

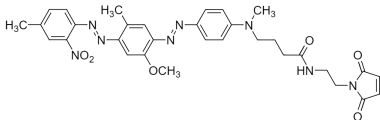
## AzoDye-1 Maleimide

Catalog number: 2404  
Unit size: 1 mg

### Product Details

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

### Chemical Properties

Appearance	Dark red Solid
Molecular Weight	626.67
Soluble In	DMSO
Chemical Structure	

### Applications

AzoDye-1 maleimide is an excellent building block for labeling thiol-containing 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.