

# PRODUCT INFORMATION



## Calcein Deep Red™

Item No. 20634

**FW:** 373.7  
**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 Deep Red™ is supplied as a solid. A stock solution may be made by dissolving the Calcein Deep Red™ in the solvent of choice, which should be purged with an inert gas. Calcein Deep Red™ is soluble in the organic solvent DMSO.

### Description

Calcein Deep Red™ is a membrane-impermeant fluorescent dye. It is a deep red-emitting variant of the green fluorophore calcein (Item No. 16221), which has been entrapped in liposomes and similar vesicles for studying membrane characteristics.<sup>1,2</sup> Calcein Deep Red™ displays excitation/emission maxima of 646/659 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, 12/01/2022

#### 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](http://WWW.CAYMANCHEM.COM)