

5,5'-Dimethyl BAPTA, AM Ester

BAPTA and its derivatives are calcium chelators that are commonly used to form calcium buffers with well-defined calcium concentrations. By injecting the chelators into cells or by incubating cells with the AM ester form of the chelators, one can control the cytosolic calcium concentration, an important means to study the roles of calcium.



Product attributes

CAS number	147504-94-7
Cell permeability	Membrane permeant

Product Description

BAPTA and its derivatives are calcium chelators that are commonly used to form calcium buffers with well-defined calcium concentrations. By injecting the chelators into cells or by incubating cells with the AM ester form of the chelators, one can control the cytosolic calcium concentration, an important means to study the roles of calcium. Key advantages of these calcium chelators include relative insensitivity toward intracellular pH change and fast release of calcium. Biotium offers several BAPTA chelators with calcium dissociation constants covering the biologically significant range from 10^{-7} to 10^{-2} M.

Dimethyl BAPTA AM ester is a membrane permeable form of Dimethyl BAPTA that can be loaded into cells by incubation. Dimethyl BAPTA AM ester itself does not bind calcium, but once inside the cell is converted into Dimethyl BAPTA by cytoplasmic esterases. See catalog no. [50008](#) for information on Dimethyl BAPTA.

- White to off-white solid soluble in DMSO
- Store desiccated at -20°C
- $\text{C}_{36}\text{H}_{44}\text{N}_2\text{O}_{18}$
- MW: 792.75
- [147504-94-7]

