Glowing products for science

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

Bietun

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 <a href="Tech Tip: Multicolor Fluorescence Imaging using Tyramide Amplification Kits">Tech Tip: Multicolor Fluorescence Imaging using Tyramide Amplification Kits</a>.

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. <u>Tyramide Amplification Buffer</u>, <u>CF® Dye Tyramides</u>, and other tyramide compounds are also available separately.

Also learn about TyraMaxTM 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

Call us: 800-304-5357

Storage Conditions

Store at -10 to -35  $^{\circ}\text{C},$  Protect dye component from light

Email: btinfo@biotium.com

## **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	<u>33001</u>
		Streptavidin HRP	33002
CF®543	541/560 nm	Goat anti-mouse HRP	33003
		Goat anti-rabbit HRP	<u>33004</u>
		Streptavidin HRP	<u>33005</u>
CF®568	562/583 nm	Goat anti-mouse HRP	<u>33006</u>
		Goat anti-rabbit HRP	<u>33007</u>
		Streptavidin HRP	33008
CF®594	593/614 nm	Goat anti-mouse HRP	<u>33009</u>
		Goat anti-rabbit HRP	<u>33010</u>
		Streptavidin HRP	<u>33011</u>
CF®640R	642/662 nm	Goat anti-mouse HRP	<u>33012</u>
		Goat anti-rabbit HRP	<u>33013</u>
		Streptavidin HRP	<u>33014</u>
CF®680R	680/701 nm	Goat anti-mouse HRP	<u>33015</u>
		Goat anti-rabbit HRP	<u>33016</u>
		Streptavidin HRP	<u>33017</u>
Biotin-XX	N/A	Goat anti-mouse HRP	<u>33018</u>
		Goat anti-rabbit HRP	<u>33019</u>
		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 November 15, 2025 at 04:15:13 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/tyramide-amplification-kit-with-hrp-streptavidin-and-cfbiotin-tyramide/">https://biotium.com/product/tyramide-amplification-kit-with-hrp-streptavidin-and-cfbiotin-tyramide/</a>