

PRODUCT INFORMATION



Calcein Red™ (sodium salt)

Item No. 20633

FW: 825.4
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Light sensitive

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

Laboratory Procedures

Calcein Red™ (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the Calcein Red™ (sodium salt) in the solvent of choice. Calcein Red™ (sodium salt) is soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Calcein Red™ is a membrane-impermeant fluorescent dye. It is a red-emitting variant of the green fluorophore calcein (Item No. 16221), which has been entrapped in liposomes and similar vesicles for studying membrane characteristics.^{1,2} Calcein Red™ displays excitation/emission maxima of 560/574 nm, respectively.

References

1. Sine, J., Urban, C., Thayer, D., *et al.* Photo activation of HPPH encapsulated in "Pocket" liposomes triggers multiple drug release and tumor cell killing in mouse breast cancer xenografts. *Int. J. Nanomedicine* **10**, 125-145 (2014).
2. Sudji, I.R., Subburaj, Y., Frenkel, N., *et al.* Membrane disintegration caused by the steroid saponin digitonin is related to the presence of cholesterol. *Molecules* **20(11)**, 20146-20160 (2015).

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, 11/02/2022

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