

Rat Ceruloplasmin ELISA Kit (CERP) ab108819

1 图像

概述

产品名称

检测方法

精确度

大鼠Ceruloplasmin ELISA试剂盒(CERP)

Colorimetric

批次内

样品	n	Mean	SD	CV%
Overall				4.8%

批次间

样品	n	Mean	SD	CV%
Overall				8.3%

样品类型

检测类型

灵敏度

范围

回收率

检测时间

实验步骤

种属反应性

产品概述

Cell culture supernatant, Urine, Serum, Plasma, Cell Lysate

Sandwich (quantitative)

= 74 ng/ml

1 ng/ml - 10 ng/ml

95 %

4h 00m

Multiple steps standard assay

与反应: Rat

Abcam's Ceruloplasmin (CERP) rat *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Ceruloplasmin in urine and cell culture supernatants.

A Ceruloplasmin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Ceruloplasmin specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Ceruloplasmin captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

平台 Microplate

性能

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

组件	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Rat Ceruloplasmin Antibody	1 x 120µl
Ceruloplasmin Microplate (12 x 8 well strips)	1 unit
Ceruloplasmin Standard	1 vial
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

功能 Ceruloplasmin is a blue, copper-binding (6-7 atoms per molecule) glycoprotein. It has ferroxidase activity oxidizing Fe(2+) to Fe(3+) without releasing radical oxygen species. It is involved in iron transport across the cell membrane.

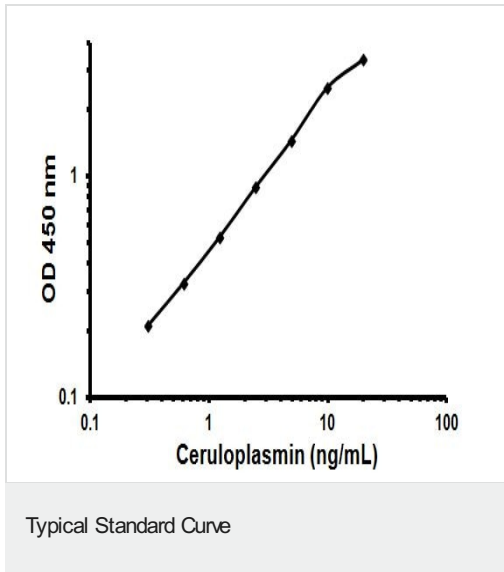
组织特异性 Expressed by the liver and secreted in plasma.

疾病相关 Defects in CP are the cause of aceruloplasminemia (ACERULOP) [MIM:604290]. It is an autosomal recessive disorder of iron metabolism characterized by iron accumulation in the brain as well as visceral organs. Clinical features consist of the triad of retinal degeneration, diabetes mellitus and neurological disturbances.
Note=Ceruloplasmin levels are decreased in Wilson disease, in which copper cannot be incorporated into ceruloplasmin in liver because of defects in the copper-transporting ATPase 2.

序列相似性 Belongs to the multicopper oxidase family.
Contains 3 F5/8 type A domains.
Contains 6 plastocyanin-like domains.

细胞定位 Secreted.

图片



Representative Standard Curve using ab108819.

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