

AATOM™ 532 BCN

Catalog Number: 70557

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 N/A

Soluble In DMSO

Spectral Properties

Excitation Wavelength 531 nm

Emission Wavelength 552 nm

Applications

AATOM™ 532 BCN is a clickable derivative of AATOM™ 532, a green fluorescent dye designed for labeling peptides, oligonucleotides, and other biomolecules. AATOM™ 532 is known for its strong absorption, high fluorescence quantum yield, and excellent water solubility. It is optimally excited within the 515-545 nm range, with the 532 nm output of a frequency-doubled Nd laser serving as an ideal excitation source.

To improve conjugation performance, AATOM™ 532 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.