

## DABCYL acid [4-((4-(Dimethylamino)phenyl)azo)benzoic acid] \*CAS 6268-49-1\*

Catalog Number: 2001

Unit Size: 5 g

### Product Details

---

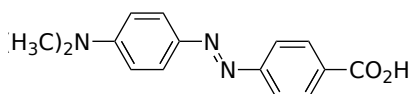
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Solid red
Molecular Weight	269.3
Soluble In	DMSO

Chemical Structure



### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

DABCYL acid is one of the most popular acceptors for developing FRET-based nucleic acid probes such as Molecular Beacon-based PCR probes. It is also used to prepare a variety of FRET peptides that are often used for the detection of various proteases. It is predominantly paired with EDANS (as fluorescence donor) in FRET-based fluorescent probes.