

Human Cytoglobin ELISA ab277460

画像数 1

製品の概要

製品名	Human Cytoglobin ELISA			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			

サンプル	N	平均値	SD	CV%
Overall				< 10%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall				< 12%

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	0.3 ng/ml
検出範囲	0.287 ng/ml - 70 ng/ml

添加回収試験	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
Serum	90.59	82% - 101%
Plasma	87.98	84% - 94%
Cell culture media	74.86	69% - 87%

ステップ	Multiple steps standard assay
------	-------------------------------

種交差性	交差種: Human
------	------------

製品の概要	Human Cytoglobin ELISA Kit (ab277460) is an <i>in-vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human Cytoglobin in serum, plasma and cell culture supernatants.
-------	--

This assay employs an antibody specific for Human Cytoglobin coated on a 96-well plate.

Standards and samples are pipetted into the wells and Human Cytochrome b5 present in a sample is bound to the wells by the immobilized antibody. The wells are washed, and biotinylated anti-Human Cytochrome b5 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 Human Cytochrome b5 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

製品の特長

保存方法

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

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Biotinylated Anti-Human Cytochrome b5 Antibody	2 vials
Anti-Human Cytochrome b5 coated Microplate	1 unit
Human Cytochrome b5 standard protein (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

機能

May have a protective function during conditions of oxidative stress. May be involved in intracellular oxygen storage or transfer.

組織特異性

Ubiquitously expressed. Highest expression in heart, stomach, bladder and small intestine.

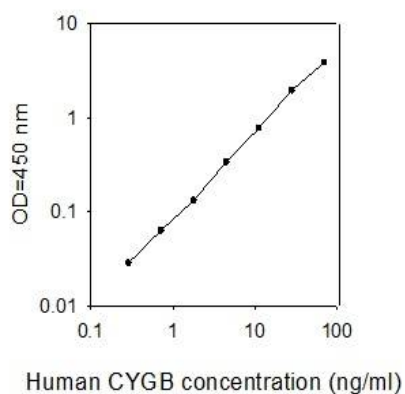
配列類似性

Belongs to the globin family.

細胞内局在

Cytoplasm.

画像



These standard curves are for demonstration only. A standard curve must be run with each assay.

Example data

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