

BAPTA, AM ester

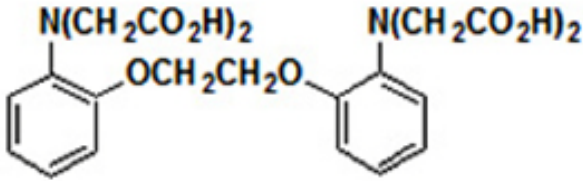
Catalog Number	Packaging Size
C203	25 mg

Storage upon receipt: -20°C

Introduction

BAPTA, AM ester is a cell-permeant chelator, which is highly selective for Ca^{2+} over Mg^{2+} , and it can be used to control the level of intracellular 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:	$\text{C}_{34}\text{H}_{40}\text{N}_2\text{O}_{18}$	
Molecular Weight:	764.69	
CAS Number:	126150-97-8	
Product Size:	100 mg	
Storage Conditions:	-20°C	
Shipping Condition:	Room Temperature	

Applications

Calcium chelator

References:

[Tonotopic variations of calcium signalling in turtle auditory hair cells.](#)

Ricci AJ, Gray-Keller M, Fettiplace R
J Physiol (2000) 524 Pt 2:423-436

[Cadherins mediate intercellular mechanical signaling in fibroblasts by activation of stretch-sensitive calcium-permeable channels.](#)

Ko KS, Arora PD, McCulloch CA
J Biol Chem (2001) 276:35967-35977

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

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