

Human BCA1 ELISA Kit ab99981

2 Images

Overview

Product name	Human BCA1 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 1.5 pg/ml
Range	1.37 pg/ml - 1000 pg/ml
Recovery	

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	95.62	84% - 105%
Serum	95.79	82% - 105%
Plasma	93.68	83% - 104%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's BCA1 (BLC) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human BCA1 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human BCA1 coated on a 96-well plate. Standards and samples are pipetted into the wells and BCA1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human BCA1 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 BCA1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Human BCA1 ELISA Kit (CXCL13) ([ab179881](#)) from our SimpleStep ELISA® range.

Notes Optimization may be required with urine samples.

Properties

Storage instructions

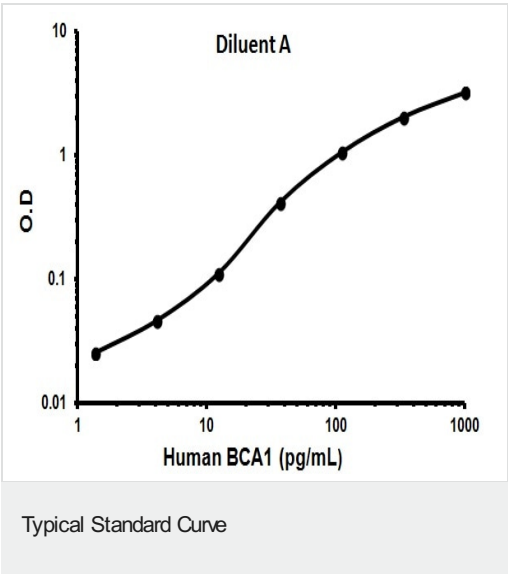
Store at -20°C. Please refer to protocols.

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
BCA1 Microplate (12 x 8 wells)	1 unit
Biotinylated anti-Human BCA1	2 vials
Recombinant Human BCA1 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

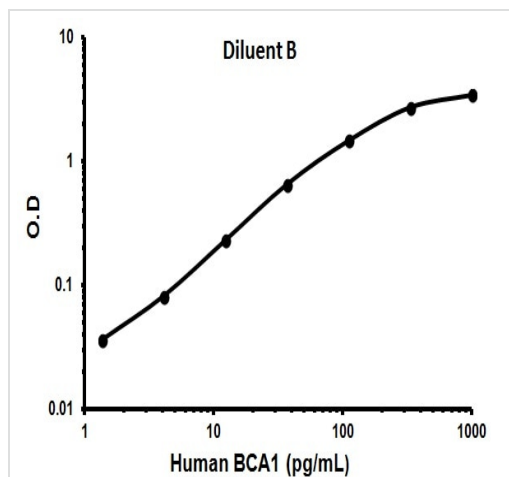
Relevance

B lymphocyte chemoattractant, independently cloned and named Angie, is a CXC chemokine strongly expressed in the follicles of the spleen, lymph nodes, and Peyer's patches. It preferentially promotes the migration of B lymphocytes (compared to T cells and macrophages), apparently by stimulating calcium influx into, and chemotaxis of, cells expressing Burkitt's lymphoma receptor 1 (BLR1). It may therefore function in the homing of B lymphocytes to follicles.

Images



Representative Standard Curve using ab99981.



Representative Standard Curve using ab99981.

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