

Product Information Sheet

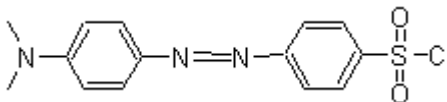
Ordering Information

Product Number: 2030
Product Name: DABSYL chloride [4-Dimethylaminoazobenzene-4-sulfonyl chloride]
Unit Size: 1 g
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 323.8

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 430 nm
Emission Wavelength: NA

Application Notes

DABSYL is widely used for labeling amino acids, also used as an acceptor for developing FRET-based nucleic acid probes and protease substrates (often paired with EDANS).