



Glowing products for science

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

## 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 [CF@ dye tyramides](#), [Ready-to-Use Tyramide Amplification Buffer](#), [Tyramide Amplification Buffer Plus](#) (an improved formulation for enhanced TSA sensitivity), and [CF@ Dye Tyramide Amplification Kits](#).

Learn more about [Tyramide Signal Amplification](#).

Also see [CF@ Dye Tyramides](#), available with a wide selection of our bright and photostable CF@ Dyes.

### Product attributes

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

This datasheet was generated on January 21, 2026 at 11:40:17 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cyanine-555-tyramide/>