

Human Chemerin ELISA Kit ab155430

★★★★★ 1 Abreviews 6 References 画像数 1

製品の概要

製品名	Human Chemerin ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	overall			CV%
				< 10%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	overall			CV%
				< 12%
サンプルの種類	Serum, Plasma, Cell culture media			
アッセイタイプ	Sandwich (quantitative)			
検出感度	< 0.5 ng/ml			
検出範囲	0.51 ng/ml - 50 ng/ml			
添加回収試験	特定サンプルでの回収試験			

サンプルの種類	平均 %	測定範囲
Serum	= 126.3	107% - 145%
Plasma	= 134.6	105% - 144%
Cell culture media	= 136.9	120% - 146%

ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<p>Abcam's Chemerin Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay designed for the quantitative measurement of Human Chemerin in serum, plasma and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human Chemerin coated on a 96-well plate. Standards and samples are pipetted into the wells and Chemerin present in a sample is bound to</p>

the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Chemerin 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 Chemerin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

#### 特記事項

Optimisation may be required with urine samples.

#### 試験プラットフォーム

Microplate

#### 製品の特性

#### 保存方法

Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
700X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent C	1 x 30ml
Biotinylated anti-Human Chemerin	2 vials
Chemerin Microplate (12 x 8 wells)	1 unit
Recombinant Human Chemerin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

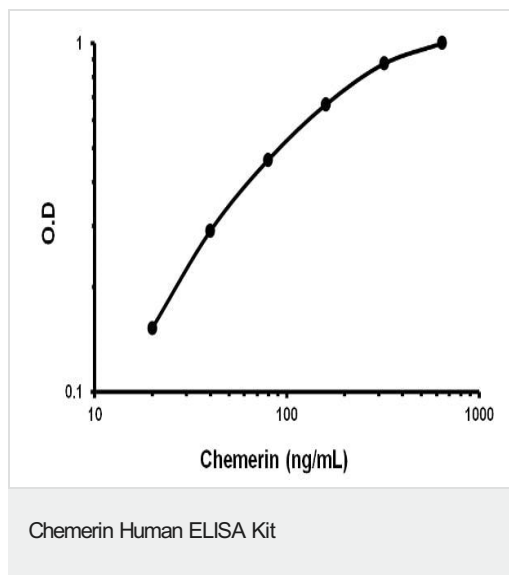
#### 関連性

This gene encodes a secreted chemotactic protein that initiates chemotaxis via the ChemR23 G protein-coupled seven-transmembrane domain ligand. Expression of this gene is upregulated by the synthetic retinoid tazarotene and occurs in a wide variety of tissues. The active protein has several roles, including that as an adipokine, and is truncated on both termini from the proprotein.

#### 細胞内局在

Secreted

#### 画像



Representative standard curve using ab155430

**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.co.jp/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors