

**AzoDye-3 PEG10 Alkyne**Catalog number: 2485  
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 | 1138.29 |
| Soluble In       | DMSO    |

**Applications**

---

AzoDye-3 PEG10 alkyne is an excellent building block for labeling azide-containing molecules. It contains a PEG10 spacer that significantly improves its water solubility compared to AzoDye-3 alkyne (#2484) although both AzoDye-3 alkyne (#2484) and AzoDye-3 PEG10 alkyne (#2485) bear the same extremely hydrophobic BHQ-2 chromophore. 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.