

# AATOM™ 514 NHS ester

Catalog Number: 2863

Unit Size: 1 mg

## Product Details

---

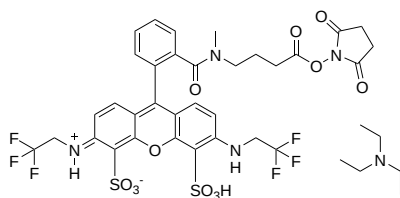
|                    |  |
|--------------------|--|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
| Expiration Date    | 12 months upon receiving                   |

## Chemical Properties

---

|                  |           |
|------------------|-----------|
| Appearance       | Solid red |
| Molecular Weight | 951.91    |
| Soluble In       | DMSO      |

Chemical Structure



## Spectral Properties

---

|                       |        |
|-----------------------|--------|
| Excitation Wavelength | 510 nm |
| Emission Wavelength   | 531 nm |

## Applications

---

AATOM™ 514, a hydrophilic rhodamine-based fluorophore, is known for its excellent water solubility, strong absorption, high fluorescence quantum yield, and thermal and photo-stability. It is well-suited for single-molecule detection and high-resolution microscopy techniques like PALM, dSTORM, and STED. Additionally, it finds applications in flow cytometry (FACS), fluorescence in-situ hybridization (FISH), and other scientific methodologies. AATOM™ 514 exhibits optimal fluorescence efficiency when excited within the 510-535 nm range, making it a match for the 514 nm line of an Argon-Ion laser. AATOM™ 514 NHS ester is a popular tool used to label the primary amines of proteins, peptides, and amino-modified oligonucleotides. This product is manufactured by AAT Bioquest and is not affiliated with ATTO-TEC GmbH.