

Human Superoxide Dismutase 1 ELISA Kit ab119520

8 References 1 图像

概述

产品名称 人Superoxide Dismutase 1 ELISA试剂盒

检测方法 Colorimetric

批次内				
样品	n	Mean	SD	CV%
Overall				5.1%

批次间				
样品	n	Mean	SD	CV%
Overall				5.8%

样品类型 Cell culture supernatant, Urine, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 0.04 ng/ml

范围 0.08 ng/ml - 5 ng/ml

回收率 98 %

特定样本回收率		
样品类型	平均%	范围
Serum	98	89% - 108%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's Human Superoxide Dismutase 1 *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Human Cu/Zn Superoxide Dismutase concentrations in cell culture supernatant, serum and plasma (EDTA, citrate, heparin), amniotic fluid, fetal umbilical vein blood and urine.

Cu/Zn Superoxide Dismutase specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a Cu/Zn Superoxide Dismutase HRP-conjugated detection antibody and the microplate is then incubated at room temperature.

After the removal of unbound proteins by washing, TMB is added and catalyzed by HRP to produce a blue color product that changes to yellow after addition of an acidic stop solution. The density of yellow coloration is directly proportional to the Cu/Zn Superoxide Dismutase amount of sample captured in plate.

说明 ab119520 Human Superoxide Dismutase 1 ELISA kit has been shown to detect both SOD1 and SOD3.

平台 Microplate

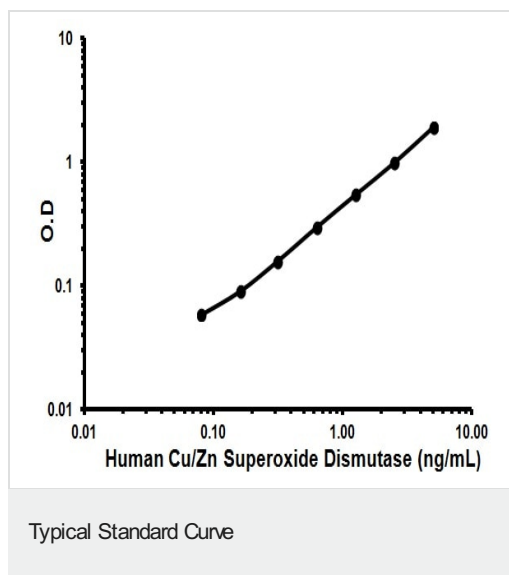
性能

存放说明 Store at +4°C. Please refer to protocols.

组 件	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	2 units
HRP-Conjugate anti-Human Cu/Zn Superoxide Dismutase monoclonal antibody	2 x 20µl
Human Cu/Zn Superoxide Dismutase Standard (5 ng/mL)	2 x 500µl
Microplate coated with anti Human Cu/Zn SOD monoclonal antibody (12 x 8 wells)	1 unit
Phosphate Buffered Saline Concentrate (PBS) (20x)	1 x 5ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
TMB Substrate Solution	1 x 15ml

细胞定位 Cytoplasmic. Note: The pathogenic variants ALS1 Arg-86 and Ala-94 gradually aggregates and accumulates in mitochondria.

图片



Representative Standard Curve using ab119520.

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