

AzoDye-2 Carboxylic Acid

Catalog number: 2420

Unit size: 25 mg

Product Details

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

Expiration Date 12 months upon receiving

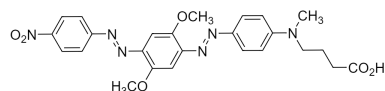
Chemical Properties

Appearance Solid

Molecular Weight 506.52

Soluble In DMSO

Chemical Structure



Applications

AzoDye-2 Carboxylic Acid is an excellent building block for labeling amino-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.