BAPTA, Tetrapotassium Salt

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 73630-08-7

Cell permeability Membrane impermeant

Call us: 800-304-5357 Email: btinfo@biotium.com

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 Mg²): 0.59 uM, Kd (1 mM Mg²): 0.70 uM
- White solid soluble in water
- Store at 4°C
- C₂₃H₂₄K₄N₂O₁₀
- MW: 656.85
- [73630-08-7]

This datasheet was generated on November 9, 2025 at 08:19:38 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/baptatetrapotassium-salt/