

**AzoDye-3 Azide**Catalog number: 2482  
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	714.75
Soluble In	DMSO

**Applications**

---

AzoDye-3 azide is an excellent building block for labeling alkyne-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, Cy5 derivatives such as Cy5, Cy5.5, DyLight 647, Alexa Fluor® 647 and iFluor® 647. This may reflect the AzoDye-3's ability to form stable, non-fluorescent complexes with a fluorophore.