

DABCYL succinimidyl ester [4-((4-(Dimethylamino)phenyl)azo)benzoic acid, succinimidyl ester] *CAS 146998-31-4*

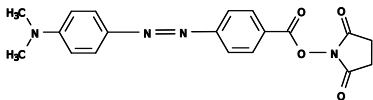
Catalog Number: 2004, 2005

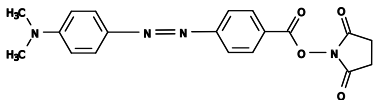
Unit Size: 1 g, 5 g

Product Details

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

Chemical Properties

Appearance	Solid red
Molecular Weight	366.37
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

DABCYL, SE is the amino-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.