

PRODUCT INFORMATION



Quin-2 AM

Item No. 20422

CAS Registry No.: 83104-85-2

Formal Name: N-[2-[(acetyloxy)methoxy]-2-oxoethyl]-
N-[2-[[8-[bis[2-[(acetyloxy)methoxy]-
2-oxoethyl]amino]-6-methoxy-2-
quinolinyl]methoxy]-4-methylphenyl]-
glycine(acetyloxy)methyl ester

Synonym:

Quin-2 Acetoxymethyl ester

MF:

C₃₈H₄₃N₃O₁₈

FW:

829.8

Purity:

≥90%

Supplied as:

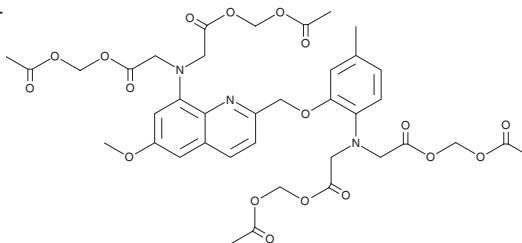
A solid

Storage:

-20°C

Stability:

As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

Quin-2 AM is supplied as a solid. A stock solution may be made by dissolving the quin-2 AM in the solvent of choice. Quin-2 AM is soluble in DMSO, which should be purged with an inert gas. We do not recommend storing the aqueous solution for more than one day.

Description

Quin-2 AM is a cell-permeant acetoxymethyl ester of the high affinity fluorescent calcium indicator quin-2 (Item No. 20421). As quin-2 AM enters cells, it is hydrolyzed by intracellular esterases to produce quin-2. Quin-2 binds calcium with a K_d value of 115 nM.¹ It displays high selectivity for calcium, as it is not affected by sodium gradients, membrane potential, or intracellular pH.¹ High affinity probes like quin-2 are ideal for monitoring low levels of calcium, as are found in resting cells.^{1,2} Loadings of up to 2 mM quin-2 are without serious toxic effects, so quin-2 may be used to buffer intracellular calcium transients. Excitation and emission maxima for quin-2 are 339 and 492 nm, respectively.

References

1. Tsien, R. Y., Pozzan, T., and Rink, T. J. Calcium homeostasis in intact lymphocytes: Cytoplasmic free calcium monitored with a new, intracellularly trapped fluorescent indicator *J. Cell. Biol.* **947(2)**, 325-334, (1982).
2. Minta, A., Kao, J. P., and Tsien, R. Y. Fluorescent indicators for cytosolic calcium based on rhodamine and fluorescein chromophores. *J. Bio. Chem.* **264(14)**, 8171-8178 (1989).

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, 09/15/2016

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