

BAPTA Alkyne

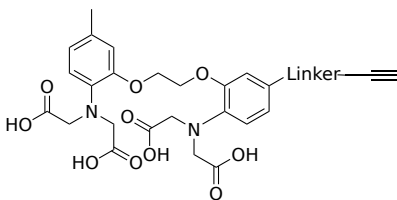
Catalog Number: 20424

Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	571.54
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

BAPTA Alkyne is an excellent building block to introduce BAPTA calcium chelator into a biological molecule that contains an azido group such as azido-modified antibodies, proteins, peptides and oligos via the well-known click chemistry. BAPTA (1,2-bis(o-aminophenoxy)ethane-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. BAPTA is commonly used to chelate Ca²⁺, similarly to EGTA.