



Glowing products for science

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

Tyramide Amplification Kit with HRP Streptavidin and CF® Dye or Biotin Tyramide



Kits for increasing immunofluorescence sensitivity using HRP-catalyzed tyramide signal amplification; with your choice of next-generation fluorescent CF® dyes or biotin.

Product Description

This kit contains reagents for performing tyramide signal amplification (TSA) using streptavidin HRP conjugate and fluorescent CF® Dyes or biotin to increase immunofluorescence signal.

Tyramide amplification, sometimes called Catalyzed Reporter Deposition (CARD), is an enzyme-mediated detection method that uses the catalytic activity of horseradish peroxidase (HRP) to generate high density labeling of a target protein or nucleic acid sequence *in situ*. The tyramide amplification method has been reported to increase the sensitivity by up to 100-fold compared to conventional avidin-biotin detection. Multiple rounds of tyramide signal amplification can be performed for multiplex detection; to learn more, see our [Tech Tip: Multicolor Fluorescence Imaging using Tyramide Amplification Kits](#).

These kits include your choice of 6 CF® dye tyramides or biotin tyramide. CF® dyes are Biotium's line of next-generation fluorescent dyes, with advantages in brightness and photostability compared to other dyes. Learn more about [CF® dyes](#).

Also see our [Tyramide Amplification Kits with HRP Goat Anti-Mouse](#) and [Tyramide Amplification Kits with HRP Goat Anti-Rabbit](#).

Each kit contains reagents for staining 50-150 slides; the actual number of samples that can be stained will depend on staining volumes used. [Tyramide Amplification Buffer](#), [CF® Dye Tyramides](#), and other tyramide compounds are also available separately.

Also learn about [TyrMax™ 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 [Ready-to-Use Tyramide Amplification Buffer](#) and [Tyramide Amplification Buffer Plus](#) (an improved formulation for enhanced TSA sensitivity). Learn more about [Tyramide Signal Amplification](#).

Product attributes

Storage Conditions	Store at -10 to -35 °C, Protect dye component from light

Tyramide Signal Amplification Kits

Tyramide Label	Ex/Em	Secondary conjugate	Catalog no.
CF@488A	490/515 nm	Goat anti-mouse HRP	33000
Goat anti-rabbit HRP	33001		
Streptavidin HRP	33002		
CF@543	541/560 nm	Goat anti-mouse HRP	33003
Goat anti-rabbit HRP	33004		
Streptavidin HRP	33005		
CF@568	562/583 nm	Goat anti-mouse HRP	33006
Goat anti-rabbit HRP	33007		
Streptavidin HRP	33008		
CF@594	593/614 nm	Goat anti-mouse HRP	33009
Goat anti-rabbit HRP	33010		
Streptavidin HRP	33011		
CF@640R	642/662 nm	Goat anti-mouse HRP	33012
Goat anti-rabbit HRP	33013		
Streptavidin HRP	33014		
CF@680R	680/701 nm	Goat anti-mouse HRP	33015
Goat anti-rabbit HRP	33016		
Streptavidin HRP	33017		
Biotin-XX	N/A	Goat anti-mouse HRP	33018
Goat anti-rabbit HRP	33019		
Streptavidin HRP	33020		

*Note: Older versions of this kit contain Ready-to-Use Tyramide Amplification Buffer (component 22027-25mL) without a separate vial of hydrogen peroxide. Download the product information sheet for detailed protocols. The SDS for the older version of the kit can be downloaded here: [SDS-Tyramide-Amplification-Kit with 22027](#).
[Back to top](#)

This datasheet was generated on January 14, 2026 at 03:07:11 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/tyramide-amplification-kit-with-hrp-streptavidin-and-cfbiotin-tyramide/>

