

Human CEACAM1 ELISA Kit (BGP) ab99993

[2 Images](#)

Overview

| | |
|-------------------------|---|
| Product name | Human CEACAM1 ELISA Kit (BGP) |
| Detection method | Colorimetric |
| Sample type | Cell culture supernatant, Serum, Plasma |
| Assay type | Sandwich (quantitative) |
| Sensitivity | < 20 pg/ml |
| Range | 20.58 pg/ml - 15000 pg/ml |
| Recovery | > 83 % |

Sample specific recovery

| Sample type | Average % | Range |
|--------------------------|-----------|------------|
| Cell culture supernatant | 83.49 | 75% - 90% |
| Serum | 96.92 | 82% - 112% |
| Plasma | 86.5 | 70% - 105% |

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's CEACAM1 (BGP) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human CEACAM1 in serum, plasma, cell culture supernatant.

This assay employs an antibody specific for Human CEACAM1 coated on a 96-well plate. Standards and samples are pipetted into the wells and CEACAM1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human CEACAM1 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 CEACAM1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples.

Platform Microplate

Properties

Storage instructions

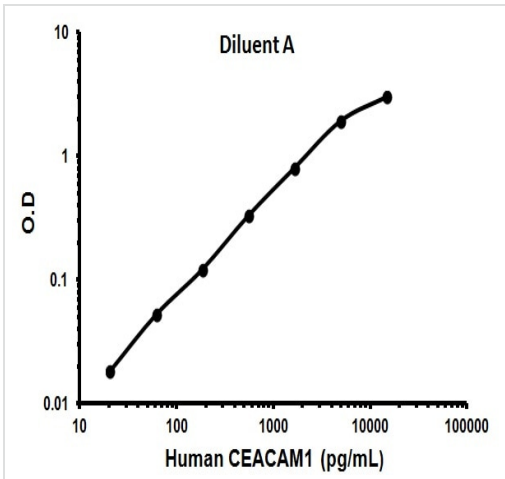
Store at -20°C. Please refer to protocols.

| Components | 1 x 96 tests |
|--|--------------|
| 400X 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 CEACAM1 | 2 vials |
| CEACAM1 Microplate (12 x 8 well strips) | 1 unit |
| Recombinant Human CEACAM1 Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

Relevance

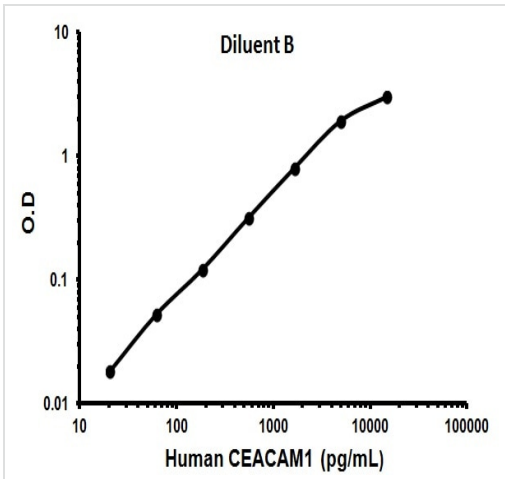
CEA-related cell adhesion molecules (CEACAM) belong to the carcinoembryonic antigen (CEA) family. It consists of seven CEACAM (CEACAM 1, CEACAM 3-CEACAM 8) and 11 pregnancy-specific glyco-protein (PSG 1-PSG 11) members. The CEA family proteins belong to the immunoglobulin (Ig) superfamily and are composed of one Ig variable-like (IgV) and a varying number (0-6) of Ig constant-like (IgC) domains. CEACAM molecules are membrane-bound either via a transmembrane domain or a glycosyl phosphatidyl inositol (GPI) anchor. CEACAM molecules are differentially expressed in epithelial cells or in leucocytes. Over-expression of CEA/CEACAM 5 in tumors of epithelial origin is the basis of its wide-spread use as a tumor marker. The function of CEACAM family members varies widely: they function as cell adhesion molecules, tumor suppressors, regulators of lymphocyte and dendritic cell activation, receptors of Neisseria species and other bacteria.

Images



Representative Standard Curve using ab99993.

Typical Standard Curve



Representative Standard Curve using ab99993.

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