



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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.

Product attributes

Storage Conditions
Store at 2 to 8 °C, Protect from light

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.

Also learn about [TyraMax™ Amplification Dyes and Kits](#), Biotium's next generation tyramide dyes that offer brighter signal compared to the original [CF® Dye Tyramides](#), and have advantages in brightness, photostability, and working solution stability compared to other TSA dyes. We also offer [CF® Dye Tyramide Amplification Kits](#). Learn more about [Tyramide Signal Amplification](#).

This datasheet was generated on June 15, 2026 at 01:08:00 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/ready-to-use-tyramide-amplification-buffer-1x/>