

PRODUCT INFORMATION



Calcein Deep Red™ AM

Item No. 30575

Synonym:	Calcein Deep Red™ Acetoxymethyl ester
FW:	1,000.0
Purity:	≥90%
Ex./Em. Max:	643/663 nm
Supplied as:	A solid
Storage:	-20°C
Stability:	≥4 years

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

Laboratory Procedures

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

Description

Calcein Deep Red™ AM is a fluorogenic probe. As the acetoxymethyl ester, Calcein Deep Red™ AM is non-fluorescent but cell-permeant and is hydrolyzed to the fluorescent probe Calcein Deep Red™ (Item No. 20634) by intracellular esterases. Calcein Deep Red™ is retained by live cells and displays excitation/emission maxima of 643/663 nm, respectively.

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, 12/07/2022

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