

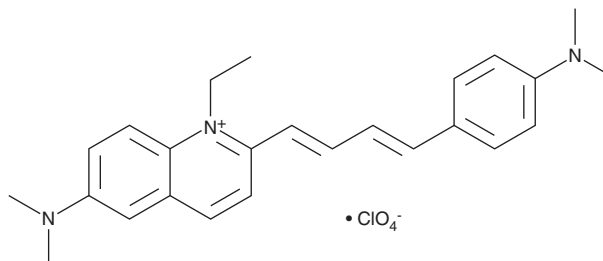
PRODUCT INFORMATION



LDS 751

Item No. 26859

CAS Registry No.: 181885-68-7
Formal Name: 6-(dimethylamino)-2-[4-[4-(dimethylamino)phenyl]-1,3-butadien-1-yl]-1-ethyl-quinolinium, monoperchlorate
MF: $C_{25}H_{30}N_3 \cdot ClO_4$
FW: 472.0
Ex./Em. Max: 543/712 nm
Supplied as: A powder
Storage: -20°C
Stability: ≥ 2 years



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

Description

LDS 751 is a cell-permeable fluorescent nucleic acid dye that can be used to label double-stranded DNA (dsDNA) within live or dead cells. Upon dsDNA binding, LDS 751 fluorescence is increased by approximately 20-fold, and this fluorescence can be measured using fluorescence microscopy, microplate fluorometry, or flow cytometry. LDS 751 displays excitation/emission maxima of 543/712 nm, respectively.

Assay Protocol

1. Prepare an LDS 751 stock solution by diluting to 5-10 mM in DMSO.
2. Add LDS 751 stock solution to suspension or adherent cell culture to a final concentration of 1 to 10 μ M and stain cells for 15 to 60 minutes.
3. Observe cells using fluorescence technique of choice.

Note: 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, 02/01/2019

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