

PLAQPRO™ Lp-PLA₂ Assay

Enzyme activity kit to measure Lp-PLA₂ activity in serum samples

The PLAQPRO™ Lp-PLA₂ Assay measures the enzyme activity of Lp-PLA₂ (lipoprotein-associated phospholipase A2) in serum samples. The assay is designed to run on automated clinical chemistry analyzers.

Approximately two-thirds of heart attacks and strokes are caused by the rupture of an unstable, rupture-prone atherosclerotic plaque. Unstable plaques are characterized by high levels of inflammatory cells and inflammatory mediators.

Lipoprotein-associated phospholipase A2 (Lp-PLA₂) is produced by inflammatory cells and circulates primarily bound to low-density lipoprotein (LDL) and is to a lesser extent associated with high-density lipoprotein (HDL) in human plasma. LDL oxidation is an early key event in the pathogenesis of atherosclerosis. Elevated Lp-PLA₂ levels have been found in atherosclerotic plaques and rupture lesions.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-KIT179-0200	200 tests
-----------------	-----------

Manuals, SDS & CofA

[View Online »](#)

- Cost-efficient alternative to competitor's Lp-PLA₂ detection system
- Wide analytical measurement range from 0 - 400.0 nmol/min/mL
- Analytical sensitivity of 0.236 nmol/min/mL
- High specificity and reproducibility, lot-after-lot
- Ready-to-use calibration standards and QC controls are provided
- Designed to run on automated clinical chemistry analyzers



Handling & Storage

Use/Stability	Store all components, with the exception of Reagent 1, at -80°C until the kit's expiration date. Reagent 1 may be stored at or below -20°C. Once thawed, the components can be stored at 2-8°C for up to 4 weeks.
Handling	Avoid freeze/thaw cycles.
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Colorimetric detection, HTS
Application Notes	For the measurement of enzyme activity of Lp-PLA ₂ (lipoprotein-associated phospholipase A2) in serum samples.
Contents	Lp-PLA ₂ assay buffer Lp-PLA ₂ assay substrate 5 calibrator standards QC low control QC high control
Sensitivity	0.236 nmol/min/mL
Species Reactivity	Human
Technical Info / Product Notes	<p>The PLAQPRO™ Lp-PLA₂ Assay is a colorimetric readout enzyme activity assay. During the test, Lp-PLA₂ in serum samples hydrolyze the S_N2 position of the substrate, 1-myristoyl-2-(4-nitrophenylsuccinyl)phosphatidylcholine. The resulting colored reaction product, 4-nitrophenol, is measured spectrophotometrically and the rate of formation of 4-nitrophenol is determined over successive measurements.</p> <p>The assay kit includes five Lp-PLA₂ calibrator standards for use in generating a standard curve fit of change in absorbance versus Lp-PLA₂ activity level in nmol/min/mL. The activity of Lp-PLA₂ in serum samples is interpolated from the calibration curve.</p>
UniProt ID	Q13093



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)