

## AzoDye-3 Amine

Catalog number: 2476  
Unit size: 5 mg

### Product Details

---

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

### Chemical Properties

---

Appearance	Solid
Molecular Weight	816.81
Soluble In	DMSO

### Applications

---

AzoDye-3 amine is an excellent building block for labeling carbonyl-containing molecules. AzoDye-3 dyes are nearly non-fluorescent with moderate extinction coefficients. These dark quenchers are most typically used in Molecular Beacon probes in which FRET has minimal contribution to their fluorescence quenching, thus dependence on donor-acceptor spectral overlap is not significant as to other FRET probes such as FRET peptides. AzoDye-3 dyes can quench a variety of fluorophores, in particular, fluorescein derivatives such as FAM, HEX, TET, NED and JOE. This may reflect the AzoDye-3's ability to form stable, non-fluorescent complexes with a fluorophore.