

Mouse DPPIV ELISA Kit (CD26) ab155455

1 图像

概述

产品名称

检测方法

精确度

小鼠DPPIV ELISA试剂盒(CD26)

Colorimetric

批次内

样品	n	Mean	SD	CV%
Overall				< 10%

批次间

样品	n	Mean	SD	CV%
Overall				< 12%

样品类型

检测类型

灵敏度

范围

回收率

Cell culture supernatant, Serum, Plasma

Sandwich (quantitative)

< 0.45 ng/ml

0.41 ng/ml - 100 ng/ml

> 68 %

特定样本回收率		
样品类型	平均%	范围
Serum	101.4	94% - 109%
Plasma	88.06	77% - 97%
Cell culture media	77.89	68% - 89%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Mouse
产品概述	Abcam's DPPIV (CD26) Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay designed for the quantitative measurement of mouse DPPIV (CD26) in serum, plasma and cell culture supernatants. This assay employs an antibody specific for DPPIV coated on a 96-well plate. Standards and

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

说明 Optimization may be required with urine samples

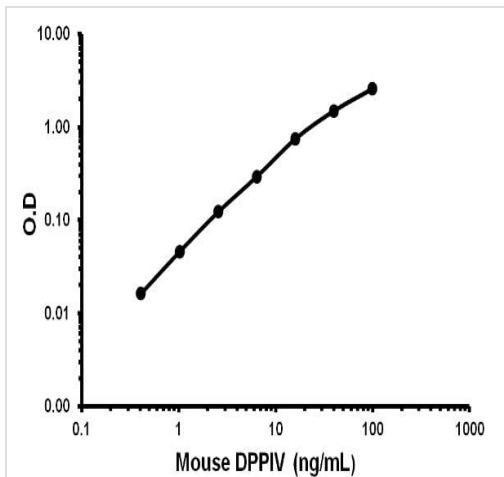
平台 Microplate

性能

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

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

图片



Representative standard curve using ab155455

DPPIV (CD26) Mouse ELISA Kit

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