

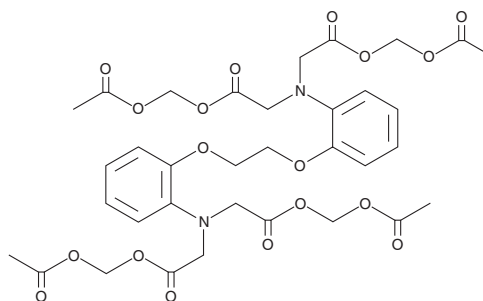
# PRODUCT INFORMATION



## BAPTA AM

Item No. 15551

**CAS Registry No.:** 126150-97-8  
**Formal Name:** N,N'-[1,2-ethanediylbis(oxy-2,1-phenylene)]  
bis[N-[2-[(acetyloxy)methoxy]-2-oxoethyl]-  
1,1'-bis[(acetyloxy)methyl] ester-glycine  
**Synonym:** BAPTA Acetoxymethyl ester  
**MF:** C<sub>34</sub>H<sub>40</sub>N<sub>2</sub>O<sub>18</sub>  
**FW:** 764.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 213, 248, 286 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

BAPTA AM is supplied as a crystalline solid. A stock solution may be made by dissolving the BAPTA AM in the solvent of choice. BAPTA AM is soluble in organic solvents such as DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of BAPTA AM in these solvents is approximately 20 mg/ml.

BAPTA AM is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BAPTA AM should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BAPTA AM has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

BAPTA (Item No. 11706) is membrane-impermeable calcium chelator that binds extracellular calcium ions ( $K_d = 0.11 \mu\text{M}$ ), with selectivity over magnesium ions or protons.<sup>1,2</sup> BAPTA and its derivatives can also be used as calcium indicators, since the absorption maximum for BAPTA changes when it is complexed with calcium (absorption maxima free/complexed = 254/274 nm, emission maxima free/complexed = 363/363 nm).<sup>3</sup> BAPTA AM is a cell permeable analog of BAPTA that binds calcium only after the acetoxymethyl group is removed by cytoplasmic esterases.<sup>4</sup> It is commonly used at 10-100  $\mu\text{M}$  to evaluate the role of intracellular calcium in cell signaling.<sup>4-7</sup> It can also be used in animals.<sup>4</sup> BAPTA AM also inhibits voltage-gated potassium ( $K_v$ ) channels, including  $K_v1.3$ ,  $K_v1.5$ , and  $K_v11.1$  ( $K_i = 1.45$ , 1.23, and 1.30  $\mu\text{M}$ , respectively).<sup>8</sup>

### References

1. Sabnis, R.W. Handbook of biological dyes and stains: Synthesis and industrial applications. John Wiley & Sons, Inc., Hoboken, NJ, USA (2010).
2. Gryniewicz, G., Poenie, M., and Tsien, R.Y. *J. Biol. Chem.* **260**(6), 3440-3450 (1985).
3. Tsien, R.Y. *Annu. Rev. Biophys. Bioeng.* **12**, 91-116 (1983).
4. Tymianski, M., Spigelman, I., Zhang, L., et al. *J. Cereb. Blood Flow Metab.* **14**(6), 911-923 (1994).
5. Verrière, V., Higgins, G., Al-Alawi, M., et al. *PLoS One* **7**(5), e37746 (2012).
6. Chang, M.G., Chang, C.Y., de Lange, E., et al. *Biophys. J.* **102**(12), 2706-2714 (2012).
7. Bray, G.E., Ying, Z., Baillie, L.D., et al. *J. Neurosci.* **33**(19), 8202-8215 (2013).
8. Tang, Q., Jin, M.W., Xiang, J.Z., et al. *Biochem. Pharmacol.* **74**(11), 1596-1607 (2007).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/29/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM