

## AzoDye-2 Succinimidyl Ester

Catalog number: 2422  
Unit size: 5 mg

### Product Details

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

### Chemical Properties

|                    |   |
|--------------------|---|
| Appearance         | Dark purple Solid   |
| Molecular Weight   | 603.59  |
| Soluble In         | DMSO  |
| Chemical Structure |  |

### Applications

AzoDye-2 Succinimidyl Ester 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.