

PRODUCT INFORMATION



Rhod-FF AM

Item No. 20444

Formal Name: N-(9-(4-(bis(2-(acetoxymethoxy)-2-oxoethyl)amino)-3-(2-(6-(bis(2-(acetoxymethoxy)-2-oxoethyl)amino)-2,3-difluorophenoxy)ethoxy)phenyl)-6-(dimethylamino)-3H-xanthen-3-ylidene)-N-methylmethanaminium bromide

Synonym: Rhod-FF Acetoxymethyl ester

MF: C₅₁H₅₅F₂N₄O₁₉ • Br

FW: 1,145.9

Purity: ≥90%

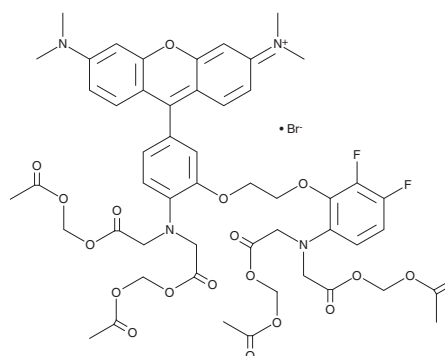
Supplied as: A solid

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

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

Description

Rhod-FF AM is a cell-permeant acetoxymethyl ester of the fluorescent calcium indicator rhod-FF (Item No. 20433). As rhod-FF AM enters cells, it is hydrolyzed by intracellular esterases to produce rhod-FF. Rhod-FF is a difluorinated analog of the cell-impermeant calcium indicator rhod-2. It has a very low affinity for calcium ($K_d = 320 \mu\text{M}$).¹ Low affinity calcium indicators are particularly useful for studying compartments with high concentrations of calcium, such as endoplasmic reticulum, where high affinity dyes will be insensitive to luminal fluctuations. Rhod-FF has excitation/emission maxima of approximately 552 and 580 nm, respectively.

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, 11/22/2022

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