

Mouse PF4 ELISA Kit (CXCL4) ab100735

★★★★★ 1 Abreviews 2 References 4 Images

Overview

| | |
|------------------|---|
| Product name | Mouse PF4 ELISA Kit (CXCL4) |
| Detection method | Colorimetric |
| Sample type | Cell culture supernatant, Serum, Plasma |
| Assay type | Sandwich (quantitative) |
| Sensitivity | < 6 pg/ml |
| Range | 4.12 pg/ml - 3000 pg/ml |
| Recovery | 100 % |

Sample specific recovery

| Sample type | Average % | Range |
|--------------------------|-----------|-------------|
| Cell culture supernatant | 73.99 | 67% - 83% |
| Serum | 118.2 | 104% - 127% |
| Plasma | 109.3 | 92% - 129% |

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview

Abcam’s PF4 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse PF4 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse PF4 coated on a 96-well plate. Standards and samples are pipetted into the wells and PF4 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse PF4 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 PF4 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Microplate

Properties

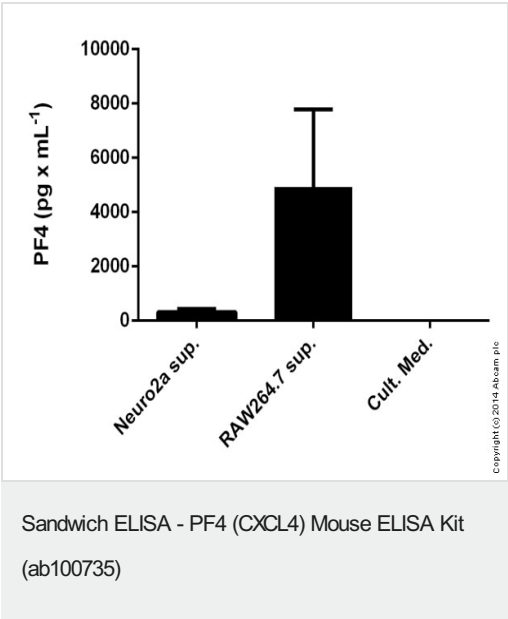
Storage instructions

Store at -20°C. Please refer to protocols.

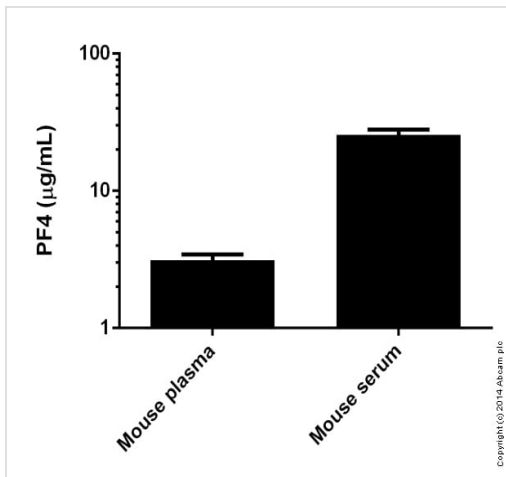
| Components | 1 x 96 tests |
|--|--------------|
| 20X Wash Buffer | 1 x 25ml |
| 5X Assay Diluent B | 1 x 15ml |
| 500X HRP-Streptavidin Concentrate | 1 x 200µl |
| Assay Diluent A | 1 x 30ml |
| Biotinylated anti-mouse PF4 | 2 vials |
| PF4 Microplate (12 x 8 wells) | 1 unit |
| Recombinant Mouse PF4 Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

| | |
|----------------------------------|---|
| Function | Released during platelet aggregation. Neutralizes the anticoagulant effect of heparin because it binds more strongly to heparin than to the chondroitin-4-sulfate chains of the carrier molecule. Chemotactic for neutrophils and monocytes. Inhibits endothelial cell proliferation, the short form is a more potent inhibitor than the longer form. |
| Sequence similarities | Belongs to the intercrine alpha (chemokine CxC) family. |
| Post-translational modifications | Binds non-covalently to a proteoglycan molecule. |
| Cellular localization | Secreted. |

Images

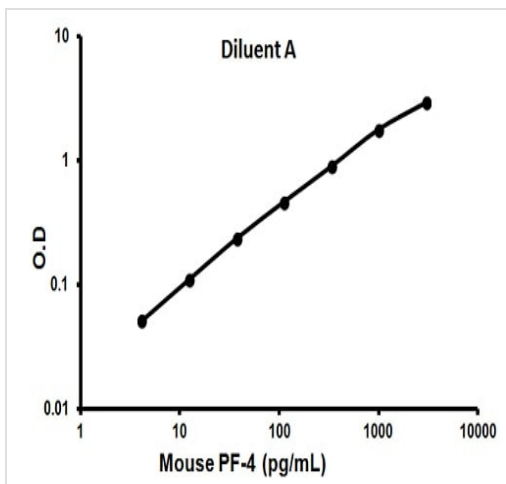


PF4 measured in cell supernatanses showing quantity (pg) per mL tested sample



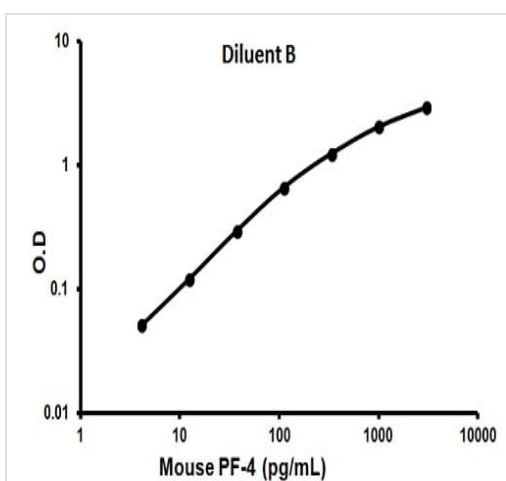
PF4 measured in biofluids showing quantity (microgram) per mL of tested sample

Sandwich ELISA - PF4 (CXCL4) Mouse ELISA Kit (ab100735)



Representative standard curve using ab100735

Typical standard curve



Representative standard curve using ab100735

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.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