

# AATOM™ 594 TCO

Catalog Number: 2843

Unit Size: 1 mg

## Product Details

---

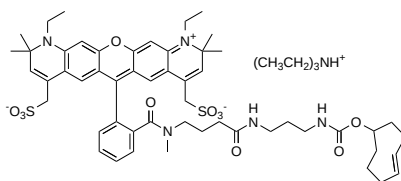
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

## Chemical Properties

---

Appearance	Solid purple
Molecular Weight	1115.46
Soluble In	DMSO

Chemical Structure



## Spectral Properties

---

Excitation Wavelength	602 nm
Emission Wavelength	621 nm

## Applications

---

AATOM™ 594 is a bright, red fluorescent dye characterized by a strong absorption, high fluorescence quantum yield, and exceptional thermal and photostability. The dye exhibits superior water solubility and hydrophilicity, facilitating its use in various aqueous environments. AATOM™ 594 is optimally excited within the 560-615 nm range, making it compatible with both 561 nm and 594 nm laser lines commonly used in advanced fluorescence imaging systems. Upon conjugation to biomolecules, AATOM™ 594 becomes anionic, carrying a net charge of -1, which may influence its binding characteristics and performance in assays. Its photostability and brightness make it particularly suited for high-resolution techniques like single-molecule detection and super-resolution microscopy, including PALM, dSTORM, and STED. Additionally, AATOM™ 594 is highly compatible with flow cytometry (FACS), fluorescence in situ hybridization (FISH), and various other fluorescence-based assays, supporting its broad utility in complex biological studies.

AATOM™ 594 TCO is particularly useful for labeling tetrazine-modified biomolecules under copper-free conditions. It reacts with tetrazine-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.