

SLPI (human), ELISA kit

Rapid ELISA kit to quantify SLPI, a secretory leukocyte peptidase inhibitor and unique cancer biomarker.

The SLPI ELISA kit is a complete, colorimetric, immunometric immunoassay kit for the quantitative determination of SLPI in serum with results in just 2.5 hours.

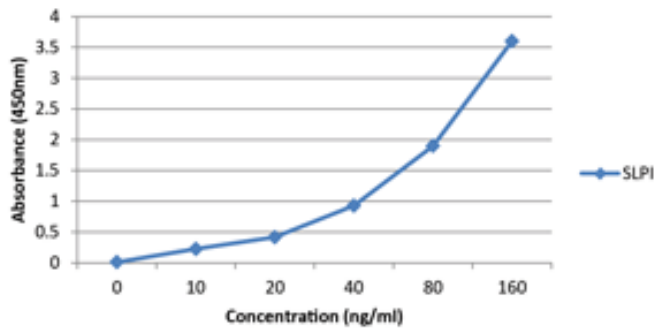
The ALP (antileukoproteinase) or SLPI (secretory leukocyte peptidase inhibitor) protein is a secreted inhibitor which protects epithelial tissues from serine proteases. It is found in various secretions including seminal plasma, cervical mucus, and bronchial secretions, and has affinity for trypsin, leukocyte elastase, and cathepsin G. Its inhibitory effect contributes to the immune response by protecting epithelial surfaces from attack by endogenous proteolytic enzymes. The ALP protein is also thought to have broad-spectrum antibiotic activity.

Manuals, SDS & CofA

[View Online »](#)

- Rapid – results in <2.5 hours from up to 40 samples in duplicate
- Convenient – user-friendly alternative to mass spectrometry
- Easy-to-use – simple procedure, liquid color-coded reagents and pre-coated plates to reduce error
- Quantitative – fully quantitative results surpass semi-quantitative Western blot analysis
- Sensitive – measure as little as 1 ng/ml of SLPI (ALP)

SLPI Standard Curve



Handling & Storage

Use/Stability	Store all components at +4°C
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

 RUO - Research Use Only

Product Details

Alternative Name	Secretory leukocyte proteinase inhibitor, ALP, Antileukoproteinase
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of human SLPI (ALP, Antileukoproteinase) in serum.
Assay Time	2.5 hours
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	Microtiter Plate, Standards, Calibrators, Antibody-conjugate, TMB Reagent, Stop Solution, Wash Buffer Concentrate
GenBank ID	758101
Gene/Protein Identifier	6590 (Gene ID), OTTHUMP00000031775 (Protein ID)
Sensitivity	1 ng/ml (range 10-160ng/ml)
Species Reactivity	Human
UniProt ID	P03973
Wavelength	450 nm



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Last modified: May 29, 2024