

AATOM™ 565 BCN

Catalog Number: 70559

Unit Size: 1 mg

Product Details

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

Expiration Date

Chemical Properties

Appearance Solid

Molecular Weight N/A

Soluble In DMSO

Spectral Properties

Excitation Wavelength 562 nm

Emission Wavelength 589 nm

Applications

AATOM™ 565 BCN is a clickable derivative of AATOM™ 565, an orange fluorescent dye designed for labeling peptides, oligonucleotides, and other biomolecules. AATOM™ 565 is known for its strong absorption, high fluorescence quantum yield, and excellent thermal and photo-stability. This dye exhibits moderate hydrophilicity, with an optimal excitation range of 545-575 nm. AATOM™ 565 is well-suited for advanced applications in single-molecule detection and high-resolution microscopy techniques, including PALM, dSTORM, and STED microscopy. It is also compatible with flow cytometry (FACS), fluorescence in situ hybridization (FISH), and a variety of other biological assays.

To improve conjugation performance, AATOM™ 565 BCN incorporates a PEG spacer, which reduces steric hindrance and minimizes potential interference with target binding sites. This design maximizes conjugation efficiency while preserving the biological activity of the resulting conjugate. The bicyclononyne (BCN) moiety enables strain-promoted azide-alkyne cycloaddition (SPAAC) with azido groups, forming stable triazole linkages under catalyst-free conditions. In addition, unlike dibenzocyclooctyne (DBCO), BCN also reacts efficiently with tetrazines through an inverse electron-demand Diels-Alder (IEDDA) reaction. This reaction is rapid, selective, and bioorthogonal, allowing labeling of biomolecules under physiological conditions without the need for metal catalysts or disruption of native biological processes.

This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.