

IGF-1 (human), ELISA kit

Highly sensitive ELISA, measuring as little as 34.2 pg/ml of IGF-1 in just 4 hours.

The IGF-1 (human), ELISA kit is a colorimetric immunometric enzyme-linked immunosorbent assay kit with results in 4 hours. Absorbance is read at 450 nm. This kit offers ultra-sensitive measurement of as little as 34.2 pg/ml of IGF-1.

Insulin-like growth factor-I (IGF-I) is a polypeptide hormone similar in structure to insulin. It is involved in regulation of neuronal growth and development in the central and peripheral nervous system. IGF-1 is a potent inhibitor of apoptosis in neurons, and activates the Akt signaling pathway via the IGF-1R and Insulin receptors. The IGF-1R acts as a receptor for IGF-1 and the related IGF-2, and is subject to autophosphorylation on multiple tyrosine residues that regulate receptor activation and assembly of signal transduction complexes.

Citations: 21

[View Online »](#)

Ordering Information

[Order Online »](#)

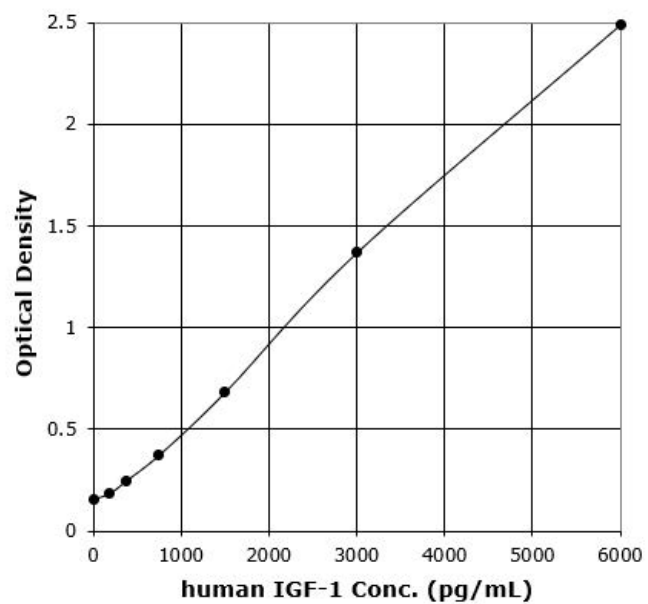
ADI-900-150

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive, detecting as low as 34.2 pg/ml IGF-1
- Broad dynamic range suitable for use with a variety of samples
- High throughput format with results for up to 40 samples in duplicate in just 4 hours
- Fully quantitative results that surpass semi-quantitative Western blot analysis



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Insulin-like growth factor 1

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human IGF-1 in plasma and serum. Cited sample types include culture supernatants and milk.

Assay Time 4 hours

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

Contents Microtiter plate, Conjugate, Antibody, Assay buffer 10 concentrate, Standard, Wash buffer concentrate, TMB substrate, Stop solution 2, Neutralizing reagent 2, 2N HCl

Crossreactivity Human IGF-1 (100%), human IGF-2 (0.11%) and <0.1%: IGFBP-2, IGFBP-3, IGFBP-4, mouse IGF-1, mouse IGF-2, HGH, HGF, PDGF-BB, human insulin

Sensitivity 34.2 pg/ml (range 187 – 6,000 pg/ml)

Species Reactivity Human

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

UniProt ID P05019

Wavelength



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-fr@enzolifesciences.com

450 nm

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: August 22, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.