

Human PR3 ELISA Kit ab226902

Recombinant SimpleStep ELISA®

1 References 6 Images

Overview

Product name	Human PR3 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Supernatant	8			3.7%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Supernatant	3			11%
Sample type	Cell culture supernatant, Cell culture extracts				
Assay type	Sandwich (quantitative)				
Sensitivity	150 pg/ml				
Range	1.2 ng/ml - 150 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %	Range	
	Cell culture supernatant		101	90% - 117%	
	Cell culture extracts		89	87% - 93%	
Assay time	1h 30m				
Assay duration	One step assay				
Species reactivity	Reacts with: Human				
Product overview	Human PR3 ELISA Kit (ab226902) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of PR3 protein in cell culture extracts and cell culture supernatant. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human PR3 with 150 pg/ml sensitivity.				

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Notes

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform

Pre-coated microplate (12 x 8 well strips)

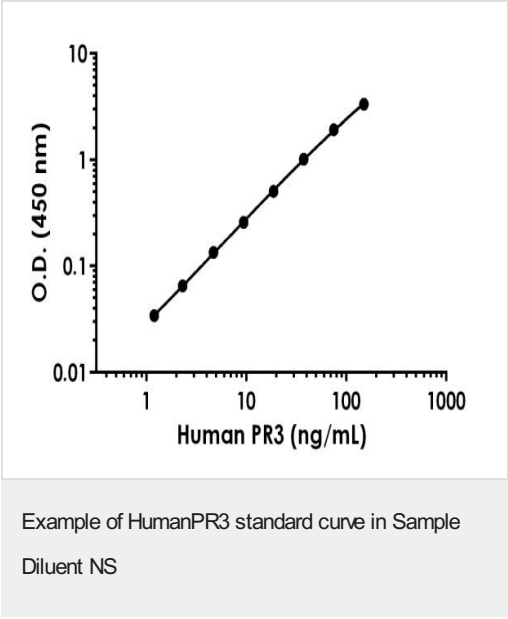
Properties

Storage instructions Store at +4°C. Please refer to protocols.

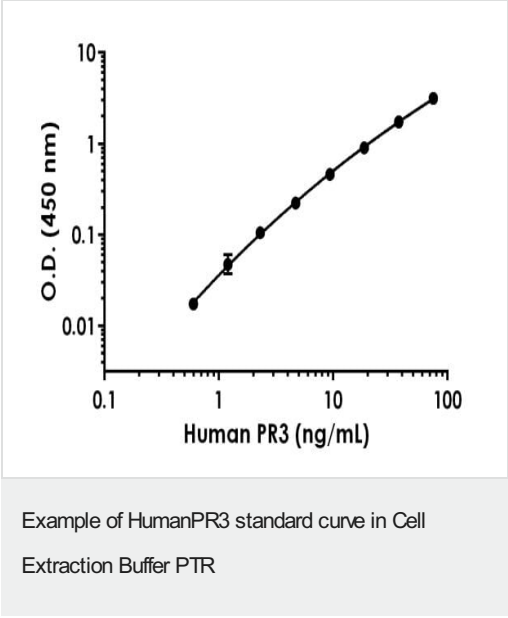
Components	1 x 96 tests
10X Human PR3 Capture Antibody	1 x 600µl
10X Human PR3 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent 4BI	1 x 6ml
Human PR3 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

Function	Polymorphonuclear leukocyte serine protease that degrades elastin, fibronectin, laminin, vitronectin, and collagen types I, III, and IV (in vitro) and causes emphysema when administered by tracheal insufflation to hamsters.
Sequence similarities	Belongs to the peptidase S1 family. Elastase subfamily. Contains 1 peptidase S1 domain.

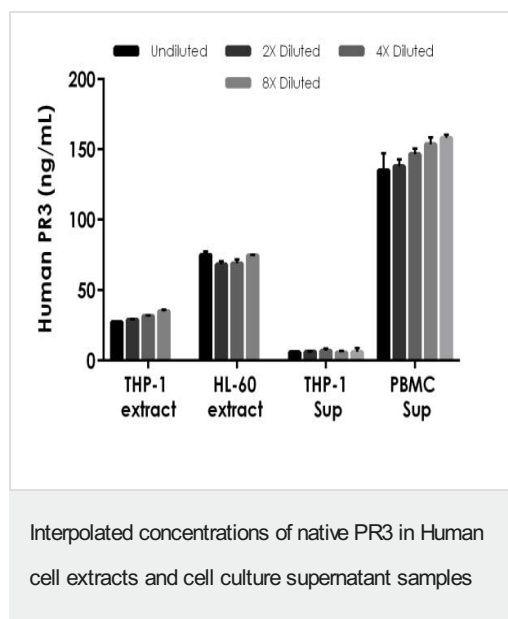
Images



Background-subtracted data values (mean +/- SD) are graphed.



Background-subtracted data values (mean +/- SD) are graphed.



The concentrations of PR3 were measured in duplicates, interpolated from the PR3 standard curves and corrected for sample dilution. Undiluted samples are as follows: THP-1 cell extract (100 μ g/ml); HL-60 (50 μ g/ml); untreated THP-1 cell supernatant (100 %) and untreated Human PBMC supernatant (50%). The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean PR3 concentration was determined to be 24.62 ng/mL in THP-1 cell extract, 72.65 ng/mL in HL-60 cell extract and 6.16 ng/mL in THP-1 supernatant and 146.21 ng/ml in PBMC supernatant.

Powered by
recombinant antibodies

Research with confidence
Consistent and reproducible results

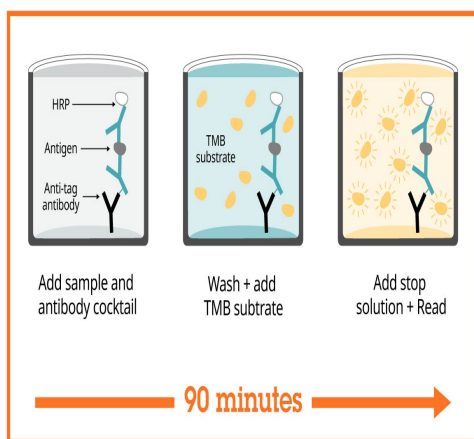
Long-term and scalable supply
Recombinant technology

Success from the first experiment
Confirmed specificity

Ethical standards compliant
Animal-free production

Sandwich ELISA - Human PR3 ELISA Kit
(ab226902)

To learn more about the advantages of recombinant antibodies see [here](#).



Sandwich ELISA - Human PR3 ELISA Kit
(ab226902)

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Get more done with SimpleStep ELISA



Easy to use
Single-wash 90-minute protocol



Flexible
Matched antibody pairs available



Precision antibodies
High sensitivity, specificity and reproducibility



Scalable
Now in 10-pack and 384-well formats

Sandwich ELISA - Human PR3 ELISA Kit
(ab226902)

To learn more about the advantages of SimpleStep ELISA® kits see [here](#).

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