

Human Procalcitonin ELISA Kit ab100630

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

Overview

Product name	Human Procalcitonin ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 20 pg/ml
Range	27.43 pg/ml - 20000 pg/ml
Recovery	83 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	100.5	80% - 111%
Plasma	72.35	65% - 88%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Procalcitonin Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Procalcitonin in plasma and cell culture supernatants. (Human Procalcitonin concentration is pretty low in normal plasma, it may not be detected in this assay).

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

We have not been able to detect endogenous Human Procalcitonin in normal serum with ab100630, only in serum spiked with Human Procalcitonin.

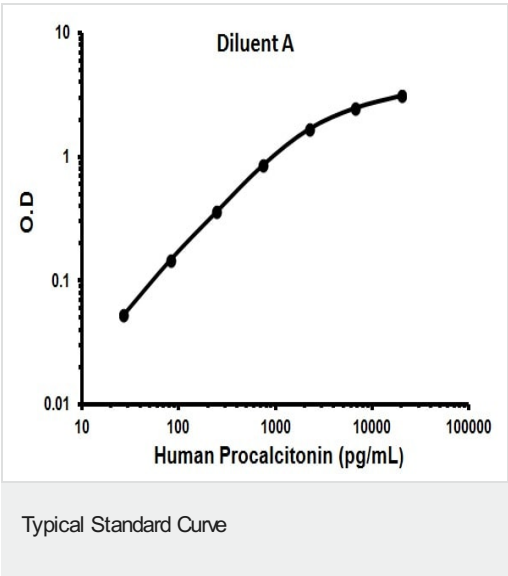
Notes	Optimization may be required with urine samples.
Platform	Microplate

Properties

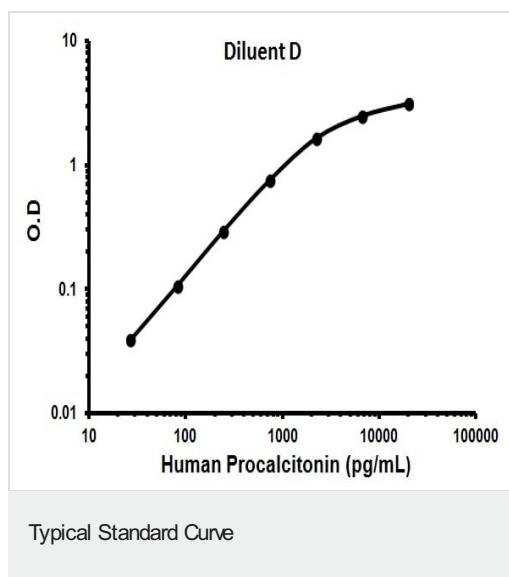
Storage instructionsStore at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
700X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human Procalcitonin	2 vials
Procalcitonin Microplate (12 x 8 wells)	1 unit
Recombinant Human Procalcitonin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Images



Representative Standard Curve using ab100630.



Representative Standard Curve using ab100630.

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