

Product Information Sheet

Ordering Information

Product Number: V100080

Product Name: Anti-Platelet-activating factor acetylhydrolase (Lp-PLA2) antibody *Mouse anti-human, monoclonal IgG1*

Unit Size: 50 ug

Storage Conditions: Freeze (<-15 °C)

Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Appearance: Solid

Molecular Weight: N/A

Soluble In: Water

Excitation Wavelength: N/A

Emission Wavelength: N/A

Application Notes

Lipoprotein-associated phospholipase A2 (Lp-PLA2), or platelet-activating factor acetylhydrolase (PAF-AH), is a 45 kDa phospholipase enzyme produced by inflammatory cells such as mature macrophages and activated platelets. Lp-PLA2 is typically found present in the blood coupled to lipoproteins: 80% to LDL and 20% to HDL. It is essential for the hydrolysis of oxidized phospholipids, yielding pro-inflammatory lysophospholipids and oxidized fatty acids. In clinically biology, Lp-PLA2 is a significant biomarker associated with the development of atherosclerosis and cardiac disease. Anti-Lp-PLA2 monoclonal antibodies are highly specific for the detection of human Lp-PLA2 immunogens. They can be used for a broad range of immunoassays such as chemiluminescent immunoassay (CLIA), latex enhanced turbidimetric immunoassay (LETIA), ELISA and rapid sandwich immunoassays.