

# PRODUCT INFORMATION

## Cysteinyl Leukotriene ELISA Standard

Item No. 400394



### Overview

---

**Contents:** This vial contains 500 µl of CysLT as a solution in ethanol.  
**Storage:** -80°C  
**Stability:** ≥1 year

### Laboratory Procedures

---

Cysteinyl Leukotriene ELISA Standard has been tested and formulated to work exclusively with Cayman's Cysteinyl Leukotriene ELISA Kit (Item No. 500390). Please visit [www.caymanchem.com/catalog/kit number 500390](http://www.caymanchem.com/catalog/kit%20number%20500390) for the kit protocol, procedures, and product handling.

### Precaution

---

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, 02/17/2022

**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](http://WWW.CAYMANCHEM.COM)