

# BAPTA, tetrapotassium salt \*CAS 73630-08-7\*

Catalog Number: 21003

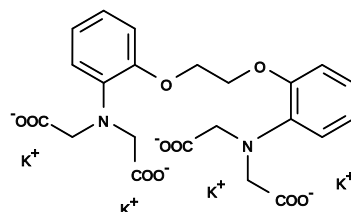
Unit Size: 100 mg

## Product Details

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

## Chemical Properties

Appearance	Solid off-white
Molecular Weight	628.79
Soluble In	Water
Chemical Structure	



## Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

## Applications

BAPTA (1,2-bis(o-aminophenoxy)ethane-N,N,N',N'-tetraacetic acid) is a water-soluble and cell-impermeable calcium chelator. It has high selectivity to calcium. BAPTA is a calcium-specific aminopolycarboxylic acid. The presence of four carboxylic acid functional groups makes possible the binding of two calcium ions. The extensive flexibility of the carboxylate ligands is critical to the coordination of calcium and other metal ions.