer	100 mL	4°C
Stabilized 3% Hydrogen peroxide	100 µL	4°C

Protocol summary

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

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

Styramide (PSA) stock solution (100X):

Add 100 µL of DMSO into a vial of iFluor® dye-labeled Styramide conjugate (for example: ab326081, ab326082, ab326083, ab326084 and ab326085) to make 100X Styramide stock solution.

Note: Make single-use aliquots, and store unused 100X stock solution at 2-8 °C in a dark place and avoid repeat freeze-thaw cycles.

Hydrogen peroxide stock solution:

Add 10 µL of 3% hydrogen peroxide to 90 µL of ddH₂O.

Note: Prepare the 100X H₂O₂ solution fresh on the day of use.

Preparation of Working Solutions:

Styramide (PSA) working solution (1X):

Every 1 mL of PSA working solution requires 1 mL TSA/PSA buffer, 10 µL of Styramide stock solution, and 10 µL of H₂O₂ stock solution.

Note: 1 mL is enough for 10 tests based on 100 µL of Styramide working solution needed per coverslip or per well in a 96-well microplate.

Note: The Styramide working solution must be used within 2 hours after preparation and avoid direct exposure to light.

Technical Support

For technical support contact information, visit:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)