

Bovine FGF1 ELISA Kit ab273195

1 Image

Overview

Product name	Bovine FGF1 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Cell culture supernatant, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	0.2 ng/ml				
Range	0.205 ng/ml - 50 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		127.4		104% - 126%
	Plasma		124.3		74% - 147%
	Cell culture media		136.9		121% - 147%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Cow				
Product overview	Bovine FGF1 ELISA kit is an <i>in vitro</i> enzyme linked immunosorbent assay for the quantitative measurement of bovine FGF1 in serum (bovine FGF1 concentration is low in normal serum/plasma, and may not be detectable in this assay), plasma and cell culture supernatants.				

This assay employs an antibody specific for bovine FGF1 coated on a 96-well plate. Standards

and samples are pipetted into the wells and FGF1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-bovine FGF1 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 FGF1 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.

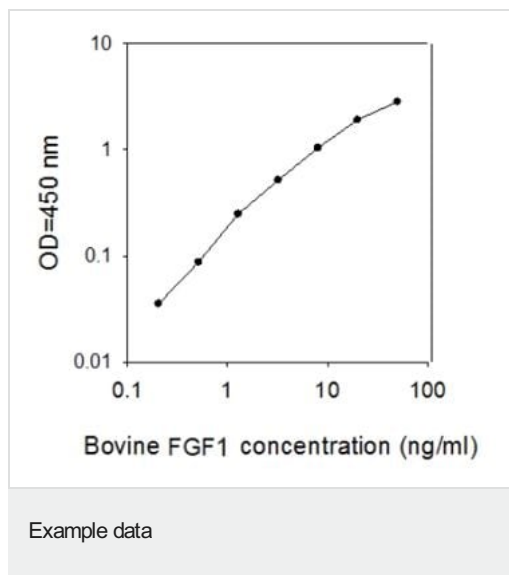
Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
500X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Anti-Bovine FGF1 coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Bovine FGF1 Antibody	2 vials
Bovine FGF1 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

Function The heparin-binding fibroblast growth factors play important roles in the regulation of cell survival, cell division, angiogenesis, cell differentiation and cell migration. They are potent mitogens in vitro.

Sequence similarities Belongs to the heparin-binding growth factors family.

Cellular localization Secreted. Cytoplasm. Cytoplasm > cell cortex. Lacks a cleavable signal sequence. Within the cytoplasm, it is transported to the cell membrane and then secreted by a non-classical pathway that requires Cu(2+) ions and S100A13. Secreted in a complex with SYT1.

Images



This standard curve is 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