

AATOM™ 532 PEG3 azide

Catalog Number: 2824

Unit Size: 1 mg

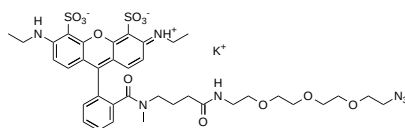
Product Details

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

Chemical Properties

Appearance	Solid red
Molecular Weight	884.03
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	531 nm
Emission Wavelength	552 nm

Applications

AATOM™ 532 is a rhodamine-based fluorescent dye notable for its high molar absorptivity and fluorescence quantum yield (0.90), providing robust signal intensity in fluorescence applications. Its combination of photostability, aqueous solubility, and sufficient Stokes shift makes it suitable for single-molecule detection and high-resolution microscopy techniques, including SIM and STED microscopy. AATOM™ 532 is also effective in flow cytometry, FISH, and various biological assays, offering flexibility for diverse fluorescence-based experimental protocols. The dye is optimally excited within the 515-545 nm range, with a frequency-doubled Nd:YAG laser at 532 nm serving as an ideal excitation source.

The azide derivative of AATOM™ 532 is widely used for labeling terminal alkynes on peptides, antibodies, and other biomolecules via click chemistry. It participates in copper-catalyzed azide-alkyne cycloaddition (CuAAC) with alkyne-containing molecules and strain-promoted alkyne-azide cycloaddition (SPAAC) with DBCO- or BCN-containing molecules. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.