

Mouse GREM2 ELISA kit (PDRC) ab267575

1 Image

Overview

Product name	Mouse GREM2 ELISA kit (PDRC)				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Serum, Plasma, Cell culture media				
Assay type	Sandwich (quantitative)				
Sensitivity	3 pg/ml				
Range	3.28 pg/ml - 800 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		107		85% - 146%
	Plasma		77		68% - 113%
	Cell culture media		94		89% - 102%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	Mouse GREM2 ELISA Kit is designed for the quantitative determination of GREM2 in cell culture supernatants, plasma and serum samples.				

This assay employs an antibody specific for mouse GREM2 coated on a 96-well plate. Standards and samples are pipetted into the wells and GREM2 present in a sample is bound to the wells by

the immobilized antibody. The wells are washed and biotinylated anti-mouse GREM2 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of GREM2 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

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

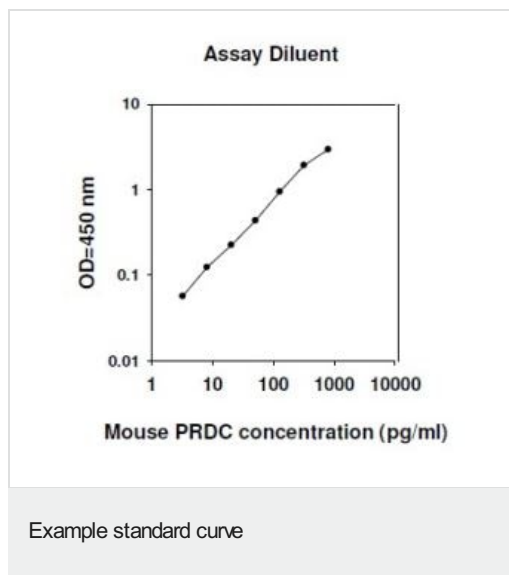
Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Anti-Mouse GREM2 coated Microplate (12 x 8 wells)	1 unit
Assay Diluent C	1 x 30ml
Biotinylated Anti-Mouse GREM2 Detection Antibody	2 vials
Mouse GREM2 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

**Function** Cytokine that inhibits the activity of BMP2 and BMP4 in a dose-dependent manner. Antagonized BMP4-induced suppression of progesterone production in granulosa cells.

**Sequence similarities** Belongs to the DAN family.  
Contains 1 CTCK (C-terminal cystine knot-like) domain.

**Cellular localization** Secreted.

## Images



These standard curves are for demonstration only. A standard curve must be run with each assay.

**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