

Product datasheet

Human IL-8 ELISA Set (without plates) ab48481

2 References 1 图像

概述

产品名称	人IL-8 ELISA试剂盒(without plates)
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 25 pg/ml
范围	2000 pg/ml - 62.5 pg/ml
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>Abcam's IL-8 Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of IL-8 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>The IL-8 ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>This ELISA set is intended to measure human IL-8 in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>Sensitivity: < 29 pg / ml</p>
经测试应用	适用于: Sandwich ELISA
平台	Reagents

性能

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

组件	标识符	5 x 96 tests	10 x 96 tests	15 x 96 tests
Detection biotinylated anti-IL8 antibody		1 vial	2 vials	3 vials
IL-8 Capture antibody		1 x 100µl	2 x 100µl	3 x 100µl

组件	标识符	5 x 96 tests	10 x 96 tests	15 x 96 tests
IL-8 Standard	Yellow	5 vials	10 vials	15 vials
Ready-to-use TMB		2 vials	4 vials	6 vials
Streptavidin-HRP		1 vial	2 vials	3 vials

相关性 IL-8 is a chemotactic factor that attracts neutrophils, basophils, and T-cells, but not monocytes. It is also involved in neutrophil activation. It is released from several cell types in response to an inflammatory stimulus. IL-8(6-77) has a 5-10-fold higher activity on neutrophil activation, IL-8(5-77) has increased activity on neutrophil activation and IL-8(7-77) has a higher affinity to receptors CXCR1 and CXCR2 as compared to IL-8(1-77), respectively.

细胞定位 Secreted

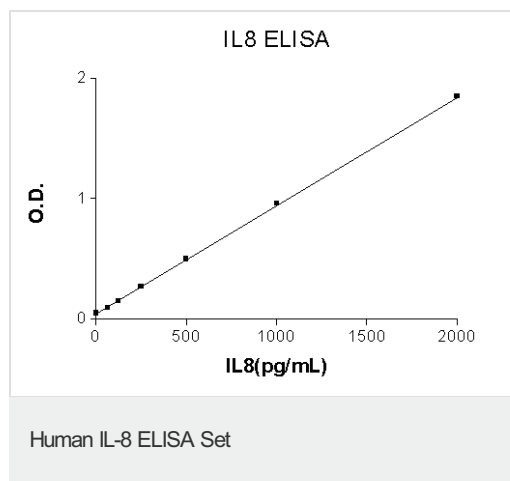
应用

Our [Abpromise guarantee](#) covers the use of **ab48481** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

图片



IL-8 Standard curve ranging from 62.5 to 2000 pg/mL

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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