

PRODUCT INFORMATION



FFA Assay Cofactor 3

Item No. 700305

Overview

Contents: The vial contains 15 mg of coenzyme A.
Stability: ≥1 year at -20°C

Laboratory Procedures

Prior to use in the Free Fatty Acid Fluorometric Assay Kit (Item No. 700310), weigh out 5 mg of FFA Assay Cofactor 3 into a separate test tube to evaluate 60 wells. Dissolve with 3 ml of FFA Assay Buffer (Item No. 700311). Add this solution to the reconstituted vial of FFA Assay Cofactor Mixture 1 (Item No. 700312) and FFA Assay Cofactor 2 (Item No. 700304). Gently mix FFA Assay Cofactor Mixture 1, FFA Assay Cofactor 2, and FFA Assay Cofactor 3 by inversion. The combined Cofactor Mixture is stable for three hours. If additional wells are being utilized, weigh out additional FFA Assay Cofactor 3.

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 Free Fatty Acid Fluorometric Assay Kit from Cayman (Item No. 700310). 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