

# PRODUCT INFORMATION



## FFA Assay Cofactor 2

Item No. 700304

### Overview

---

**Contents:** The vial contains a lyophilized powder of ATP.  
**Stability:** ≥1 year at -20°C

### Laboratory Procedures

---

Prior to use in the Free Fatty Acid Fluorometric Assay Kit (Item No. 700310), reconstitute one vial with 3 ml of FFA Assay Buffer (Item No. 700311). The reconstituted mixture is stable for three hours on ice. Cofactor 2 will be combined with the reconstituted FFA Assay Cofactor Mixture 1 (Item No. 700312) and FFA Assay Cofactor 3 (Item No. 700305) prior to utilizing in the assay. One vial is enough to assay 60 wells. If additional wells are being utilized, then reconstitute a second vial.

### 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