

AzoDye-1 CPG (1000 A)

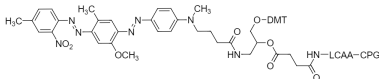
Catalog number: 2410

Unit size: 100 mg

Product Details

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

Chemical Properties

Appearance	Dark brown Solid
Molecular Weight	N/A
Soluble In	MeCN
Chemical Structure	

Applications

AzoDye-1 CPG is an excellent building block for preparing FRET oligo molecules. AzoDye-1 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-1 dyes can quench a variety of fluorophores, in particular, fluorescein derivatives such as FAM, HEX, TET, NED and JOE. This may reflect the AzoDye-1's ability to form stable, non-fluorescent complexes with a fluorophore.