

AATOM™ Rho12 Azide

Catalog Number: 70373

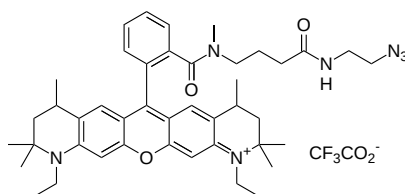
Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	831.97
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, super-resolution microscopy, and other fluorescence-based biological applications.

The azide derivative of AATOM™ Rho12 is designed for click chemistry labeling of terminal alkyne-containing biomolecules. It participates in copper-catalyzed azide-alkyne cycloaddition (CuAAC) reactions with alkyne-functionalized molecules and strain-promoted azide-alkyne cycloaddition (SPAAC) reactions with DBCO- or BCN-containing compounds. The resulting conjugates are suitable for fluorescence microscopy, flow cytometry, immunocytochemistry (ICC), high-content screening (HCS), and super-resolution imaging applications including PALM, dSTORM, and STED microscopy. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.