

AATOM™ 565 NHS ester

Catalog Number: 2828

Unit Size: 5 mg

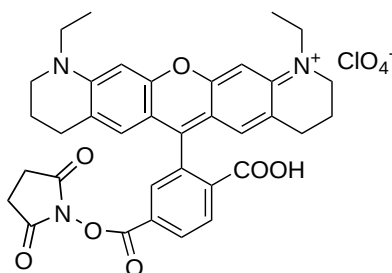
Product Details

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

Chemical Properties

Appearance	Solid purple
Molecular Weight	708.12
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	562 nm
Emission Wavelength	589 nm

Applications

AATOM™ 565 NHS ester is a commonly used amine-reactive rhodamine dye. It might be used for labeling peptides, proteins and other amino-containing compounds such as amino-modified oligos. It readily reacts with compounds containing amino groups, forming a chemically stable amide bond between the dye and amino compounds. The optimum pH range for NHS-ester coupling is pH 8.0 – 9.0. At this pH, amino groups of proteins (such as the ε-amino groups of lysines) are unprotonated enough for fast coupling. AATOM™ 565 NHS ester has strong absorption, high fluorescence quantum yield and high thermal/photostability. However, it has low water solubility. For the applications that require high water solubility, you may consider iFluor® 568 succinimidyl ester that has almost identical fluorescence properties with greatly enhanced water solubility. In addition, the protein conjugates made from iFluor® 568 succinimidyl ester are much brighter than those from AATOM™ 565 NHS ester. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.