

AATOM™ 514 PEG4 DBCO

Catalog Number: 2847

Unit Size: 1 mg

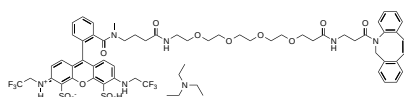
Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	1360.44
Soluble In	DMSO

Chemical Structure

**Spectral Properties**

Excitation Wavelength	510 nm
Emission Wavelength	531 nm

Applications

AATOM™ 514 is a hydrophilic fluorescent dye characterized by excellent aqueous solubility, strong molar absorptivity, high fluorescence quantum yield, and notable thermal and photostability. These properties make it well-suited for advanced imaging techniques, including single-molecule detection and super-resolution microscopy methods such as PALM, dSTORM, and STED. AATOM™ 514 is also compatible with flow cytometry (FACS), fluorescence in situ hybridization (FISH), and other fluorescence-based assays. The dye is optimally excited within the 510–535 nm range, with the 514 nm line of an Argon-Ion laser providing efficient excitation. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.