

Cyanine 555 Tyramide

Cyanine 555 tyramide is used for tyramide signal amplification (TSA), an horseradish peroxidase (HRP)-catalyzed high-density staining method for increasing immunofluorescence sensitivity.



Product Description

Cyanine 555 tyramide conjugate is used for TSA in multicolor immunocytochemistry (ICC), immunohistochemistry (IHC), or in situ hybridization (ISH). The HRP-catalyzed method enables high-density labeling of a target protein or nucleic acid with Cyanine 555. Multiple tyramide conjugates are deposited onto tyrosines adjacent to the target for enhanced fluorescent signal.

- Orange-red fluorescent TSA reagent
- High-density labeling of a target protein or nucleic acid for enhanced immunofluorescence sensitivity
- Especially suited for the detection of low abundance targets
- Detection sensitivity of over 100-fold compared to conventional procedures
- Enables multiplex multicolor detection

We also offer CF®555 tyramide, a superior alternative to Cyanine 555 with better photostability and brightness.

See also our other TSA reagents including <u>CF®</u> <u>dye tyramides</u>, <u>Ready-to-Use Tyramide Amplification Buffer, Tyramide Amplification Buffer Plus</u> (an improved formulation for enhanced TSA sensitivity), and <u>CF®</u> <u>Dye Tyramide Amplification Kits</u>.

Learn more about Tyramide Signal Amplification.

Also see CF® Dye Tyramides, available with a wide selection of our bright and photostable CF® Dyes.

This datasheet was generated on November 22, 2025 at 11:38:09 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cyanine-555-tyramide/

Product attributes

Call us: 800-304-5357

Excitation/Emission	555/565 nm
Chemical reactivity (reacts with)	Tyrosine residues
Functional group	Tyramide
Assay type/options	Tissue staining
Storage Conditions	Store at -10 to -35 °C, Protect from light
Detection method/readout	Fluorescence microscopy
Molecular weight	~851 g/mol