

Product datasheet

Human Furin ELISA Kit ab113322

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

Overview

Product name	Human Furin ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 15 pg/ml
Range	13.72 pg/ml - 10000 pg/ml
Recovery	84 %

Sample specific recovery

Sample type	Average %	Range
Serum	89.57	80% - 98%
Plasma	77.74	69% - 85%
Cell culture media	84.69	75% - 93%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Furin Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Furin in serum, plasma (collect plasma using heparin and EDTA as an anticoagulant. Citrate is not recommended) and cell culture supernatants.

This assay employs an antibody specific for Human Furin coated on a 96-well plate. Standards and samples are pipetted into the wells and Furin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Furin 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 Furin 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

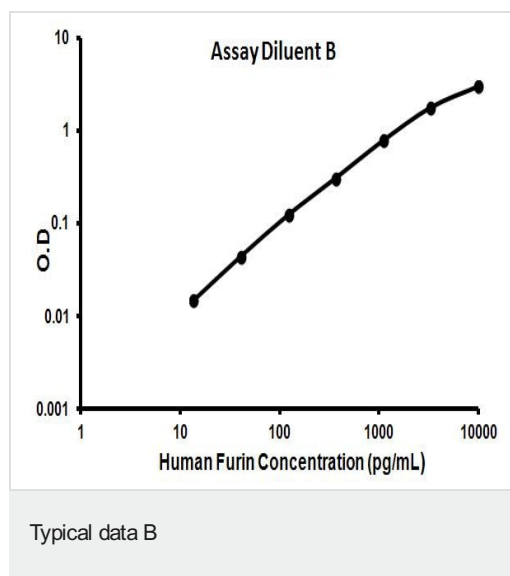
Properties

Storage instructions

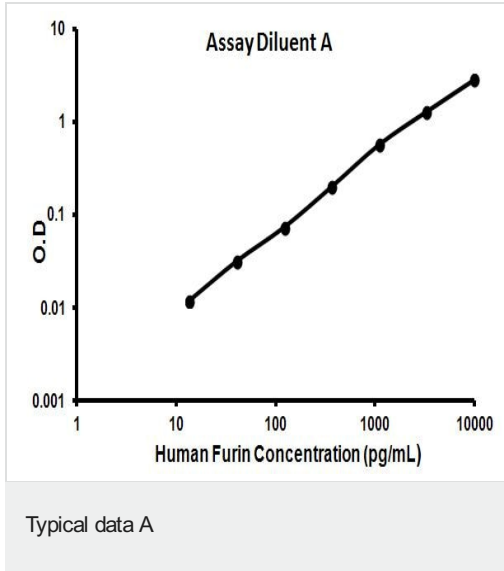
Store at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests
20X Wash Buffer		1 x 25ml
300X HRP-Streptavidin Concentrate		1 x 200µl
5X Assay Diluent B		1 x 15ml
Assay Diluent A		1 x 30ml
Biotinylated anti-Human Furin	Item F	2 vials
Furin Microplate (12 x 8 wells)	Item A	1 unit
Recombinant Human Furin Standard (lyophilized)		2 vials
Stop Solution		1 x 8ml
TMB One-Step Substrate Reagent		1 x 12ml

Images



Representative standard curve using ab113322



Representative standard curve using ab113322

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