

Human DKK4 ELISA Kit (Dickkopf-4) ab100503

画像数 1

製品の概要

製品名	Human DKK4 ELISA Kit (Dickkopf-4)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 10 pg/ml
検出範囲	13.72 pg/ml - 10000 pg/ml
添加回収試験	93 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	109.6	97% - 121%
Serum	95.59	85% - 105%
Plasma	82.62	69% - 91%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's DKK4 (Dickkopf-4) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human DKK4 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human DKK4 coated on a 96-well plate. Standards and samples are pipetted into the wells and DKK4 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human DKK4 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 DKK4 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

特記事項 Optimisation may be required with urine samples.

試験プラットフォーム Microplate

製品の特性

保存方法

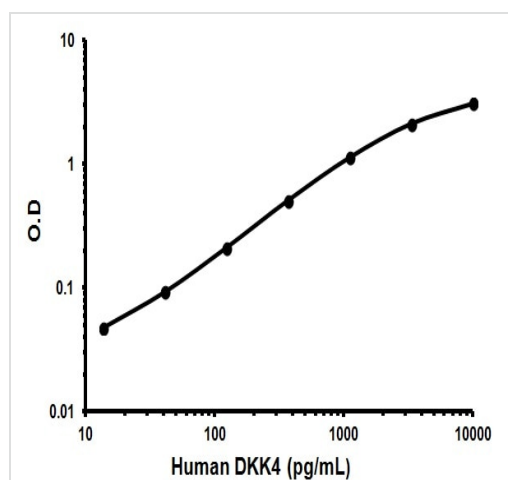
Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Biotinylated anti-Human DKK4	2 vials
DKK4 Microplate (12 x 8 well strips)	1 unit
Recombinant Human DKK4 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

関連性

DKK4, like DKK1, DKK2, and DKK3, possesses an N terminal signal peptide and 2 conserved cysteine-rich domains, which are separated by a linker region and contain 10 cysteine residues each. The second cysteine region has a putative lipid-binding function that may facilitate WNT/DKK interactions at the plasma membrane. The linker region contains 50 to 55 amino acids in DKK1, DKK2, and DKK4, whereas in DKK3 it contains only 12 amino acids. All DKKs have several potential sites for cleavage by furin-type proteases. Northern blot analysis detected no expression of DKK4, but RT-PCR analysis detected DKK4 expression in cerebellum, T cell, esophagus, and lung cDNA libraries. DKK4 blocks Xenopus Wnt8, Wnt3a, and Wnt2b, but not Dsh or Fz8, induction of a secondary axis in frog embryos, indicating that DKKs antagonize WNT function upstream of WNT receptors.

画像



Typical Standard Curve

Representative Standard Curve using ab100503.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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