

5,5'-Dimethyl BAPTA, Tetrapotassium Salt



Product attributes

CAS number	73630-09-8
Cell permeability	Membrane impermeant

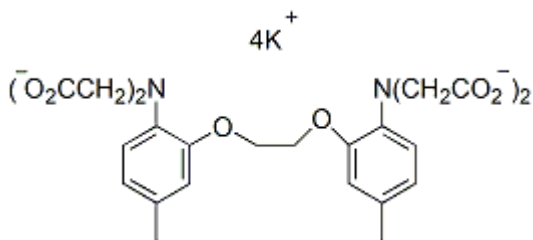
BAPTA and its derivatives are calcium chelators that are commonly used to form calcium buffers with well-defined calcium concentrations. By injecting the chelators into cells or by incubating cells with the AM ester form of the chelators, one can control the cytosolic calcium concentration, an important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

Product 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.

- Kd (no Mg^{2+}): 0.16 μ M, Kd (1 mM Mg^{2+}): 0.44 μ M
- White solid soluble in water
- Store at $-20^{\circ}C$
- $C_{24}H_{24}K_4N_2O_{10}$
- MW: 657
- [73630-09-8]



This datasheet was generated on June 15, 2026 at 12:48:25 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/55-dimethyl-baptatetrapotassium-salt/>