

PRODUCT INFORMATION



Indo-1 AM

Item No. 20417

CAS Registry No.: 112926-02-0
Formal Name: 2-[4-[bis[2-[(acetyloxy)methoxy]-2-oxoethyl]amino]-3-[2-[2-[bis[2-[(acetyloxy)methoxy]-2-oxoethyl]amino]-5-methylphenoxy]ethoxy]phenyl]-1H-Indole-6-carboxylic acid, (acetyloxy)methyl ester

Synonym: Indo-1 Acetoxymethyl ester

MF: C₄₇H₅₁N₃O₂₂

FW: 1,009.9

Purity: ≥90%

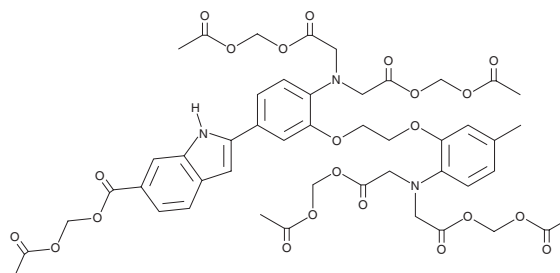
Supplied as: A solid

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

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

Description

Indo-1 AM is a cell-permeant acetoxymethyl ester of the ratiometric fluorescent calcium indicator indo-1 (Item Nos. 20418 | 20419). As indo-1 AM enters cells, it is hydrolyzed by intracellular esterases to produce indo-1. Indo-1 is ideal for analyses using flow cytometry, as it uses a single excitation source, typically 349-364 nm light from an argon-ion laser. The emission maximum shifts from 475-485 nm without calcium to 400-410 nm when indo-1 binds calcium.^{1,2} Indo-1 is prone to photobleaching, which limits its usefulness in methods involving microscopy.¹

References

1. Paredes, R.M., Etzler, J.C., Watts, L.T., *et al.* Chemical calcium indicators. *Methods* **46(3)**, 143-151 (2008).
2. Grynkiewicz, G., Poenie, M., and Tsien, R.Y. A new generation of Ca²⁺ indicators with greatly improved fluorescence properties. *J. Biol. Chem.* **260(6)**, 3440-3450 (1985).

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