

# PRODUCT INFORMATION



## FFA Assay Buffer

Item No. 700311

### Overview

---

**Contents:** The vial contains 40 ml of 50 mM Tris-HCl, pH 8.6, containing 5 mM magnesium chloride.  
**Stability:** ≥1 year at -20°C

### Laboratory Procedures

---

Prior to use in the Free Fatty Acid Fluorometric Assay Kit (Item No. 700310), thaw the Assay Buffer at room temperature. This buffer should be used to reconstitute the FFA Assay Cofactor Mixture 1 (Item No. 700312). The thawed buffer is stable at 4°C for at least six months.

### Cautions

---

For laboratory research use only. This reagent is not for human or veterinary use. 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, 09/08/2015

#### 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