

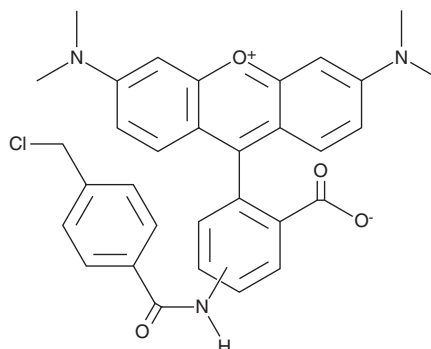
PRODUCT INFORMATION



CytoTrace™ Orange CMTMR

Item No. 20696

CAS Registry No.: 323192-14-9
Formal Name: 9-[2-carboxy-4(or 5)-[[4-(chloromethyl)benzoyl]amino]phenyl]-3,6-bis(dimethylamino)-xanthylium, inner salt
Synonym: CellTracker™ Orange CMTMR
MF: C₃₂H₂₈ClN₃O₄
FW: 554.0
Purity: ≥90%
Ex./Em. Max: 541/565 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Light sensitive



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

Laboratory Procedures

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

Description

CytoTrace™ Orange CMTMR is a rhodamine-based, cell-permeant, vital fluorescent dye that is retained by cells for prolonged periods (>72 hours) and is passed to daughter cells. Cells loaded with long-lasting dyes, including CMTMR, are used to follow cell trafficking and cell-to-cell interactions *in vivo*.¹⁻³ They have also been used to estimate the growth of implanted tumor spheroids.⁴ CytoTrace™ Orange CMTMR displays excitation/emission maxima of 541/565 nm, respectively.

References

1. Brooks, A.E.S., Parsamand, T., Kelly, R.W., *et al.* An improved quantitative method to assess adhesive properties of *Trichomonas vaginalis* to host vaginal ectocervical cells using flow cytometry. *J. Microbiol. Methods* **92**(1), 73-78 (2013).
2. Buehler, C., Dong, C.Y., So, P.T.C., *et al.* Time-resolved polarization imaging by pump-probe (stimulated emission) fluorescence microscopy. *Biophys J.* **79**(1), 536-549 (2000).
3. Sato, K., Imai, Y., and Irimura, T. Contribution of dermal macrophage trafficking in the sensitization phase of contact hypersensitivity. *J. Immunol.* **161**(12), 6835-6844 (1998).
4. Borgström, P., Hillan, K.J., Sriramarao, P., *et al.* Complete inhibition of angiogenesis and growth of microtumors by anti-vascular endothelial growth factor neutralizing antibody: Novel concepts of angiostatic therapy from intravital videomicroscopy. *Cancer Res.* **56**(17), 4032-4039 (1996).

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, 01/06/2023

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