

PRODUCT INFORMATION



CytoCalcein™ Violet 450

Item No. 20694

FW: 600.0
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥6 months
Special Conditions: Light sensitive

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

Laboratory Procedures

CytoCalcein™ Violet 450 is supplied as a solid. A stock solution may be made by dissolving the CytoCalcein™ Violet 450 in the solvent of choice. CytoCalcein™ Violet 450 is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

CytoCalcein™ Violet 450 is a fluorogenic dye used to assess cell viability.^{1,2} It is a cell-permeable dye that, upon entering cells, is cleaved by intracellular esterases. CytoCalcein™ Violet 450 is retained by living cells and displays excitation/emission maxima of 405/450 nm, respectively.

References

1. Kramer, C.E.M., Wiechert, W., and Kohlheyer, D. Artificial fluorogenic substrates in microfluidic devices for bacterial diagnostics in biotechnology. *J. Flow Chem.* **6(1)**, 3-7 (2016).
2. Price, T.O., Sheibani, N., and Shah, G.N. Regulation of high glucose-induced apoptosis of brain pericytes by mitochondrial CA VA: A specific target for prevention of diabetic cerebrovascular pathology. *Biochem. Biophys. Acta.* **1863(4)**, 929-935 (2017).

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, 05/01/2017

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