## 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 precoated plates reduce errors and save your time

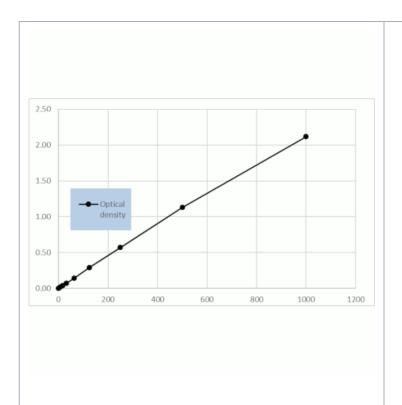


Figure 1. Typical Standard Curve for ADI-900-156A

## **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

Last modified: May 29, 2024

