

Human BMP7 ELISA Kit ab99985

★★★★★ 1 Abreviews 1 References 2 Images

Overview

Product name	Human BMP7 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 10 pg/ml
Range	8.23 pg/ml - 6000 pg/ml
Recovery	96 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	97.11	84% - 105%
Serum	96.2	85% - 104%
Plasma	95.37	86% - 104%

Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	<p>Abcam's BMP7 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human BMP7 in serum, plasma, and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human BMP7 coated on a 96-well plate. Standards and samples are pipetted into the wells and BMP7 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated antiHuman BMP7 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 BMP7 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
Notes	Optimization may be required with urine samples.
Platform	Microplate

Properties

Storage instructions

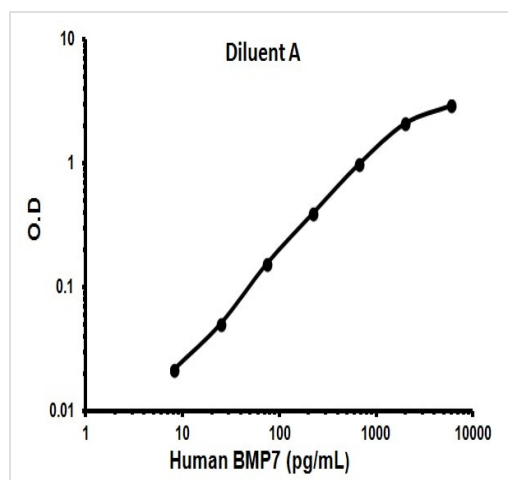
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human BMP7	2 vials
BMP7 Microplate (12 x 8 wells)	1 unit
Recombinant Human BMP7 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Relevance

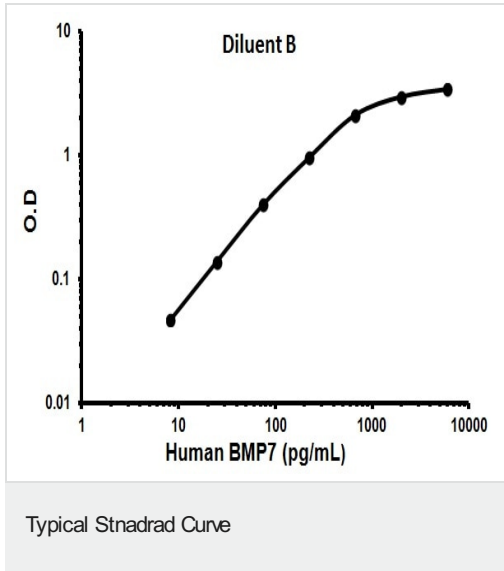
Bone Morphogenetic Protein 7 (BMP7), also known as osteogenic protein 1 (OP1), is produced from a DNA sequence encoding the human BMP2 signal peptide and human BMP2 propeptide (amino acid residues 1 to 282) fused to the human BMP7 mature subunit (amino acid residues 293 to 431). Mature human BMP7, generated after the proteolytic removal of the signal peptide and the propeptide, is a disulfide linked homodimeric protein, comprised of two 139 amino acid residue subunits. Bone Morphogenetic Proteins (BMP) are members of the TGF beta superfamily of cytokines that affect bone and cartilage formation. Similar to other TGF beta family proteins, BMPs are highly conserved across animal species.

Images



Typical Standard Curve

Representative Standard Curve using ab99985



Representative Standard Curve using ab99985

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