

AATOM™ Rho12 NHS ester

Catalog Number: 70371

Unit Size: 1 mg

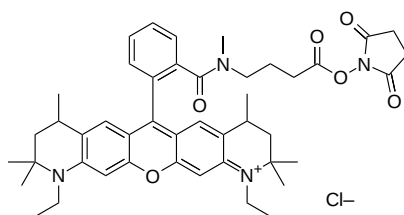
Product Details

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

Chemical Properties

Appearance	Solid purple
Molecular Weight	783.40
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	575 nm
Emission Wavelength	600 nm

Applications

AATOM™ Rho12 is a rhodamine-based fluorescent dye characterized by its bright fluorescence, good photostability, and strong absorption properties. It exhibits excitation and emission maxima of 575/600 nm and can be efficiently excited using commonly available yellow-green excitation sources. These properties make AATOM™ Rho12 suitable for fluorescence imaging, biomolecule labeling, flow cytometry, and other fluorescence-based biological applications.

AATOM™ Rho12 NHS ester reacts selectively and efficiently with primary amines (such as the side chains of lysine residues or aminosilane-coated surfaces) at pH 7-9, forming stable covalent amide bonds. This property makes AATOM™ Rho12 NHS ester suitable for labeling peptides, proteins, antibodies, and amine-modified oligonucleotides. The resulting conjugates can be used in fluorescence microscopy, immunocytochemistry (ICC), flow cytometry, high-content screening (HCS), and related fluorescence-based applications. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.