

## ATTO 532 maleimide

Catalog Number: 2823

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	532 nm
Emission Wavelength	552 nm

### Applications

---

ATTO 532 is a rhodamine-based fluorescent labeling dye known for its favorable spectral qualities. With a high molar extinction coefficient, a fluorescence quantum yield of 0.90, 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 is specifically optimized for excitation using a frequency-doubled Nd:YAG-Laser. ATTO 532 maleimide is used for labeling thiol-containing molecules, such as proteins, antibodies, ligands, and oligonucleotide thiophosphates.