

XFD532 tyramide

Catalog Number: 11072

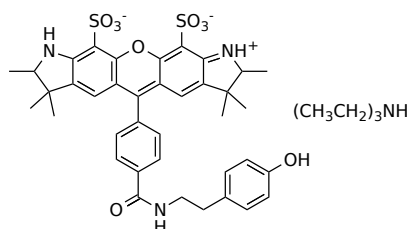
Unit Size: 200 Slides

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid red
Molecular Weight	847.06
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	534 nm
Emission Wavelength	553 nm

Applications

Tyramide signal amplification (TSA) has proven to be a particularly versatile and powerful enzyme amplification technique with improved assay sensitivity. TSA is based on the ability of HRP, in the presence of low concentrations of hydrogen peroxide, to convert labeled tyramine-containing substrate into an oxidized, highly reactive free radical that can covalently bind to tyrosine residues at or near the HRP. The signal amplification conferred by the turnover of multiple tyramide substrates per peroxidase label translates ultrasensitive detection of low-abundance targets and the use of smaller amounts of antibodies and hybridization probes. In immunohistochemical applications, sensitivity enhancements derived from TSA method allow primary antibody dilutions to be increased to reduce nonspecific background signals, and can overcome weak immunolabeling caused by suboptimal fixation procedures or low levels of target expression. XFD 532 tyramide contains the Alexa Fluor® 532 fluorophore that can be readily detected with the Alexa Fluor® 532/ATTO 532 filter set (Alexa Fluor® is the trademark of ThermoFisher). XFD 532 tyramide has intense yellow-orange fluorescence color.