

## BAPTA Amine

Catalog Number: 20420

Unit Size: 5 mg

### Product Details

---

Storage Conditions Freeze (&lt; -15 °C), Minimize light exposure

Expiration Date 24 months upon receiving

### Chemical Properties

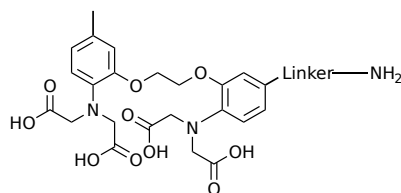
---

Appearance Solid

Molecular Weight 633.53

Soluble In DMSO

Chemical Structure



### Spectral Properties

---

Excitation Wavelength N/A

Emission Wavelength N/A

### Applications

---

BAPTA Amine is an excellent building block to introduce BAPTA calcium chelator into a biological molecule that contains a carbonyl group such as COOH or CHO. BAPTA (1,2-bis(o-aminophenoxy)ethane-N,N,N',N'-tetraacetic acid) is a common calcium-specific chelator. The presence of four carboxylic acid functional groups makes BAPTA possibly binding two calcium ions. The extensive flexibility of the carboxylate ligands is critical to the coordination of calcium and other metal ions. It is commonly used to chelate Ca<sup>2+</sup>, similarly to EGTA.