

CELLULAR STAINING PROTOCOL

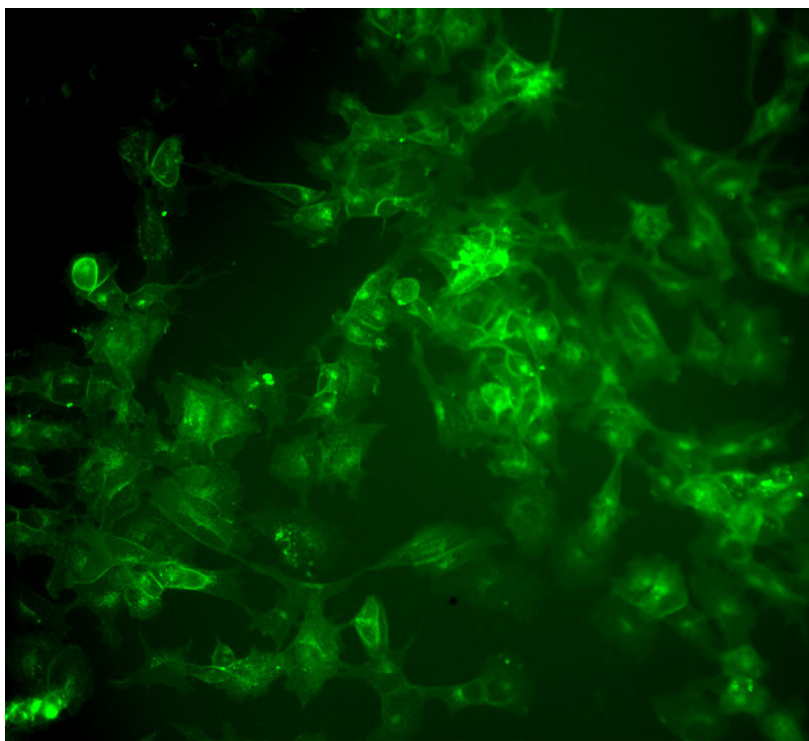


C3 BODIPY Ganglioside G_{M1} (d18:1/3:0) (ammonium salt) Item No. 33799

Staining SH-SY5Y Cells with C3-BODIPY-G_{M1}

1. Dissolve C3:0-BODIPY-G_{M1} (100 µg) in 64 µl DMSO:methanol:water (2:1:0.2) and aliquot. Store aliquots at -20°C until use.
2. Plate cells at density of 10,000/well in a 96-well plate in media with 10% FBS and allow to attach overnight.
3. The following day, remove media and replace with 2.5-10 µM C3:0-BODIPY-G_{M1} in phenol red-free media. Return cells to incubator for 30 minutes.
4. Remove media and replace with fresh phenol red-free media. Image using a fluorescence microscope with a FITC/GFP laser.

Note that complexing fluorescent lipids with bovine serum albumin (BSA) facilitates cell labeling by eliminating the need for organic solvents to dissolve the lipophilic probe. A BSA-complexed probe can be directly dissolved in water. However, gangliosides are unique lipids that are soluble in aqueous systems.



SH-SY5Y neuroblastoma cells stained with 10 µM C3-BODIPY G_{M1}.

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, 07/30/2021

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

TECHSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM