

PRODUCT INFORMATION



BTC AM

Item No. 20424

CAS Registry No.: 176767-94-5
Formal Name: N-[2-[(acetyloxy)methoxy]-2-oxoethyl]-N-[3-(2-benzothiazolyl)-6-[2-[2-[bis[2-[(acetyloxy)methoxy]-2-oxoethyl]amino]-5-methylphenoxy]ethoxy]-2-oxo-2H-1-benzopyran-7-yl]-glycine(acetyloxy) methyl ester

Synonym: BTC Acetoxymethyl ester

MF: C₄₅H₄₅N₃O₂₀S

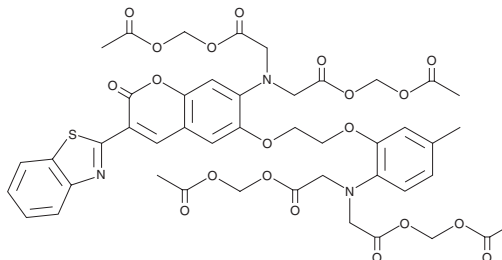
FW: 979.9

Purity: ≥90%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

BTC AM is supplied as a solid. A stock solution may be made by dissolving the BTC AM in the solvent of choice. BTC 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

BTC AM is a cell-permeable acetoxy-methyl ester of BTC (Item No. 20423). As BTC AM enters cells, it is hydrolyzed by intracellular esterases to produce BTC. BTC is a low affinity calcium indicator ($K_d = 7-26 \mu\text{M}$) that displays excitation/emission spectra of 401/529 nm, respectively. It exhibits a shift in the excitation maximum from approximately 480 to 401 nm upon calcium binding, enabling determination of ratiometric calcium measurements.¹ BTC is suitable for detecting elevated calcium levels associated with activation of smooth muscle, neurons, and intracellular calcium stores.^{1,2}

References

1. Sabnis, R. W. Handbook of biological dyes and stains: Synthesis and industrial applications. (2010).
2. Hyrc, K. L., Bownik, J. M., and Goldberg, M. P. Neuronal free calcium measurement using BTC/AM, a low affinity calcium indicator. *Cell Calcium* **24**(3), 165-175 (1998).

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, 09/22/2017

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