



Glowing products for science

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

## Tyramide Amplification Buffer Plus

An optimized buffer with enhanced sensitivity for tyramide signal amplification of immunostaining in cells or tissue sections.



### Product attributes

Storage Conditions

Store at 2 to 8 °C, Protect from light

## Product Description

Tyramide Amplification Buffer Plus is an optimized buffer with enhanced sensitivity for tyramide signal amplification of immunostaining in cells or tissue sections. It provides significantly improved brightness, specificity, and sensitivity compared to our original Ready-to-Use Tyramide Amplification Buffer ([22027](#)).

Tyramide Amplification Buffer Plus was developed as a new and improved version of our original Ready-to-Use Tyramide Amplification Buffer ([22027](#)). It is also a component in all of our Tyramide Amplification Kits. To learn more about tyramide signal amplification, including our Tyramide Amplification Kits with a selection of bright and photostable CF® dyes, see the [Tyramide Amplification Technology Page](#).

This datasheet was generated on June 18, 2026 at 10:55:49 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/tyramide-amplification-buffer-plus/>