

Human JAM-A ELISA Kit (Junctional Adhesion Molecule) ab277079

Recombinant SimpleStep ELISA[®]

5 Images

Overview

Product name	Human JAM-A ELISA Kit (Junctional Adhesion Molecule)				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Serum	8			2.5%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Serum	3			6.4%
Sample type	Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	1.9 pg/ml				
Range	15.625 pg/ml - 1000 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Cell culture supernatant		94		86% - 103%
	Serum		102		86% - 107%
	Hep Plasma		91		88% - 96%
	EDTA Plasma		85		82% - 90%
	Cit plasma		96		81% - 113%

Assay time	1h 30m
Assay duration	One step assay
Species reactivity	Reacts with: Human
Product overview	Human JAM-A ELISA kit (ab277079) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Human JAM-A protein in human serum, plasma, and cell culture media. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human JAM-A with 1.9 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	Junctional Adhesion Molecule (JAM-A), a type I transmembrane protein, appears to play a role in epithelial tight junction formation. JAM-A emerges early in primordial forms of cell junctions and recruits PARD3. The association of the PARD6-PARD3 complex may prevent the interaction of PARD3 with JAM-A, thereby preventing tight junction assembly. JAM-A plays a role in regulating monocyte transmigration involved in integrity of epithelial barrier. JAM-A shares 68.4%, 71.8%, 74.6%, and 97.0% sequence homology with mouse, rat, bovine, and monkey, respectively.
--------------	---

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 JAM-A (Junctional Adhesion Molecule) Capture Antibody	1 x 600µl
10X Human JAM-A (Junctional Adhesion Molecule) Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BI	1 x 6ml
Human JAM-A (Junctional Adhesion Molecule) Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit

Components	1 x 96 tests
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

Seems to play a role in epithelial tight junction formation. Appears early in primordial forms of cell junctions and recruits PARD3. The association of the PARD6-PARD3 complex may prevent the interaction of PARD3 with JAM1, thereby preventing tight junction assembly (By similarity). Plays a role in regulating monocyte transmigration involved in integrity of epithelial barrier. Involved in platelet activation. In case of orthoreovirus infection, serves as receptor for the virus.

Sequence similarities

Belongs to the immunoglobulin superfamily.
Contains 2 Ig-like V-type (immunoglobulin-like) domains.

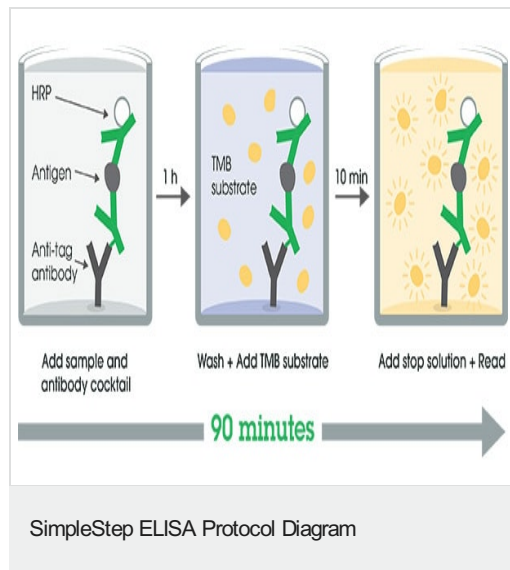
Post-translational modifications

N-glycosylated.

