

AzoDye-2 CPG (500 A)

Catalog number: 2428

Unit size: 100 mg

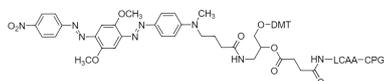
Product Details

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

Chemical Properties

| | |
|------------------|-------------------|
| Appearance | Dark purple Solid |
| Molecular Weight | N/A |
| Soluble In | MeCN |

Chemical Structure

**Applications**

AzoDye-2 CPG is an excellent building block for preparing FRET oligo molecules. AzoDye-2 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-2 dyes can quench a variety of fluorophores, in particular, fluorescein derivatives such as FAM, HEX, TET, NED and JOE. This may reflect the AzoDye-2's ability to form stable, non-fluorescent complexes with a fluorophore.