

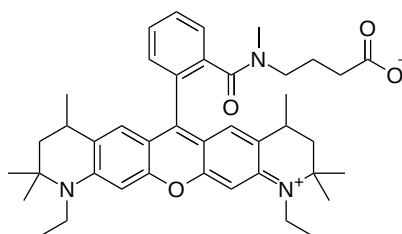
Unit Size: 5 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	649.86
Soluble In	DMSO



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 Acid is a non-reactive compound that can be employed as a reference standard in studies utilizing AATOM™ Rho12 conjugates. It is also suitable for use as a control in confocal microscopy, immunocytochemistry (ICC), high-content screening (HCS), and flow cytometry applications. Furthermore, it can be utilized in the synthesis of activated esters and can be activated for conjugation to amino-containing molecules using water-soluble carbodiimides (e.g., EDAC). This enables labeling of proteins, antibodies, amine-modified oligonucleotides, peptides, and other biomolecules. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.