

Human CD167α/DDR1 ELISA Kit ab260063

Recombinant SimpleStep ELISA®

4 Images

Overview

Product name	Human CD167a/DDR1 ELISA Kit
Detection method	Colorimetric
Precision	Intra-assay

Sample	n	Mean	SD	CV%
Supernatant	8			1.68%

Sample type	Cell culture supernatant
Assay type	Sandwich (quantitative)
Sensitivity	143.7 pg/ml
Range	312.5 pg/ml - 20000 pg/ml
Recovery	Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	89	% - %

Assay time	1h 30m
Assay duration	One step assay
Species reactivity	Reacts with: Human
Product overview	Human CD167a/DDR1 ELISA Kit (ab260063) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of CD167a/DDR1 protein in cell culture supernatant. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human CD167a/DDR1 with 20.53 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 Epithelial discoidin domain-containing receptor 1 (DDR1) is a tyrosine kinase receptor involved in cell attachment to the extracellular matrix. DDR1 is also involved in cell proliferation and differentiation. DDR1's contribution to remodeling of the extracellular matrix supports cell migration and wound healing.

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 DDR1 Capture Antibody	1 x 600µl
10X Human DDR1 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent CPI - HAMA Blocker (ab193969)	1 x 6ml
Human DDR1 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 May be involved in cell-cell interactions and recognition.

Tissue specificity Detected in T-47D, MDA-MB-175 and HBL-100 breast carcinoma cells, A431 epidermoid carcinoma cells, SW48 and SNU-C2B colon carcinoma cells and Hs 294T melanoma cells (at protein level). Expressed at low levels in most adult tissues and is highest in the brain, lung, placenta and kidney. Lower levels of expression are detected in melanocytes, heart, liver, skeletal muscle and pancreas. Abundant in breast carcinoma cell lines. In the colonic mucosa, expressed in epithelia but not in the connective tissue of the lamina propria. In the thyroid gland, expressed in the epithelium of the thyroid follicles. In pancreas, expressed in the islets of Langerhans cells, but not in the surrounding epithelial cells of the exocrine pancreas. In kidney, expressed in the epithelia of the distal tubules. Not expressed in connective tissue, endothelial cells, adipose

tissue, muscle cells or cells of hematopoietic origin.

Sequence similarities

Belongs to the protein kinase superfamily. Tyr protein kinase family. Insulin receptor subfamily.

Contains 1 F5/8 type C domain.

Contains 1 protein kinase domain.

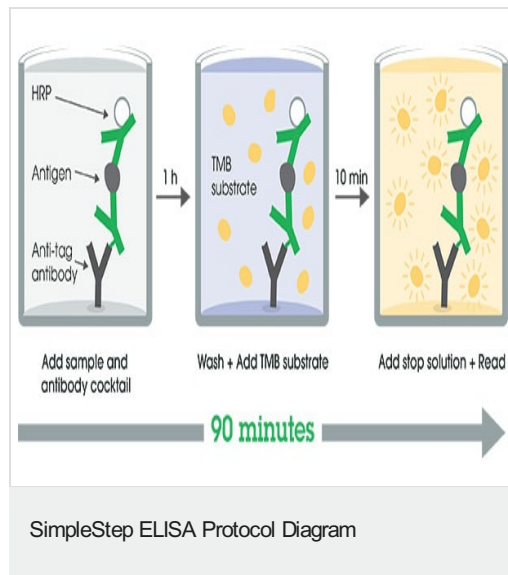
Domain

The Gly/Pro-rich domains may be required for an unusual geometry of interaction with ligand or substrates.

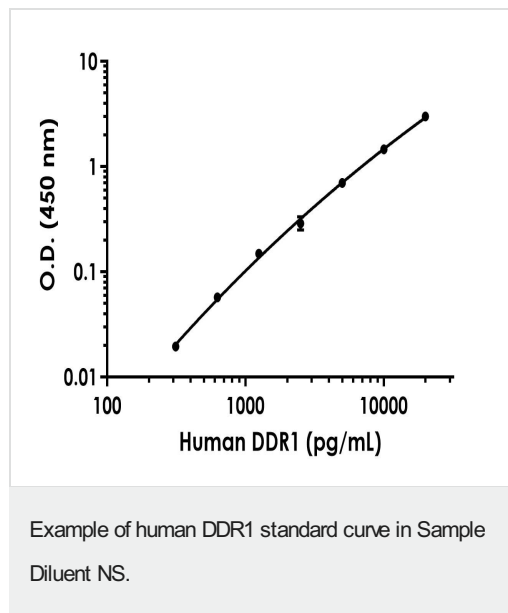
Cellular localization

Secreted and Membrane.

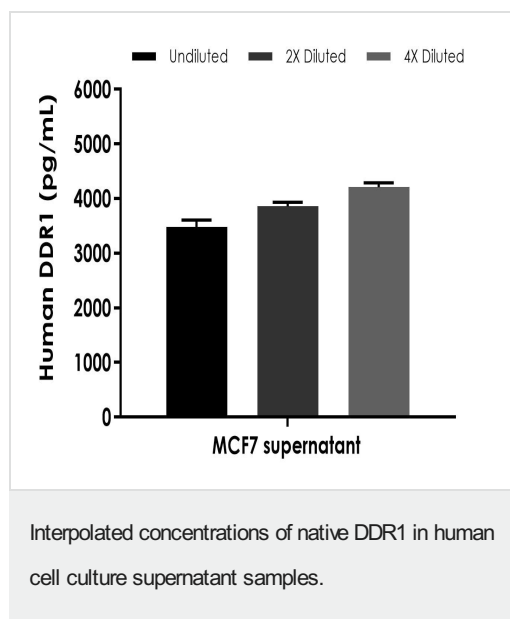
Images



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.



The DDR1 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean \pm SD) are graphed.



The concentrations of DDR1 were measured in duplicates, interpolated from the DDR1 standard curves and corrected for sample dilution. Undiluted samples are as follows: MCF7 supernatant 50%. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean DDR1 concentration was determined to 3,910.99 pg/mL in MCF7 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 DDR1 ELISA Kit
(ab260063)

To learn more about the advantages of recombinant antibodies 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