

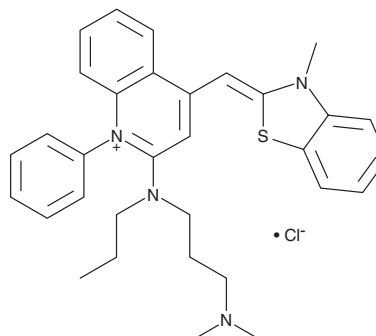
PRODUCT INFORMATION



Nucleic Acid Dye Green I

Item No. 35669

CAS Registry No.: 2225748-05-8
Formal Name: 2-[[3-(dimethylamino)propyl]propylamino]-4-[[3-methyl-2(3H)-benzothiazolylidene)methyl]-1-phenylquinolinium, monochloride
Synonyms: Green I Nucleic Acid Gel Stain, NucGel Green I Nucleic Acid Gel Stain
MF: C₃₂H₃₇N₄S • Cl
FW: 545.2
Purity: ≥95%
Ex./Em.: 497/520 nm
Supplied as: A solution in DMSO (10,000X)
Storage: -20°C
Stability: ≥4 years



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

Description

Nucleic Acid Dye Green I is a cell-impermeable fluorescent dye that has commonly been used for the quantification of dsDNA and ssDNA.¹ It displays excitation and emission maxima at 497 and 520 nm, respectively. Nucleic Acid Dye Green I has also been used in gene expression and RT-PCR assays.^{2,3}

References

1. Suzuki, T., Matsuzaki, T., Hagiwara, H., *et al.* Recent advances in fluorescent labeling techniques for fluorescence microscopy. *Acta Histochem. Cytochem.* **40(5)**, 131-137 (2007).
2. Ramos-Payán, R., Aguilar-Medina, M., Estrada-Parra, S., *et al.* Quantification of cytokine gene expression using an economical real-time polymerase chain reaction method based on SYBR® Green I. *Scand. J. Immunol.* **57(5)**, 439-445 (2003).
3. Ponchel, F., Toomes, C., Bransfield, K., *et al.* Real-time PCR based on SYBR-Green I fluorescence: An alternative to the TaqMan assay for a relative quantification of gene rearrangements, gene amplifications and micro gene deletions. *BMC Biotechnol.* **3**, 18 (2003).

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, 12/06/2023

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