

IL-8 (human), ELISA kit

Ultrasensitive ELISA, measuring as low as 1 pg/ml of IL-8 in just 3 hours

The IL-8 (human), ELISA kit is a colorimetric immunometric enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 450 nm. Ultrasensitive measurement of < 1 pg/mL IL-8 with excellent lot-to-lot reproducibility.

Interleukin-8 (IL-8) is a cytokine produced by monocytes, macrophages, neutrophils, lymphocytes, dermal fibroblasts, keratinocytes, vascular endothelial cells, melanocytes, hepatocytes and various tumor cell lines. It is also a member of the CXC family of chemokines, displaying chemotactic activity for neutrophils and T cells.

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-900-156A

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Ultrasensitive measurement, < 1 pg/ml IL-8
- Fast results < 3 hours from up to 37 samples in duplicate
- Reproducible and fully validated in-house in a variety of sample types
- Ready-to-use reagents and pre-coated plates reduce errors and save your time

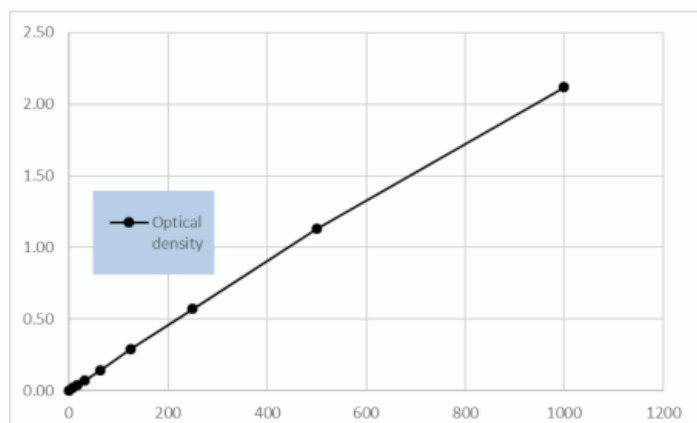


Figure 1. Typical Standard Curve for ADI-900-156A

Handling & Storage

Use/Stability Store all components at +4°C, except standard at -20°C.

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-8

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human IL-8 in culture supernatants, plasma, and serum.

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter Plate, Conjugate, Antibody, Assay Buffer 13, Wash Buffer Concentrate, Standard, TMB Substrate, Stop Solution 2, Assay Layout Sheet, Plate Sealer

Sensitivity 0.72 pg/ml (range 7.8-1,000 pg/ml)

Species Reactivity Human

Technical Info / Product Notes **Cited samples:**
[Cytokine ELISAs Cited Samples](#)

UniProt ID P10145

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)