

Echelon Biosciences Inc.

Certificate of Analysis

Product Name:	Human Lysosomal Phospholipase A2, active
Catalog Number:	E-7000
Size:	10 µg, Lyophilized Enzyme
Lot #	XCM072017L-21
Description:	Human β Lysosomal Phospholipase A2 (hLPLA2) with C-terminal 6-His tag was expressed in Hek 293 cells and purified using nickel-NTA chromatography.
Formulation:	Contains HEPES, Trehalose, Dextrose, NaCl, and Tween 20.
Purity:	$\geq 90\%$ by SDS-PAGE
Storage/Stability:	Store enzyme at -80°C . The lyophilized enzyme is stable for at least six months from date of arrival. Rehydrate in 50 μL ddH ₂ O on ice, Aliquot to avoid repeated freeze/thaw cycles. Store aliquots at -80°C .
Specific Activity:	8358 nmol/min/mg at 37°C . Specific activity is determined using 1 $\mu\text{g/mL}$ hLPLA2 enzyme in reaction buffer containing 35 mM sodium citrate, pH 4.5, 10 $\mu\text{g/mL}$ BSA, and 200 μM para-Nitrophenyl butyrate (PNPB or 4NPB) as substrate. Absorbance for product 4-nitrophenol (4NP) is measured at 401 nm immediately after equal volume of cold 0.2 M Tris (pH 9.0 at RT) is added to reactions to bring pH up to show color.