

**AzoDye-3 CPG (500 A)**

Catalog number: 2478

Unit size: 100 mg

**Product Details**

---

Storage Conditions Freeze (&lt;math&gt;-15\text{ }^{\circ}\text{C}&lt;/math&gt;), Minimize light exposure

Expiration Date 12 months upon receiving

**Chemical Properties**

---

Appearance Solid

Soluble In MeCN

**Applications**

---

AzoDye-3 CPG is an excellent building block for preparing FRET oligo 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.