

Mouse Beta nerve growth factor ELISA Kit ab277398

1 图像

概述

产品名称 小鼠Beta nerve growth factor ELISA试剂盒

检测方法 Colorimetric

批次内				
样品	n	Mean	SD	CV%
General				< 10%

批次间				
样品	n	Mean	SD	CV%
General				< 12%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 40 pg/ml

范围 0.041 ng/ml - 100 ng/ml

特定样本回收率		
样品类型	平均%	范围
Serum	117.8	81% - 149%
Plasma	108.8	78% - 148%
Cell culture media	108.4	79% - 139%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

预测可用于: Human 

产品概述 Mouse Beta nerve growth factor ELISA Kit (ab277398) is an *in-vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse Beta nerve growth factor in serum (mouse beta-NGF concentration is low in normal serum and plasma samples, and may not be detectable in this assay), plasma and cell culture supernatants.

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

平台Pre-coated microplate (12 x 8 well strips)

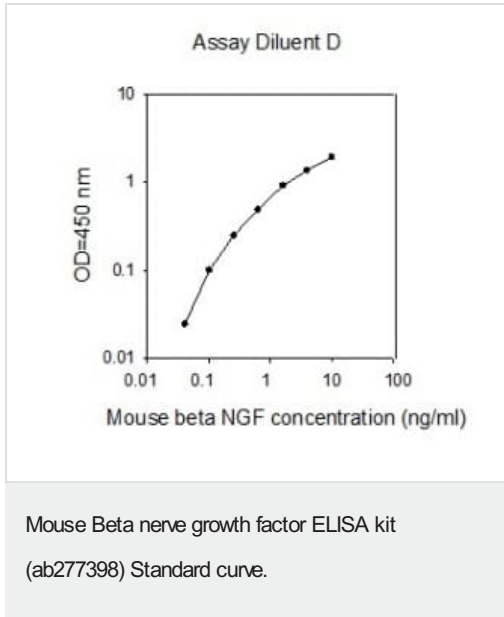
性能

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

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
700X HRP-Streptavidin Concentrate	1 x 200µl
Biotinylated Anti-Mouse Beta nerve growth factor Antibody	2 vials
Anti-Mouse Beta nerve growth factor coated Microplate	1 unit
Mouse Beta nerve growth factor standard protein (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

相关性Nerve growth factor (NGF) is one of a family of neurotrophins that induce the survival and proliferation of neurons. In cell culture, NGF induces the formation of neurite projections and, in vivo, may stimulate the innervation of tissues. NGF plays a role in the repair, regeneration, and protection of neurons, and as such could serve as a therapeutic agent in neurodegenerative conditions such as Alzheimer's disease. NGF enhances survival, growth, neurotransmitter biosynthesis of sympathetic and sensory neurons; neurotrophic factor; cutaneous innervation; growth, differentiation and survival of B lymphocytes. It also has a possible role in allergy and tissue repair. NGF is found in the hypothalamus, pituitary, thyroid gland, testes, epididymis, vascular smooth muscle cells, fibroblasts, mast cells and eosinophils. NGF is upregulated by glutamate, vitamin D3, IL6, FGF basic, astrocyte specific IL1, TNF alpha, PDGF and TGF beta. It is downregulated by GABAergic neuronal activity, glucocorticoids and Schwann cell-specific TGF beta.

细胞定位Secreted.



Typical standard curve – data provided **for demonstration purposes only**. A new standard curve must be generated for each assay performed.

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