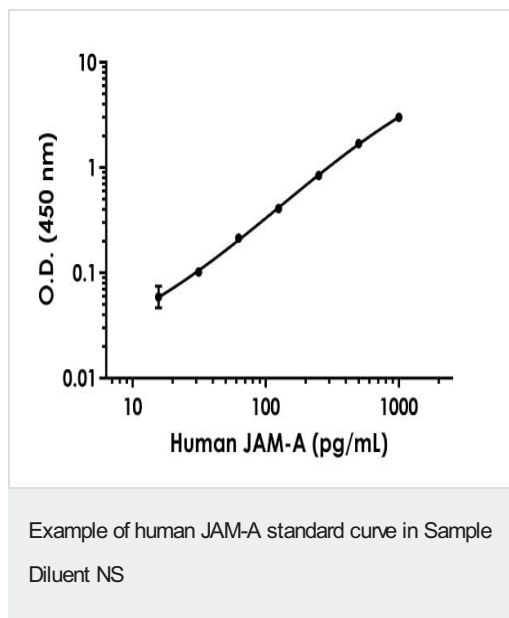
Cellular localization

Cell junction > tight junction. Cell membrane. Localized at tight junctions of both epithelial and endothelial cells.

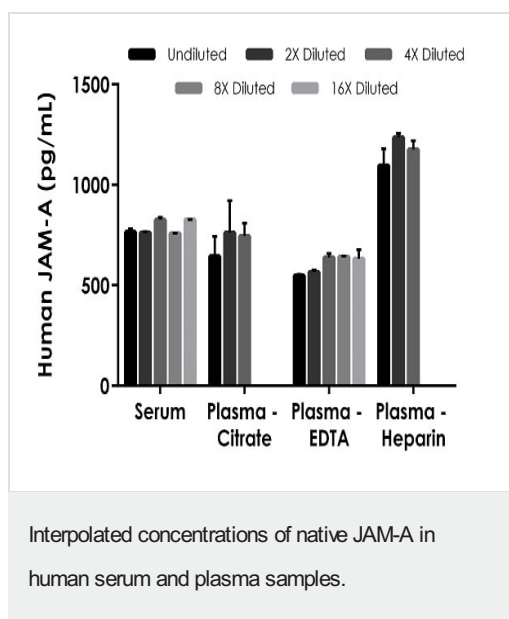
Images



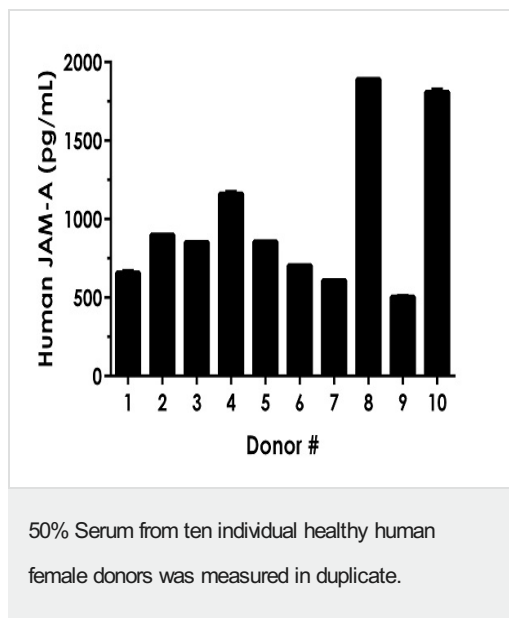
SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



The JAM-A 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 JAM-A were measured in duplicates, interpolated from the JAM-A standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 100%, plasma (citrate) 25%, plasma (EDTA) 100%, and plasma (heparin) 100%. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean JAM-A concentration was determined to be 787.60 pg/mL in serum, 716.49 pg/mL in plasma (citrate), 605.26 pg/mL in plasma (EDTA), and 1072.39 pg/mL in plasma (Heparin).



Interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean JAM-A concentration was determined to be 994.21 pg/mL with a range of 504.21 – 1890.29 pg/mL.

Powered by
recombinant antibodies

 <p>Research with confidence Consistent and reproducible results</p>	 <p>Long-term and scalable supply Recombinant technology</p>
 <p>Success from the first experiment Confirmed specificity</p>	 <p>Ethical standards compliant Animal-free production</p>

Sandwich ELISA - Human JAM-A ELISA Kit
(Junctional Adhesion Molecule)

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