

Human ErbB3 / HER3 ELISA Kit ab100511

1 References 2 Images

Overview

Product name	Human ErbB3 / HER3 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 4 pg/ml
Range	3.43 pg/ml - 2500 pg/ml
Recovery	> 90 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	100.1	78% - 113%
Serum	95.12	79% - 111%
Plasma	88.87	71% - 106%

Assay duration	Multiple steps standard assay
Species reactivity	<b>Reacts with:</b> Human
Product overview	<p>Abcam's ErbB3 (Epidermal Factor Growth Factor Receptor 3) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of human ErbB3 in serum, plasma, and cell culture supernatants.</p> <p>This assay employs an antibody specific for human ErbB3 coated on a 96-well plate. Standards and samples are pipetted into the wells and ErbB3 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-human ErbB3 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 ErbB3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
Notes	Optimisation may be required with urine samples.
Platform	Microplate

## Properties

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

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human ErbB3	2 vials
ErbB3 Microplate (12 x 8 wells)	1 unit
Recombinant Human ErbB3 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function** Binds and is activated by neuregulins and NTAK.

**Tissue specificity** Epithelial tissues and brain.

**Involvement in disease** Defects in ERBB3 are the cause of lethal congenital contracture syndrome type 2 (LCCS2) [MIM:607598]; also called Israeli Bedouin multiple contracture syndrome type A. LCCS2 is an autosomal recessive neurogenic form of a neonatally lethal arthrogryposis that is associated with atrophy of the anterior horn of the spinal cord. The LCCS2 syndrome is characterized by multiple joint contractures, anterior horn atrophy in the spinal cord, and a unique feature of a markedly distended urinary bladder. The phenotype suggests a spinal cord neuropathic etiology.

**Sequence similarities** Belongs to the protein kinase superfamily. Tyr protein kinase family. EGF receptor subfamily. Contains 1 protein kinase domain.

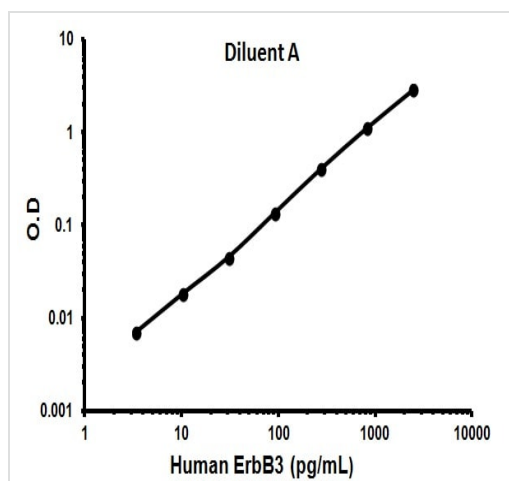
**Developmental stage** Overexpressed in a subset of human mammary tumors.

**Domain** The cytoplasmic part of the receptor may interact with the SH2 or SH3 domains of many signal-transducing proteins.

**Post-translational modifications** Ligand-binding increases phosphorylation on tyrosine residues and promotes its association with the p85 subunit of phosphatidylinositol 3-kinase.

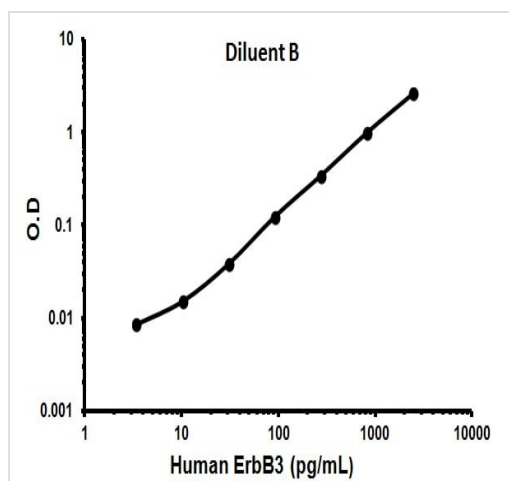
**Cellular localization** Secreted and Cell membrane.

## Images



TYPICAL STANDARD CURVE - Assay Diluent A

TYPICAL STANDARD CURVE - Assay Diluent A – Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



TYPICAL STANDARD CURVE - Assay Diluent B

TYPICAL STANDARD CURVE - Assay Diluent B – Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

**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