

# Human Junctional Adhesion Molecule 1/JAM-A ELISA Kit ab213800

[1 References](#)   [1 Image](#)

Overview

Product name	Human Junctional Adhesion Molecule 1/JAM-A ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	1	16	81pg/ml	5.83	7.2%
	2	16	325pg/ml	16.9	5.2%
	3	16	1280pg/ml	67.84	5.3%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	1	24	73pg/ml	5.32	7.3%
	2	24	310pg/ml	16.43	5.3%
	3	24	1209pg/ml	93.09	7.7%
Sample type	Cell culture supernatant, Serum				
Assay type	Sandwich (quantitative)				
Sensitivity	< 10 pg/ml				
Range	46.9 pg/ml - 3000 pg/ml				
Assay time	3h 30m				
Assay duration	Multiple steps standard assay				
Species reactivity	<b>Reacts with:</b> Human				
Product overview	The Human Junctional Adhesion Molecule 1/JAM-A Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213800) is designed for the quantitative detection of Human JAM-A in cell culture supernatants and serum.				

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A monoclonal antibody from mouse specific for JAM-A has been pre-coated onto 96-well plates. Standards (Expression system for standard: E.coli; Immunogen sequence: S28-V238) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for JAM-A is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human JAM-A amount of sample captured in plate.

## Notes

Junctional adhesion molecule A/JAM-A is a protein that in humans is encoded by the F11R gene. It is mapped to 1q23.3. This gene is an immunoglobulin-like molecule that co-localizes with tight junctions in endothelium and epithelium and is also found on blood leukocytes and platelets. JAM-A plays an important role in the regulation of tight junction assembly in epithelia. In addition, it can act as a receptor for reovirus, a ligand for the integrin LFA1, involved in leukocyte transmigration and a platelet receptor. JAM-A has a non-redundant role in controlling DC motility, trafficking to lymph nodes, and activation of specific immunity.

## Platform

Pre-coated microplate (12 x 8 well strips)

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

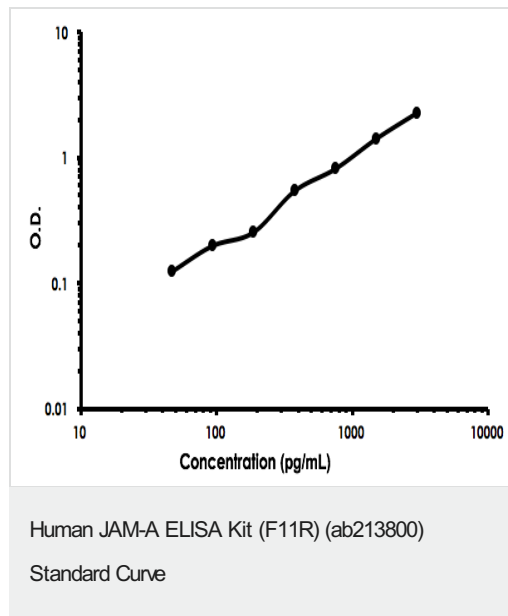
Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human JAM-A coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human JAM-A antibody		1 x 100µl
Lyophilized recombinant Human JAM-A standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

## Function

Seems to plays 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.
<b>Sequence similarities</b>	Belongs to the immunoglobulin superfamily. Contains 2 Ig-like V-type (immunoglobulin-like) domains.
<b>Post-translational modifications</b>	N-glycosylated.
<b>Cellular localization</b>	Cell junction > tight junction. Cell membrane. Localized at tight junctions of both epithelial and endothelial cells.

## Images



Human JAM-A ELISA Kit (F11R) (ab213800) Standard Curve.

**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