

BXQ-1 Succinimidyl EsterCatalog number: 2402
Unit size: 5 mg**Product Details**

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

Chemical Properties

Appearance	Solid
Molecular Weight	601.62
Soluble In	DMSO

Applications

Black Hole Quencher™ dyes (BHQ®) are non-fluorescent dyes mainly used for preparing FRET oligonucleotides (Black Hole Quencher™ and BHQ® are the trademarks of Biosearch). They 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. BHQ dyes can quench a variety of fluorophores, in particular, fluorescein derivatives such as FAM, HEX, TET, NED and JOE. This may reflect the BHQ's ability to form stable, non-fluorescent complexes with a fluorophore. BXQ-1 is chemically equivalent to BHQ®-1. BXQ-1 succinimidyl ester is an excellent building block for labeling amino-modified oligonucleotides.