

DABCYL C2 maleimide

Catalog Number: 2008

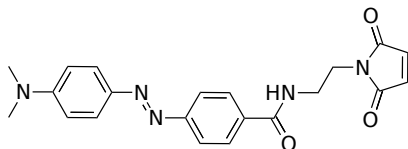
Unit Size: 25 mg

Product Details

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

Chemical Properties

Appearance	Solid orange
Molecular Weight	391.42
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

DABCYL C2 maleimide is the thiol-reactive form of DABCYL, and widely used to prepare a variety of FRET-based probes that contain DABCYL. DABCYL is one of the most popular acceptors for developing FRET-based nucleic acid probes and protease substrates. It is often paired with EDANS or a blue fluorescent donor such as coumarin dyes.