

# AATOM™ 565 acid

Catalog Number: 2827

Unit Size: 25 mg

## Product Details

---

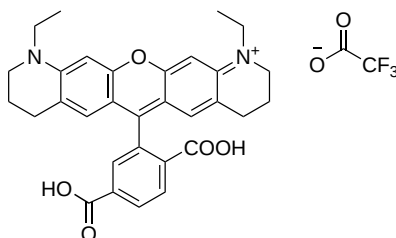
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

## Chemical Properties

---

Appearance	Solid
Molecular Weight	624.61
Soluble In	DMSO

Chemical Structure



## Spectral Properties

---

Excitation Wavelength	562 nm
Emission Wavelength	589 nm

## Applications

---

AATOM™ 565 acid is a commonly used amine-reactive rhodamine dye. It might be used for labeling peptides, proteins and other amino-containing compounds such as amino-modified oligos. It reacts with compounds containing amino groups in the presence of a coupling reagent such as EDC. It forms a chemically stable amide bond between the dye and amino compounds. AATOM™ 565 acid has strong absorption, high fluorescence quantum yield and high thermal/photostability. However, it has low water solubility. For the applications that require high water solubility, you may consider iFluor® 568 acid. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.