

CERTIFICATE of ANALYSIS



PLA2 (human, recombinant Group XV)

Item No. 43839 • Batch No. 0811746 • Expires 27APR2027

Overview

Formulation:	50 mM HEPES, pH 7.5, 125 mM sodium chloride
Concentration:	2.2 mg/mL
Storage/stability:	-80°C/1yr
Applications:	enzyme activity assay, WB
Specific activity:	3,699.00 U/mg
Unit definition:	One unit is defined as the amount of enzyme required to hydrolyze 1 nmol of diheptanoyl thio-PG (Item No. 38436) per minute at 25 °C in 25 mM Tris, pH 7.5, 100 mM potassium chloride, 0.3 mM Triton X 100 in the presence of 1 mM DTNB and 1.25 mM diheptanoyl thio-PG.
Activity:	8,064.00 U/ml
Purity:	≥80% estimated by SDS-PAGE

Reviewed and approved by: Beth Hurst

Approved date: 28APR2026

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