



Sensitive Measurement of VEGF in 2.5 Hours

VEGF (human), ELISA Kit (ENZ-KIT156)

For the quantitative determination of human VEGF in serum, plasma, and culture supernatants

Vascular Endothelial Growth Factor (VEGF), originally known as vascular permeability factor (VPF), is a family of proteins that stimulate vasculogenesis and angiogenesis. The VEGF family comprises five members: VEGF-A, VEGF-B, VEGF-C, VEGF-D and PGF (placenta growth factor) 1-3. Being the first of the VEGF family members to be discovered, VEGF-A is often simply referred to as VEGF. The VEGF (human) ELISA kit is a sensitive colorimetric immunometric enzyme immunoassay kit with results in just 2.5 hours.

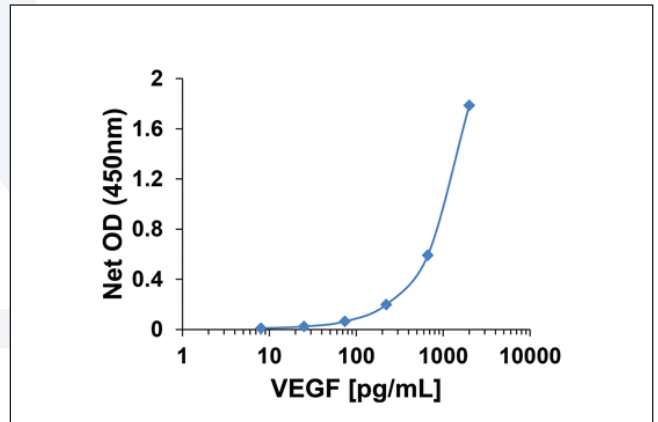
Key Features

- Sensitive Assay – Detect VEGF as low as 3.979 pg/ml
- Specific – 100% cross-reactivity to VEGF 165 with negligible cross-reactivity to related molecules including VEGF 121
- Flexible – Wide dynamic range ideal for various matrices including serum, plasma, and culture supernatants
- High-throughput – Assay up to 39 samples in duplicate with each kit

Areas of Research

- Asthma
- Autoimmune
- Cancer
- Cardiovascular
- Diabetes
- Inflammation
- Neuroscience
- Stem Cells

Highly Sensitive Measurement of VEGF



Standard curve experiment using VEGF (human), ELISA Kit (ENZ-KIT156).

RELATED PRODUCTS

PRODUCT NAME	PRODUCT #	PRODUCT NAME	PRODUCT #
Flt1 (human), (recombinant) (His-tag)	BML-SE445	VEGF monoclonal antibody (6B7)	ADI-905-164
VEGF 121 (human), (recombinant)	ENZ-PRT128	VEGF-C (human) ELISA Kit	ALX-850-306
VEGF 165 (human), (recombinant)	ENZ-PRT129	VEGFR-1 (human) ELISA Kit	ALX-850-264
VEGF A monoclonal antibody (4G3)	ADI-905-861	VEGFR2 (human), (recombinant) (GST-tag)	BML-SE404

Accelerate 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