

## DABCYL C2 amine

Catalog Number: 2006

Unit Size: 100 mg

### Product Details

---

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

### Chemical Properties

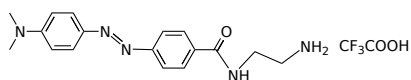
---

Appearance Solid red

Molecular Weight 425.41

Soluble In DMSO

Chemical Structure



### Spectral Properties

---

Excitation Wavelength N/A

Emission Wavelength N/A

### Applications

---

DABCYL C2 amine is the carbonyl-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.