

BAPTA, free acid

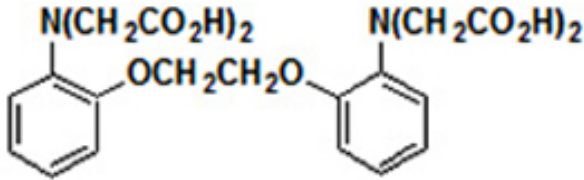
Catalog Number	Packaging Size
C200	100 mg

Storage upon receipt: Room temperature

Introduction

BAPTA, free acid is a cell-impermeant chelator, which is a highly selective for Ca^{2+} over Mg^{2+} , and it can be used to control the level of extracellular Ca^{2+} . BAPTA is more selective for Ca^{2+} than EDTA and EGTA, and its metal binding is also less sensitive to pH.

Specifications

Molecular Formula:	$\text{C}_{22}\text{H}_{24}\text{N}_2\text{O}_{10}$	
Molecular Weight:	476.44	
CAS Number:	85233-19-8	
Product Size:	100 mg	
Storage Conditions:	Room Temperature	
Shipping Condition:	Room Temperature	

Applications

Calcium chelator

References:

[A practical guide to the preparation of \$\text{Ca}^{2+}\$ buffers.](#)

Bers DM, Patton CW, Nuccitelli R
Methods Cell Biol (1994) 40:3-29

[A store-operated calcium channel in *Drosophila* S2 cells.](#)

Yeromin AV, Roos J, Stauderman KA, Cahalan MD
J Gen Physiol (2004) 123:167-182

[Role of the inositol 1,4,5-trisphosphate receptor in \$\text{Ca}^{2+}\$ feedback inhibition of calcium release-activated calcium current \(I_{crac}\).](#)

Broad LM, Armstrong DL, Putney JW
J Biol Chem (1999) 274:32881-32888