

# PRODUCT INFORMATION



## Lipid Droplets Assay Fixative (10X)

Item No. 10008981

### Overview

---

**Contents:** This vial contains 10 ml of 10X formaldehyde stock solution.  
**Storage:** Room temperature  
**Stability:** ≥2 years

### Laboratory Procedures

---

To make 10 ml of working solution (enough for one 96-well plate), add 1 ml of this stock solution to 9 ml of PBS Buffer, pH 7.2. The stock solution is hazardous and should be handled with care.

### Precaution

---

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