



Analyze and Detect IL-6

High-Sensitivity IL-6 ELISA Kit

Most Sensitive Colorimetric Kit for IL-6 Quantification

IL-6 (human), High Sensitivity ELISA Kit (ENZ-KIT178)

Interleukin-6 (IL-6) is a cytokine critical to the regulation of the immune and hematopoietic systems. Human IL-6 is expressed by T-cells, mast cells, monocytes, macrophages, fibroblasts, endothelial cells, keratinocytes, and many tumor cell lines. It plays a role in acute phase reactions and response to injury and inflammation. Human IL-6 also stimulates the differentiation of B-cells for antibody production, promotes expansion of activated T-cells, expands hematopoietic cell production, and induces the expression of acute phase proteins.

We are proud to provide a comprehensive portfolio of products to detect and modulate the complex processes of immunological disease – our IL-6 High-Sensitivity ELISA Kit can detect up to 0.057 pg/mL of IL-6 in culture supernatants, plasma, serum, and urine.

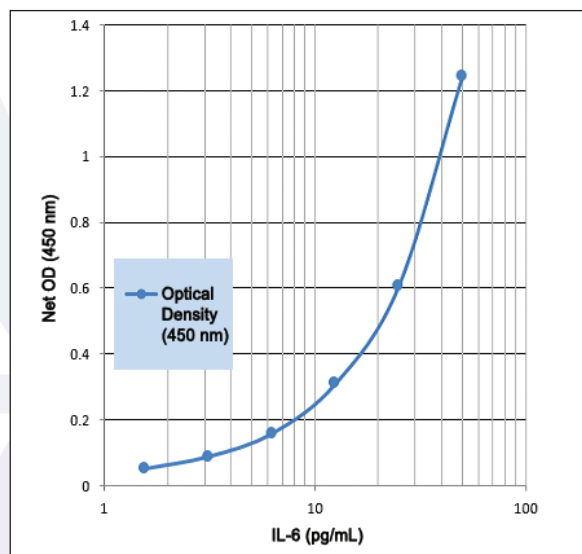
Key Features

- High-sensitivity ELISA enables quantification of low amounts of IL-6
- Low cross-reactivity with other interleukins and similar proteins
- Easy-to-use liquid color-coded reagents reduce errors
- Rapid quantitative results in 3 hours
- High-throughput immunoassay measures up to 40 samples in duplicate per 96 well kit
- Ready-to-use reagents and pre-coated plates reduce errors and save time

RELATED PRODUCTS

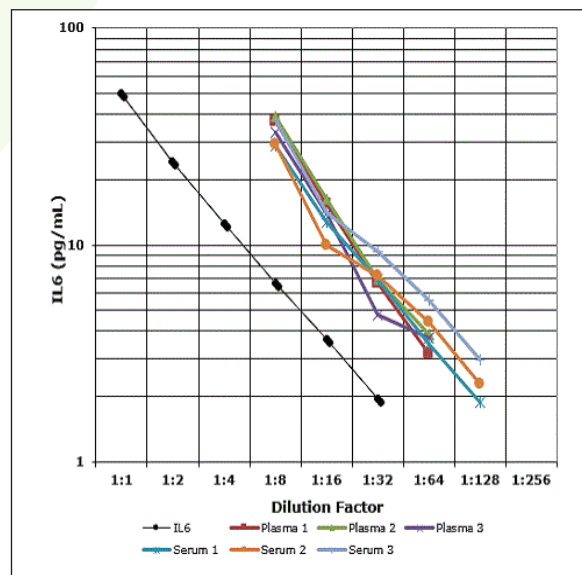
Product Name	Size	Product #
IL-6 (human) ELISA Kit	96 wells	ADI-900-033
IL-6 (human) ELISA Kit	5x96 wells	ADI-901-033
CRP (human) ELISA Kit	96 wells	ENZ-KIT102

Highly Sensitive Measurement of IL-6



Standard Curve for IL-6 High-Sensitivity ELISA Kit (ENZ-KIT178)

IL-6 ELISA Shows Consistent Results Across Sample Types



Advance Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
 10 Executive Boulevard
 Farmingdale, NY 11735
 Phone: 800.942.0430
 Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
 Industriestrasse 17
 CH-4415 Lausen, Switzerland
 Phone: +41 61 926 8989
 Fax: +41 61 926 8979
info-eu@enzolifesciences.com