

Human IL-8 ELISA Kit ab108869

1 References 1 图像

概述

产品名称	人IL-8 ELISA试剂盒																				
检测方法	Colorimetric																				
精确度	<div>批次内</div> <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td></td><td></td><td></td><td>4.1%</td></tr></table> <div>批次间</div> <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td></td><td></td><td></td><td>7.6%</td></tr></table>	样品	n	Mean	SD	CV%	Overall				4.1%	样品	n	Mean	SD	CV%	Overall				7.6%
样品	n	Mean	SD	CV%																	
Overall				4.1%																	
样品	n	Mean	SD	CV%																	
Overall				7.6%																	
样品类型	Cell culture supernatant, Serum, Plasma, Tissue																				
检测类型	Sandwich (quantitative)																				
灵敏度	> 0.01 ng/ml																				
范围	0.016 ng/ml - 1 ng/ml																				
回收率	97 %																				
检测时间	5h 00m																				
实验步骤	Multiple steps standard assay																				
种属反应性	与反应: Human																				
产品概述	Abcam's IL-8 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for																				

An IL-8 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an IL-8 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 IL-8 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 Human IL-8 Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
IL-8 Microplate (12 x 8 well strips)	1 unit
IL-8 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

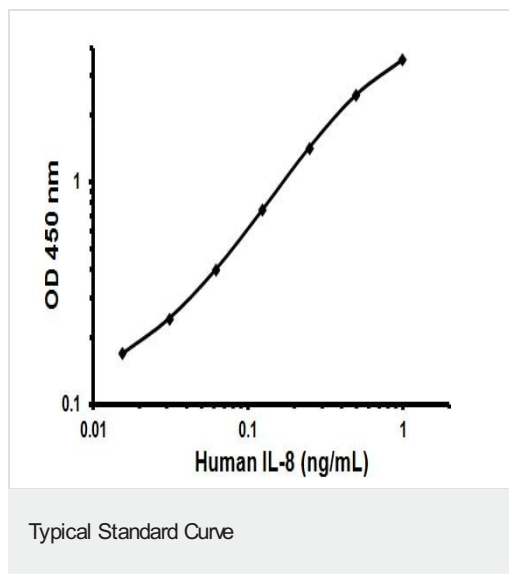
功能 IL-8 is a chemotactic factor that attracts neutrophils, basophils, and T-cells, but not monocytes. It is also involved in neutrophil activation. It is released from several cell types in response to an inflammatory stimulus. IL-8(6-77) has a 5-10-fold higher activity on neutrophil activation, IL-8(5-77) 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.

序列相似性 Belongs to the intercrine alpha (chemokine CxC) family.

翻译后修饰 Several N-terminal processed forms are produced by proteolytic cleavage after secretion from at least peripheral blood monocytes, leukocytes and endothelial cells. In general, IL-8(1-77) is referred to as interleukin-8. IL-8(6-77) is the most prominent form.

细胞定位 Secreted.

图片



Representative Standard Curve using ab108869.

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