

# PRODUCT INFORMATION



## Adipolysis Assay IBMX Solution (1,000X)

Item No. 10009948

### Overview and Properties

---

**Contents:** This vial contains 100  $\mu$ l of 1,000X IBMX stock solution.  
**Storage:** -20°C  
**Stability:**  $\geq$ 3 years

### Procedures

---

To prepare 10 ml of induction culture medium, add 10  $\mu$ l of this IBMX (1,000X) solution, along with 10  $\mu$ l of Insulin Solution (1,000X) (Item No. 10009949) and 10  $\mu$ l of Dexamethasone Solution (1,000X) (Item No. 10009950) to DMEM containing 10% FBS. For long term storage, we recommend that you make small aliquots to avoid frequent freezing and thawing of the stock solution.

### CAUTIONS

---

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work with the Adipolysis Assay Kit from Cayman (Item No. 10009381). 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, 10/25/2021

**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