

# PRODUCT INFORMATION



## Lipid Droplets Assay Dye Extraction Solution

Item No. 600046

### Overview

---

**Contents:** This vial contains 30 ml of dye extraction solution.  
**Storage:** Room temperature  
**Stability:** ≥2 years

### Laboratory Procedures

---

This dye extraction solution is supplied ready to use. For a 96-well plate, add 100 µl of the solution to each well and gently shake on an orbital plate shaker for up to 30 minutes before reading the absorbance with a plate reader. For 6-, 12-, or 24-well plates, add 1ml, 0.5 ml, and 0.25 ml of the dye extraction solution respectively. This solution is flammable and should be kept away from open flames.

### Precaution

---

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, 05/18/2021

#### 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