

AATOM™ 565 Acid, 5-Isomer

Catalog Number: 2826

Unit Size: 25 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	624.61
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	562 nm
Emission Wavelength	589 nm

Applications

AATOM™ 565 Acid, 5-Isomer is manufactured by AAT Bioquest for research and development use. AATOM™ 565 Acid, 5-Isomer is an amine-reactive rhodamine dye commonly used for labeling peptides, proteins, and other amine-containing molecules, such as amino-modified oligonucleotides. In the presence of a coupling reagent such as EDC, this dye reacts with amine groups to form a stable amide bond. AATOM™ 565 Acid, 5-Isomer demonstrates strong absorption, high fluorescence quantum yield, and good thermal and photostability. However, it has low water solubility. For applications requiring high water solubility, iFluor® 568 acid is recommended as an alternative. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.