Tyramide Amplification Kit with HRP Goat Anti-Mouse and CF® Dye or Biotin Tyramide



Product attributes

Cross adsorption Bovine, Horse, Human, Rabbit, Swine

Storage Conditions Store at -10 to -35 °C. Protect dve component from light

Call us: 800-304-5357 Email: btinfo@biotium.com

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 goat anti-mouse HRP secondary antibody 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. Learn more about tyramide signal amplification.

The HRP goat anti-mouse antibody provided in the kits has been adsorbed against bovine, horse, human, rabbit, and swine serum to minimize cross-reactivity; the antibody is not cross-adsorbed against rat. The 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 <u>Tyramide Amplification Kits with HRP Goat Anti-Rabbit and Tyramide Amplification Kits with HRP Streptavidin.</u>

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. See our full selection of tyramide products.

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	<u>33003</u>
		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	<u>33008</u>
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	<u>33020</u>

*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