

# PRODUCT INFORMATION



## Calcein Orange™ Diacetate

Item No. 20641

FW:	877.2
Purity:	≥90%
Ex./Em. Max:	525/550 nm
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 Orange™ Diacetate is supplied as a solid. A stock solution may be made by dissolving the Calcein Orange™ Diacetate in the solvent of choice, which should be purged with an inert gas. Calcein Orange™ Diacetate is soluble in the organic solvent DMSO.

### Description

Calcein Orange™ Diacetate is a fluorogenic dye that is used to assess cell viability. It is hydrolyzed to the fluorescent probe Calcein Orange™ (Item No. 20640). Calcein Orange™ is retained by living cells and displays excitation/emission maxima of 525/550 nm, respectively. Calcein Orange™ Diacetate is an orange-emitting variant of the green fluorophore Calcein AM (Item No. 14948), which has been used to evaluate cell viability and cytotoxicity.<sup>1,2</sup>

### References

1. Bratosin, D., Mitrofan, L., Pali, C., *et al.* Novel fluorescence assay using calcein-AM for the determination of human erythrocyte viability and aging. *Cytometry* **66(1)**, 78-84 (2005).
2. Lévesque, A., Paquet, A., and Pagé, M. Measurement of tumor necrosis factor activity by flow cytometry. *Cytometry* **20(2)**, 181-184 (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, 11/11/2022

#### 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