

PRODUCT INFORMATION



Lipoxygenase (potato) Screening Enzyme

Item No. 60401

Overview

Contents:	This vial contains purified potato 5-lipoxygenase (5-LO) supplied in 0.1 M Tris-HCl (pH 7.4) containing 0.1% BSA.
Storage:	-80°C (as supplied)
Stability:	As supplied, 1 year from the QC date provided on the Certificate of Analysis, when stored properly

Laboratory Procedures

5-LO is relatively unstable at room temperature; keep the stock vial of the enzyme on ice (0-4°C) at all times. This vial of 5-LO has been formulated to be used for screening 5-LO specific inhibitors using the Lipoxygenase Inhibitor Screening Assay (Item No. 760700). Dilute 180 µl of 5-LO with 8.82 ml of diluted Lipoxygenase Inhibitor Screening Assay Buffer (Item No. 760710) prior to assaying. This is enough enzyme for use in 96 wells. The diluted enzyme is stable for 1 hour when stored on ice. When following the Lipoxygenase Inhibitor Screening Assay protocol, a 90 µl aliquot of this diluted enzyme causes a final absorbance of 0.25-0.29 using linoleic acid as the substrate. 5-LO prefers the linoleic acid substrate, therefore when utilizing this enzyme with the assay do not use the arachidonic acid as a substrate.

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, 01/26/2017

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