

Human beta IG H3 ELISA Kit (TGFBI) ab155426

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

Overview

Product name Human beta IG H3 ELISA Kit (TGFBI)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
overall				< 12%

Sample type Serum, Plasma, Cell culture media

Assay type Sandwich (quantitative)

Sensitivity < 25 pg/ml

Range 0.025 ng/ml - 6 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	= 92.44	76% - 113%
Plasma	= 86.91	69% - 104%
Cell culture media	= 104.8	96% - 114%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's beta IG H3 (TGFBI) Human ELISA (Enzyme-Linked Immunosorbent Assay) Kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of Human beta IG H3 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human beta IG H3 (TGFBI) coated on a 96-well plate. Standards and samples are pipetted into the wells and beta IG H3 present in a sample is bound

to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human beta IG H3 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 beta IG H3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization maybe required with urine samples.

Platform Microplate

Properties

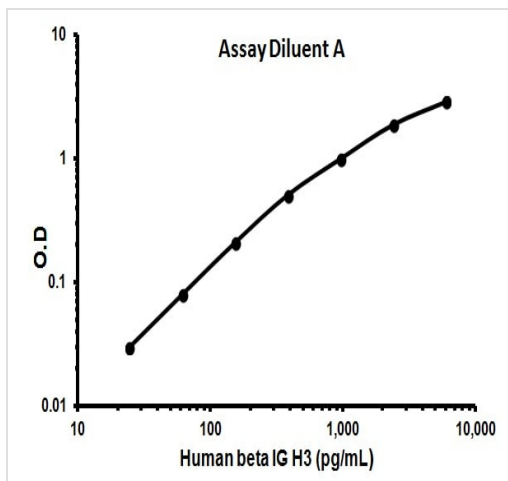
Storage instructions Store at -20°C. Please refer to protocols.

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
Assay Diluent A	1 x 30ml
beta IG H3 Microplate (12 x 8 wells)	1 unit
Biotinylated anti-Human beta IG H3	2 vials
Recombinant Human beta IG H3 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

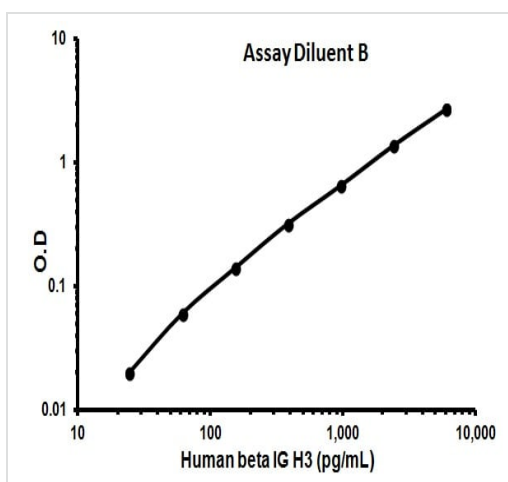
Cellular localization Secreted, extracellular space, extracellular matrix.

Images

Representative standard curve using ab155426 - Assay Diluent B



Typical Standard Curve



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