

December 11, 2018



Product Information

Ready-to-Use Tyramide Amplification Buffer, 1X

Catalog Number: 22027-25mL; 22027-100mL

Unit Size

22027-25mL: 25 mL 22027-100mL: 100 mL

Storage and Handling

Store at 4°C. Protect from light. Product is stable for at least 6 months from date of receipt when stored as recommended.

Product Description

The 1X Ready-to-use Amplification Buffer is designed for cell and tissue immunostaining with tyramide signal amplification. Simply dilute the tyramide compounds in this buffer to make a working solution; no additional H_2O_2 is needed. Although specifically optimized for our CF® Dye-, Biotin XX-, and DNP-tyramide products, the 1X Ready-to-use Amplification Buffer is also also compatible with tyramide probes from other manufacturers.

Protocol

The following protocol is optimized for fixed cells in 96-well plates, or tissue sections with size about 10 mm X 10 mm. It can be scaled up and down based on the size of sample.

- Perform fixation, blocking, and antibody labeling using methods recommended for your samples and antibodies. The sample must be labeled with antibody or streptavidin-horseradish peroxidase (HRP) conjugates. We recommend to use our tyramide amplification kits (see Related Products) for optimal blocking and secondary antibody labeling.
- After labeling with antibody-HRP conjugates, wash samples 3 times with appropriate washing buffer, incubating 5 minutes at room temperature for each wash. Based on the nature of the sample and the antibody, phosphate buffered saline (PBS) or Tris buffered saline (TBS), with or without 0.1% TWEEN® 20 can be used as washing buffer.
- 3. Make a working solution by diluting dye-tyramide probe in the 1X Ready-touse Amplification Buffer. The final concentration of dye-tyramide probe may vary depending on the dye and the application. For our CF®dye-, Biotin XX-, and DNP-tyramide, we recommend a final concentration of 2 uM. However, a titration from 0.5 uM to 10 uM may be necessary to optimize the labeling results. Make 100 uL of working solution for each well on 96-well plate or each tissue section. The working solution can be stored at room temperature, protecting from light, for up to 24 hours.
- Apply 100 uL of working solution to each sample. Incubate at room temperature for 10 minutes. Protect from light.
- 5. Wash samples 3 times with washing buffer (see Step 2), incubating 5 minutes at room temperature for each wash.
- The samples are now ready for fluorescence imaging. We recommend to cover the well or mount the coverslip with our EverBrite[™] Mounting Medium (see Related Products).

Related Products

Catalog No.	Product
33000; 33003; 33006; 33009; 33012; 33015; 33018	Tyramide Amplification Kits with HRP Goat Anti-Mouse IgG and CF® Dye/BiotinTyramide
33001; 33004; 33007; 33010; 33013; 33016; 33019	Tyramide Amplification Kits with HRP Goat Anti-Rabbit IgG and CF® Dye/BiotinTyramide
33002; 33005; 33008; 33011; 33014; 33017; 33020	Tyramide Amplification Kits with HRP Streptavidin and CF® Dye/Biotin-Tyramide
20401	HRP Goat Anti-Mouse IgG (H+L), Highly Cross-Adsorbed
20403	HRP Goat Anti-Rabbit IgG (H+L), Highly Cross-Adsorbed
20406	HRP Goat Anti-Rat IgG (H+L), Highly Cross-Adsorbed
20474	HRP Goat Anti-Chicken IgY (H+L), Highly Cross-Adsorbed
20470	HRP Goat Anti-Human IgG (H+L), Highly Cross-Adsorbed
20475	HRP Goat Anti-Llama IgG (H+L), Highly Cross-Adsorbed
92021; 92022; 92170 - 92175; 92194 - 92198	CF® Dye Tyramides
22011	Fish Gelatin Powder
22014	BSA, IgG- and protease-free, 30% Solution
22002	TWEEN® 20
23001	EverBrite™ Mounting Medium
23002	EverBrite [™] Mounting Medium with DAPI
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite [™] Hardset Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
23007	TrueBlack® Lipofuscin Autofluorescence Quencher
23008	Drop-n-Stain EverBrite™ Mounting Medium
23009	Drop-n-Stain EverBrite™ Mounting Medium with DAPI
90082	DMSO, Anhydrous
22020	10X Phosphate-Buffered Saline (PBS)
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative

Please visit our website at www.biotium.com for information on our life science research products, including fluorescent CF® Dye antibody conjugates and reactive dyes, near-infrared CF® Dye conjugates for western blotting and other applications, EvaGreen® dye and master mixes for qPCR, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.

TWEEN is a registered trademark of Uniqema Americas LLC.