

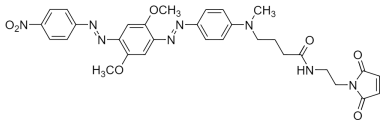
AzoDye-2 Maleimide

Catalog number: 2424
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	628.65
Soluble In	DMSO
Chemical Structure	

Applications

AzoDye-2 maleimide is an excellent building block for labeling thiol-containing 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.