

5-TAMRA ethylenediamine

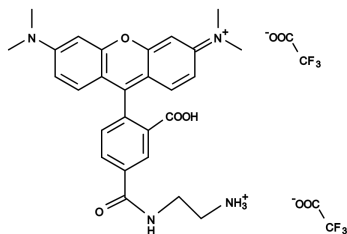
Catalog number: 358

Unit size: 5 mg

Product Details

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

Chemical Properties

| | |
|--------------------|--|
| Appearance | Red solid |
| Molecular Weight | 700.58 |
| Soluble In | DMSO |
| Chemical Structure |  |

Spectral Properties

| | |
|-----------------------|--------|
| Excitation Wavelength | 552 nm |
| Emission Wavelength | 578 nm |

Applications

5-TAMRA ethylenediamine is a building block for developing TAMRA bioconjugates. TAMRA is one of the most popular bright orange fluorophores used in various bioconjugations.