

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.

Ordering Information

[Order Online »](#)

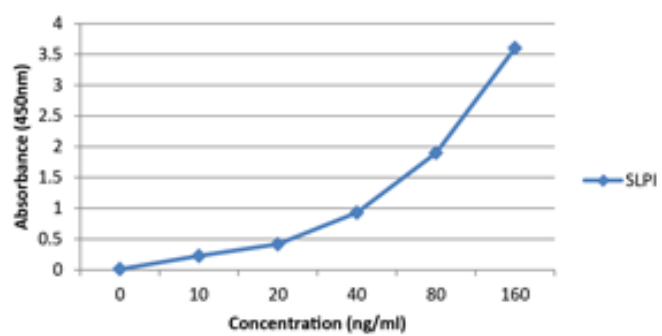
ADI-900-222-0001	96 wells
------------------	----------

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](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)