

AATOM™ 610 Tetrazine

Catalog Number: 70257

Unit Size: 1 mg

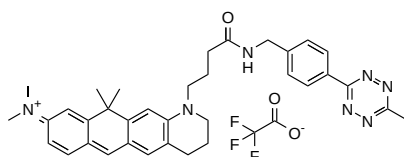
Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	687.77
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	615 nm
Emission Wavelength	632 nm

Applications

AATOM™ 610 is a carbopyronin-based fluorescent dye known for its strong absorption, high fluorescence quantum yield, and exceptional photostability and thermal stability. It is moderately hydrophilic and optimally excited at wavelengths between 595 and 625 nm. Upon coupling to a substrate, AATOM™ 610 becomes cationic, carrying a net electrical charge of +1. The dye remains stable under physiological pH conditions and in buffers with a pH of up to 8, though it gradually degrades at higher pH levels. AATOM™ 610 is ideal 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), FRET, and various other biological assays.

AATOM™ 610 tetrazine is particularly useful for labeling TCO-modified biomolecules under copper-free conditions. It reacts with TCO-functionalized molecules, forming a stable conjugate via a dihydropyrazine moiety. This click reaction is favored over others due to its extremely fast kinetics and higher yields under mild reaction conditions, making it a popular choice for researchers. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.