

PRODUCT INFORMATION



Fura-2 AM

Item No. 14591

CAS Registry No.: 108964-32-5
Formal Name: 2-[6-[bis[2-[(acetyloxy)methoxy]-2-oxoethyl]amino]-5-[2-[2-[bis[2-[(acetyloxy)methoxy]-2-oxoethyl]amino]-5-methylphenoxy]ethoxy]-2-benzofuranyl]-5-oxazolecarboxylic acid, (acetyloxy)methyl ester

Synonym: Fura-2 Acetoxymethyl ester

MF: C₄₄H₄₇N₃O₂₄

FW: 1,001.9

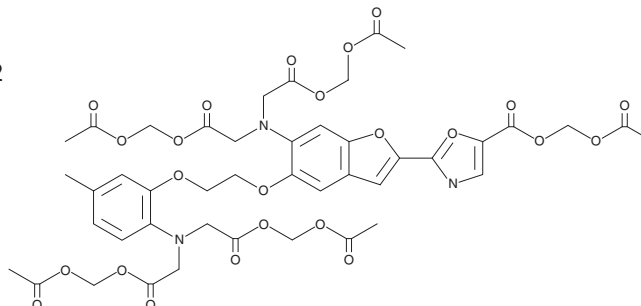
Purity: ≥95%

Ex./Em. Max: 340 and 380/510 nm

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

Fura-2 AM is supplied as a solid. A stock solution may be made by dissolving the fura-2 AM in the solvent of choice. Fura-2 AM is soluble in DMSO at a concentration of approximately 10 mM.

Description

Fura-2 AM is a membrane permeable, UV light-excitable form of Fura-2 (Item Nos. 19531 | 20414).¹ It crosses cell membranes and is converted to Fura-2 via cellular esterases.² Fura-2 is a ratiometric fluorescent calcium indicator that can be used to detect calcium in cells.³ It is a pentacarboxylate that displays excitation maxima of 340 and 380 nm at high and low calcium concentrations, respectively, when the emission is fixed at 510 nm, enabling determination of ratiometric measurements of calcium influx in live cells.

References

1. Sabnis, R.W. *Handbook of biological dyes and stains: Synthesis and industrial applications*. John Wiley & Sons, Inc., Hoboken, NJ, USA (2010).
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).
3. Kong, S.K. and Lee, C.Y. The use of fura 2 for measurement of free calcium concentration. *Biochem. Mol. Biol. Ed.* **23**(2), 97-98 (1995).

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, 05/03/2018

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