Ready-to-Use Tyramide Amplification Buffer, 1X

For simple, optimized tyramide amplification of immunostaining in cells or tissue sections. No need to add hydrogen peroxide before use.

The state of the s

Product attributes

Call us: 800-304-5357

Storage Conditions

Store at 2 to 8 °C, Protect from light

Email: btinfo@biotium.com

Product Description

Ready-to-Use Tyramide Amplification Buffer was designed for simple, optimized tyramide amplification of immunostaining in cells or tissue sections. There is no need to add hydrogen peroxide before use.

Features:

- Tyramide amplification buffer
- Ready-to-use
- No need to add hydrogen peroxide
- Simply dilute tyramide probe in buffer and proceed to signal development
- Compatible with tyramides from Biotium or other suppliers

Please note that Biotium has released a new, improved buffer for tyramide signal amplification. The Tyramide Amplification Buffer Plus (22029) provides significantly improved brightness, specificity, and sensitivity over this original buffer.

Ready-to-Use Tyramide Amplification Buffer was designed for simple, optimized signal amplification of immunostaining in cells or tissue sections. Simply mix your tyramide compound with the buffer to make a working solution for signal amplification. Although specifically optimized for our CF® Dye-, Biotin XX-, and DNP-tyramide products, Ready-to-Use Tyramide Amplification Buffer is also compatible with tyramide probes from other manufacturers.

To learn more about tyramide signal amplification, including our Tyramide Amplification Kits with a selection of bright and photostable CF® dyes, see the <u>Tyramide Amplification Technology Page</u>.

This datasheet was generated on November 7, 2025 at 11:23:22 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/ready-to-use-tyramide-amplification-buffer-1x/