

Bodi Fluor™ 576/589 acid (equivalent to BODIPY™ 576/589)

Catalog number: 660 Unit size: 10 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Solid
Molecular Weight	329.11
Soluble In	DMSO

Applications

Bodi Fluor™ 576/589 is chemically the same molecule to BODIPY™ 576/589 dye. It is a bright red fluorescent dye with high cell-permeability. It has a high extinction coefficient and fluorescence quantum yield. Its fluorescence is relatively insensitive to solvent polarity and pH change. In contrast to the highly water-soluble fluorophores such as iFluor™ and Alexa Fluor® 488 dyes, Bodi Fluor dyes have unique hydrophobic properties ideal for staining lipids, membranes, and other lipophilic compounds. Bodi Fluor 576/589 dye has a relatively long excited-state lifetime (typically 5 nanoseconds or longer), which is useful for fluorescence polarization-based assays and a large two-photon cross-section for multiphoton excitation.