

PRODUCT INFORMATION



Nuclear Blue™ DCS1

Item No. 25170

Ex./Em. Max: 350/461 nm
Supplied as: A solution in DMSO
Storage: -20°C
Stability: ≥2 years

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

Description

Nuclear Blue™ DCS1 is a DNA-selective, cell-impermeant dye that is used to label DNA content in live or fixed necrotic and apoptotic cells. Upon binding to double-stranded DNA, Nuclear Blue™ DCS1 fluoresces, and this fluorescence can be measured using fluorescence microscopy, microplate fluorometry, or flow cytometry. Nuclear Blue™ DCS1 displays excitation/emission maxima of 350/461 nm, respectively.

Assay Protocol

1. Add Nuclear Blue™ DCS1 to suspension or adherent cells to a final concentration of 2 to 10 µM, and stain the cells for 15 to 60 minutes.

Optional: Wash cells in PBS prior to fluorescence analysis.

2. Observe cells using fluorescence technique of choice.

Note 1: Aliquot and store unused Nuclear Blue™ DCS1 at -20°C. Nuclear Blue™ DCS1 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, 09/30/2022

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