

CERTIFICATE of ANALYSIS



5-Lipoxygenase (human, recombinant)

Item No. 60402 • Batch No. 0821947 • Expires 27MAR2027

Overview

Formulation:	100 mM Tris-HCl, pH 8.0, with 5 mM EGTA, 1mM calcium chloride, and 30% glycerol
Concentration:	19.3 mg/mL
Storage/stability:	-80°C/9mo
Specific activity:	43.30 U/mg
Unit definition:	One unit of enzyme consumes one nmol of oxygen per minute at 25°C in 50 mM Tris-HCl buffer, pH 7.5, with 100 µM arachidonate, 2 mM calcium chloride, 1 mM ATP, 4.6 µM 13(S)-HpODE, and 250 µM oxygen. 5-Lipoxygenase activity is measured at 25°C by monitoring oxygen consumption using a Gilson Model 5/6 H oxygraph equipped with a Clark oxygen electrode.
Activity:	835.00 U/ml
Purity:	Clarified lysate

Reviewed and approved by: Beth Hurst

Approved date: 14MAY2026

Elizabeth Hurst

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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