

PRODUCT INFORMATION



Cal Green™ 1 AM

Item No. 20400

CAS Registry No.: 186501-28-0
Formal Name: N-[2-[(acetyloxy)methoxy]-2-oxoethyl]-N-[2-[2-[5-[[[3',6'-bis(acetyloxy)-2',7'-dichloro-3-oxospiro[isobenzofuran-1(3H),9'-[9H]xanthen]-5-yl]carbonyl]amino]-2-[bis[2-[(acetyloxy)methoxy]-2-oxoethyl]amino]phenoxy]ethoxy]phenyl]-glycine, (acetyloxy)methyl ester

Synonym: Calcium Green-1 AM

MF: C₅₉H₅₃Cl₂N₃O₂₆

FW: 1,291.0

Purity: ≥90%

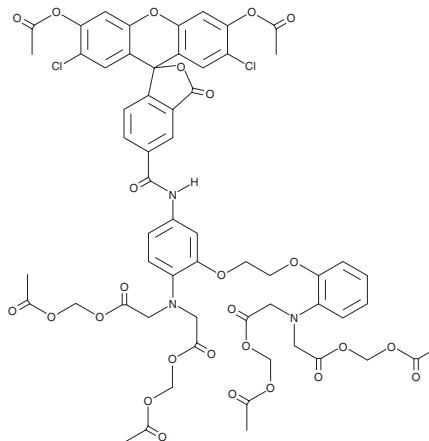
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: Protect from light

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



Laboratory Procedures

Cal Green™ 1 AM is supplied as a solid. A stock solution may be made by dissolving the Cal Green™ 1 AM in the solvent of choice, which should be purged with an inert gas. Cal Green™ 1 AM is soluble in the organic solvent DMSO.

Description

Cal Green™ 1 AM is a cell-permeant acetyloxy-methyl ester compound that is converted to the fluorescent calcium indicator Cal Green™ 1 (Item No. 20399) by intracellular esterases. This derivative is characterized by high quantum yield and low phototoxicity.¹ Its peak excitation and emission wavelengths (506 and 531 nm, respectively) are comparable to standard fluorescein dyes, making Cal Green™ 1 AM appropriate for fluorescent microscopy and flow cytometry. Cal Green™ 1 is ~5-fold brighter than fluo-3 (Item Nos. 20403 | 20404) at saturating calcium levels.¹

Reference

1. Paredes, R.M., Etzler, J.C., Watts, L.T., *et al.* Chemical calcium indicators. *Methods* **46(3)**, 143-151 (2008).

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, 12/01/2022

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