

AATOM™ 565 maleimide

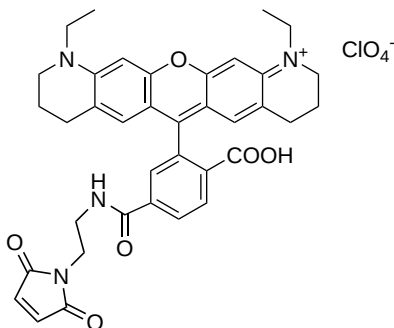
Catalog Number: 2829

Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid purple
Molecular Weight	733.17
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	562 nm
Emission Wavelength	589 nm

Applications

AATOM™ 565 Maleimide is a commonly used thiol-reactive rhodamine dye. It might be used for labeling peptides, proteins and other thiol-containing compounds such as thiol-modified oligos. It readily reacts with compounds containing thiol groups, forming a chemically stable thiol ether bond between the dye and thiol compounds. The optimum pH range for maleimide coupling is pH 5.5 – 7.0. AATOM™ 565 Maleimide 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 Maleimide that has almost identical fluorescence properties with greatly enhanced water solubility. In addition, the protein conjugates made from iFluor® 568 Maleimide are much brighter than those from AATOM™ 565 Maleimide. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.