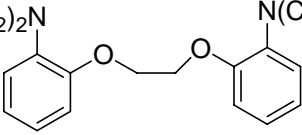


# Bapta (tetracesium salt)

O,O'-Bis(2-aminophenyl)ethyleneglycol-N,N,N',N'-tetraacetic acid, tetracesium salt, hydrate

## Instruction Manual

Catalog Number	PK-CA707-50001
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 towards intracellular pH change and fast release of calcium. PromoKine offers several BAPTA chelators with calcium dissociation constants covering the biologically significant range from $10^{-7}$ to $10^{-2}$ M (please request).
Quantity	1 g
Excitation / Emission Maxima	NA
Molecular Structure	<p style="text-align: center;">4 Cs<sup>+</sup></p> <p style="text-align: center;">(O<sub>2</sub>CCH<sub>2</sub>)<sub>2</sub>N      N(CH<sub>2</sub>CO<sub>2</sub>)<sub>2</sub></p> 
Molecular Weight / Molecular Formula	1001 Da; C <sub>23</sub> H <sub>24</sub> Cs <sub>4</sub> N <sub>2</sub> O <sub>10</sub>
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	White solid; soluble in water.
Storage & Stability	Store at 4°C.
Applications	Calcium chelator; for use in calcium assays
References	NA
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.