

Product datasheet

Human HB EGF ELISA Kit ab100531

[2 References](#) [2 Images](#)

Overview

Product name	Human HB EGF ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 20 pg/ml
Range	16.38 pg/ml - 4000 pg/ml
Recovery	90 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	81.22	73% - 90%
Plasma	93.73	84% - 103%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's HB EGF Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human HB EGF in plasma and cell culture supernatants. (Human HB EGF concentration is pretty low in normal plasma, it may not be detected in this assay).

This assay employs an antibody specific for Human HB EGF coated on a 96-well plate. Standards and samples are pipetted into the wells and HB EGF present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human HB EGF antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of HB EGF bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect the endogenous Human HB-EGF in normal serum with ab100531, only in serum spiked with Human HB-EGF.

Notes Optimization may be required with urine samples.
The standard is affinity purified.

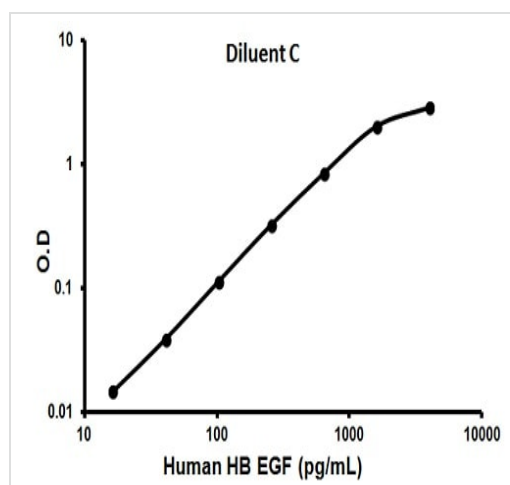
Platform Microplate

Properties

Storage instructions Store at -20°C. Please refer to protocols.

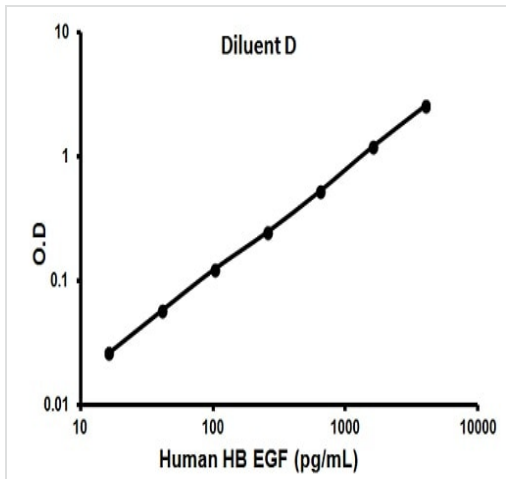
Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Human HB EGF	2 vials
HB EGF Microplate (12 x 8 wells)	1 unit
Recombinant Human HB EGF Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Images



Representative standard curve using ab100531

Typical standard curve



Representative standard curve using ab100531

Typical standard curve

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors