

AATOM™ 514 azide

Catalog Number: 2845

Unit Size: 1 mg

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

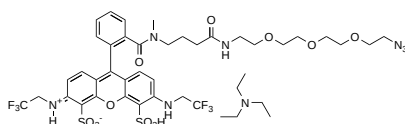
Chemical Properties

Appearance Solid

Molecular Weight 1055.07

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.