

PRODUCT INFORMATION



Cell-Based Assay IBMX Solution (1,000X)

Item No. 10008978

Overview

Contents: This vial contains 500 μ l of 1,000X IBMX stock solution.
Storage: -20°C
Stability: \geq 3 years

Laboratory Procedures

Cell-based assay IBMX Solution (1,000X) has been formulated and tested to work with Cayman's Adipogenesis Assay Kit (Item No. 10006908) and EP4 Receptor (rat) Activation Assay Kit (cAMP) (Item No. 600410). Detailed instructions for its use are contained in the respective Cayman Assay Kit booklet.

To prepare 10 ml of induction culture medium, add 10 μ l of this IBMX (1,000X) solution, along with 10 μ l of Insulin Solution (1,000X) (Item No. 10008979) and 10 μ l of Adipogenesis Assay Dexamethasone Solution (1,000X) (Item No. 10008980) to DMEM containing 10% FBS. To prepare 10 ml of 2X IBMX Solution for EP4 Receptor (rat) Activation Assay Kit (cAMP) (Item No. 600410), add 10 μ l of this stock solution to the culture medium used for the cells. To avoid frequent freeze and thaw of the stock solution, we recommend that you make small aliquots for long-term storage.

Precaution

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, 04/22/2024

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