

# AATOM™ 514 maleimide

Catalog Number: 2864

Unit Size: 1 mg

## Product Details

---

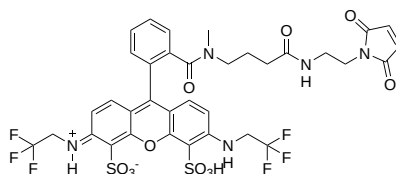
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

## Chemical Properties

---

Appearance	Solid red
Molecular Weight	875.77
Soluble In	DMSO

Chemical Structure



## Spectral Properties

---

Excitation Wavelength	510 nm
Emission Wavelength	531 nm

## Applications

---

AATOM™ 514 is a rhodamine-based fluorescent labeling dye known for its favorable spectral qualities. With a high molar extinction coefficient, a fluorescence quantum yield of 0.85, high photo-stability, and excellent water solubility, it is an ideal choice for single-molecule detection, high-resolution microscopy (e.g., SIM and STED), flow cytometry, and fluorescence in-situ hybridization (FISH). It exhibits its highest fluorescence efficiency when excited within the range of 510-535 nm, making it a perfect match for the 514 nm line of an argon-ion laser. AATOM™ 514 maleimide is used for labeling thiol-containing molecules, such as proteins, antibodies, ligands, and oligonucleotide thiophosphates. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.