

AzoDye-3 Carboxylic Acid

Catalog number: 2470

Unit size: 25 mg

Product Details

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Solid

Molecular Weight 660.70

Soluble In DMSO

Applications

AzoDye-3 Carboxylic Acid is an excellent building block for labeling amino-containing 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, Cy5 derivatives such as Cy5, Cy5.5, DyLight 647, Alexa Fluor® 647 and iFluor® 647. This may reflect the AzoDye-3's ability to form stable, non-fluorescent complexes with a fluorophore.