PRODUCT INFORMATION



EGTA AM

Item No. 20401

CAS Registry No.:	99590-86-0	
Formal Name:	3,12-bis[2-[(acetyloxy)methoxy]-	
	2-oxoethyl]-6,9-dioxa-3,12-	
	diazatetradecanedioic acid,	0 0
	1,14-bis[(acetyloxy)methyl] ester	
Synonym:	EGTA Acetoxymethyl ester	
MF:	C ₂₆ H ₄₀ N ₂ O ₁₈	
FW:	668.6	
Purity:	≥90%	$\mathcal{A} \sim \mathcal{A}$
Supplied as:	A neat oil	0 0
Storage:	-20°C	
Stability:	≥4 years	
Special Conditions: Protect from light		
1 6 1		

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

Laboratory Procedures

EGTA AM is supplied as a neat oil. A stock solution may be made by dissolving the EGTA AM in the solvent of choice, which should be purged with an inert gas. EGTA AM is soluble in DMSO.

Description

EGTA AM is a cell-permeant acetoxymethyl ester of the calcium chelator EGTA. As EGTA AM enters cells, it is hydrolyzed by intracellular esterases to produce EGTA. EGTA AM is used to evaluate the role of intracellular calcium changes in cell signaling.¹⁻³

References

- 1. Drumm, B. T., Large, R. J., Hollywood, M. A., et al. The role of Ca²⁺ influx in spontaneous Ca²⁺ wave propagation in interstitial cells of Cajal from the rabbit urethra. J. Physiol. 593(15), 3333-3350 (2015).
- 2. Jalini, S., Ye, H., Tonkikh, A. A., et al. Raised intracellular calcium contributes to ischemia-induced depression of evoked synaptic transmission. PLoS One 11(3), (2016).
- 3. Nanou, E., Sullivan, J. M., Scheuer, T., et al. Calcium sensor regulation of the CaV2.1 Ca²⁺ channel contributes to short-term synaptic plasticity in hippocampal neurons. Proc. Nat. Acad. Sci. USA 113(4), 1062-1067 (2016).

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.

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, 12/19/2022

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