

PRODUCT INFORMATION

FFA Assay Sample Buffer

Item No. 700317



Overview

Contents: The vial contains 15 ml of 50 mM Tris-HCl, pH 7.4, 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 Sample Buffer at room temperature. This buffer should be used to prepare the standard curve and to dilute samples. The thawed buffer is stable for at least six months at 4°C.

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