

Product datasheet

Human MFGE8 ELISA Kit (Lactadherin) ab213810

1 Image

Overview

Product name Human MFGE8 ELISA Kit (Lactadherin)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	100pg/ml	5.8	= 5.8%
2	16	660pg/ml	36.96	= 5.6%
3	16	1196pg/ml	90.89	= 7.6%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	101pg/ml	8.28	= 8.2%
2	24	717pg/ml	41.58	= 5.8%
3	24	1149pg/ml	102.26	= 8.9%

Sample type Cell culture supernatant, Saliva, Milk, Urine, Serum, Hep Plasma, EDTA Plasma

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 62.5 pg/ml - 4000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview The Human MFGE8 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213810) is designed for the quantitative detection of Human MFGE8 in cell culture supernatants, serum, plasma (heparin, EDTA), saliva, urine and human milk.

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

monoclonal antibody from mouse specific for MFGE8 has been precoated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: L24-C387) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for MFGE8 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 MFGE8 amount of sample captured in plate.

Notes MFGE8 (Milk Fat Globule-Egf Factor 8), also called as Lactadherin or SED1, is a protein which in humans is encoded by the MFGE8 gene. Mfge8 is secreted protein found in vertebrates, including mammals as well as birds. By fluorescence *in situ* hybridization, the MFGE8 gene was mapped to chromosome 15q25. MFGE8 is a factor that links apoptotic cells to phagocytes. It specifically bound to apoptotic cells by recognizing aminophospholipids such as phosphatidylserine. MFGE8, when engaged by phospholipids, bound to cells via its RGD motif. It bound particularly strongly to cells expressing alpha-V-beta-3 integrin. It showed that Mfge8 was expressed in intestinal lamina propria macrophages in mice. Using a wound-healing assay, they showed that Mfge8 promoted migration of intestinal epithelial cells through a PKC-epsilon (PRKCE)-dependent mechanism.

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 MFGE8 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 130µl
Biotinylated anti-Human MFGE8 antibody		1 x 130µl
Lyophilized recombinant Human MFGE8 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

Function Plays an important role in the maintenance of intestinal epithelial homeostasis and the promotion of mucosal healing. Promotes VEGF-dependent neovascularization (By similarity). Contributes to phagocytic removal of apoptotic cells in many tissues. Specific ligand for the alpha-v/beta-3 and

alpha-v/beta-5 receptors. Also binds to phosphatidylserine-enriched cell surfaces in a receptor-independent manner. Zona pellucida-binding protein which may play a role in gamete interaction. Binds specifically to rotavirus and inhibits its replication. Medin is the main constituent of aortic medial amyloid.

Tissue specificity

Mammary epithelial cell surfaces and aortic media. Overexpressed in several carcinomas.

Sequence similarities

Contains 1 EGF-like domain.
Contains 2 F5/8 type C domains.

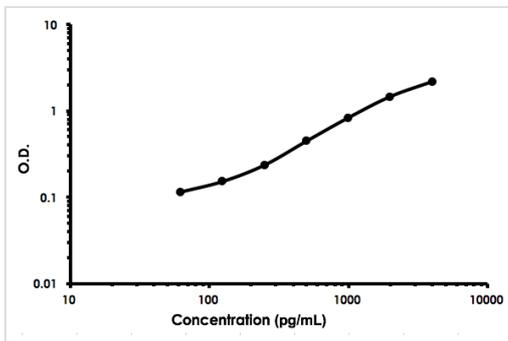
Post-translational modifications

Medin has a ragged N-terminus with minor species starting at Pro-264 and Gly-273.

Cellular localization

Membrane. Secreted.

Images



Human MFGE8 ELISA Kit (Lactadherin) (ab213810) Standard Curve

Human MFGE8 ELISA Kit (Lactadherin) (ab213810) 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