

PRODUCT INFORMATION



Methyltransferase Resorufin Buffer (10X)

Item No. 700154

Overview

Contents: This vial contains 5 ml of 1 M Tris-HCl, pH 8.0.
Stability: ≥18 months at -80°C

Laboratory Procedures

Prior to performing the Methyltransferase Fluorometric Assay Kit (Item No. 700150), dilute 4 ml of (10X) buffer with 36 ml of HPLC-grade water. This final buffer (0.1 M Tris-HCl, pH 8.0) will be used to prepare the resorufin standards (Item No. 700152). When stored at 4°C, this diluted buffer is stable for at least six months.

Cautions

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work exclusively with the Methyltransferase Fluorometric Assay Kit from Cayman (Item No. 700150). This product may not perform as described if used with other assay reagents or protocols.

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, 08/31/2015

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