

Mouse PLGF-2 ELISA Kit (PGF) ab155459

2 图像

概述				
产品名称	小鼠PLGF-2 ELISA试剂盒(PGF)			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	overall			CV% < 10%
	批次间			
	样品	n	Mean	SD
	Overall			CV% < 12%
样品类型	Cell culture supernatant, Serum, Plasma			
检测类型	Sandwich (quantitative)			
灵敏度	< 1 pg/ml			
范围	1.64 pg/ml - 400 pg/ml			
回收率	特定样本回收率			
	样品类型	平均%		范围
	Cell culture supernatant	110.3		104% - 116%
	Serum	95.3		87% - 103%
	Plasma	91.02		74% - 107%
实验步骤	Multiple steps standard assay			
种属反应性	与反应: Mouse			
产品概述	Abcam's PLGF-2 (PGF) Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay designed for the quantitative measurement of mouse PLGF-2 (PGF) in serum, plasma and cell culture supernatants.			

This assay employs an antibody specific for mouse PLGF-2 coated on a 96-well plate. Standards

and samples are pipetted into the wells and PLGF-2 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse PLGF-2 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 PLGF-2 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**说明** The standard is affinity purified.

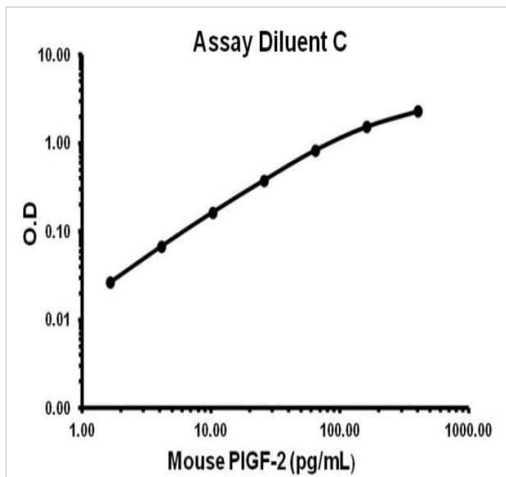
**平台** Microplate

## 性能

**存放说明** Store at -20°C. Please refer to protocols.

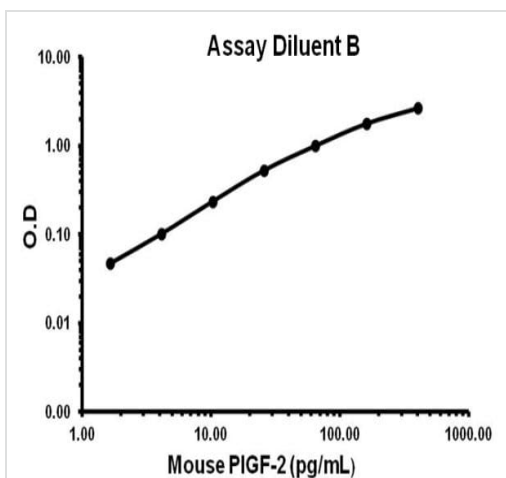
组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent C	1 x 30ml
Biotinylated anti-mouse PLGF-2	2 vials
PLGF-2 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse PLGF-2 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

## 图片



Representative standard curve using ab155459 - Assay Diluent C.

Typical Standard Curve



Representative standard curve using ab155459 - Assay Diluent B.

Typical 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.cn/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors