

PRODUCT INFORMATION



Nuclear Violet™ LCS1

Item No. 25177

Ex./Em. Max: 401/459 nm
Purity: ≥90%
Supplied as: A solution in DMSO
Storage: -20°C
Stability: ≥4 years

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

Description

Nuclear Violet™ LCS1 is a DNA-selective and cell-permeant dye that is used to label DNA content in living cells. Upon binding to double-stranded DNA, Nuclear Violet™ LCS1 fluoresces, and this fluorescence can be measured using fluorescence microscopy, microplate fluorometry, or flow cytometry. Nuclear Violet™ LCS1 displays excitation/emission maxima of 401/459 nm, respectively, and is well suited for multicolor analysis of live cells.

Assay Protocol

1. Add Nuclear Violet™ LCS1 to suspension or adherent cells to a final concentration of 0.5 to 10 µM and stain the cells for 15 to 60 minutes.
2. Wash the cells with growth medium, Hank's with 20 mM HEPES buffer, or a buffer of choice.
3. Observe cells using fluorescence technique of choice.

Note 1: Aliquot and store unused Nuclear Violet™ LCS1 at less than -15°C. Nuclear Violet™ LCS1 is light sensitive. Light exposure and repeated freeze-thaw cycles should be avoided.

Note 2: The optimal working concentration is application specific. Staining conditions may be modified according to growth medium and/or cell density. The presence of residual detergent on glassware may also affect staining.

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, 01/15/2026

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