

Rat TNF alpha ELISA Kit ab46070

★★★★★ 1 Abreviews 96 References 画像数 3

製品の概要

製品名	Rat TNF alpha 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
	A	12	543.5pg/ml	44.7
	B	10	130.1pg/ml	10.3
			CV%	
			8.2%	
			7.9%	
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	A	9	541.4pg/ml	33.2
	B	9	149.4pg/ml	9.8
			CV%	
			6.1%	
			6.5%	
サンプルの種類	Cell culture supernatant, Serum			
アッセイタイプ	Sandwich (quantitative)			
検出感度	< 15 pg/ml			
検出範囲	31.25 pg/ml - 1000 pg/ml			
全工程の試験時間	3h 45m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Rat			
製品の概要	Abcam's TNF alpha Rat <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Rat TNF alpha in cell culture supernatants, serum and plasma samples.			

A monoclonal antibody specific for TNF alpha has been coated onto the wells of the microtiter strips provided. Samples, including standards of known TNF alpha concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for TNF alpha are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is

added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of TNF alpha present in the samples.

This kit will recognize both endogenous and recombinant rat TNF alpha

## 試験プラットフォーム

Microplate

## 製品の特性

### 保存方法

Store at +4°C. Please refer to protocols.

内容	ラベル	2 x 96 tests	1 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7.5ml	1 x 7.5ml
Biotinylated anti-TNF alpha	Red	2 x 400µl	1 x 400µl	1 x 400µl
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml	1 x 11ml
HRP Diluent	Red	1 x 23ml	1 x 12ml	1 x 12ml
Plastic Plate Covers		4 units	2 units	2 units
Stop Reagent	Black	2 x 11ml	1 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl	2 x 5µl
TNF alpha Microplate (12 x 8 well strips)		2 units	1 unit	1 unit
TNF alpha Standard (Lyophilized)	Yellow	4 vials	2 vials	2 vials

### 機能

Cytokine that binds to TNFRSF1A/TNFR1 and TNFRSF1B/TNFR. It is mainly secreted by macrophages and can induce cell death of certain tumor cell lines. It is potent pyrogen causing fever by direct action or by stimulation of interleukin-1 secretion and is implicated in the induction of cachexia. Under certain conditions it can stimulate cell proliferation and induce cell differentiation.

### 関連疾患

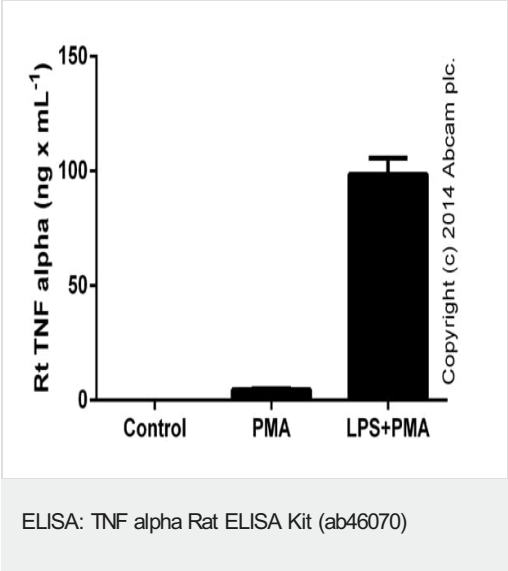
Genetic variations in TNF are a cause of susceptibility psoriatic arthritis (PSORAS) [MIM:607507]. PSORAS is an inflammatory, seronegative arthritis associated with psoriasis. It is a heterogeneous disorder ranging from a mild, non-destructive disease to a severe, progressive, erosive arthropathy. Five types of psoriatic arthritis have been defined: asymmetrical oligoarthritis characterized by primary involvement of the small joints of the fingers or toes; asymmetrical arthritis which involves the joints of the extremities; symmetrical polyarthritis characterized by a rheumatoidlike pattern that can involve hands, wrists, ankles, and feet; arthritis mutilans, which is a rare but deforming and destructive condition; arthritis of the sacroiliac joints and spine (psoriatic spondylitis).

### 配列類似性

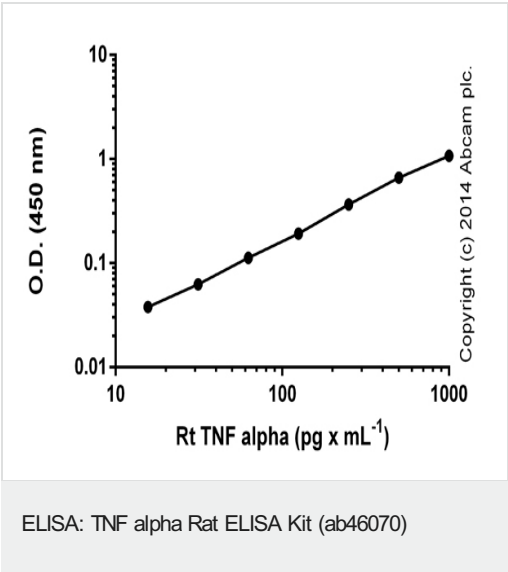
Belongs to the tumor necrosis factor family.

翻訳後修飾	The soluble form derives from the membrane form by proteolytic processing.
	The membrane form, but not the soluble form, is phosphorylated on serine residues.
	Dephosphorylation of the membrane form occurs by binding to soluble TNFRSF1A/TNFR1.
	O-glycosylated; glycans contain galactose, N-acetylgalactosamine and N-acetylneuraminic acid.
細胞内局在	Secreted and Cell membrane.

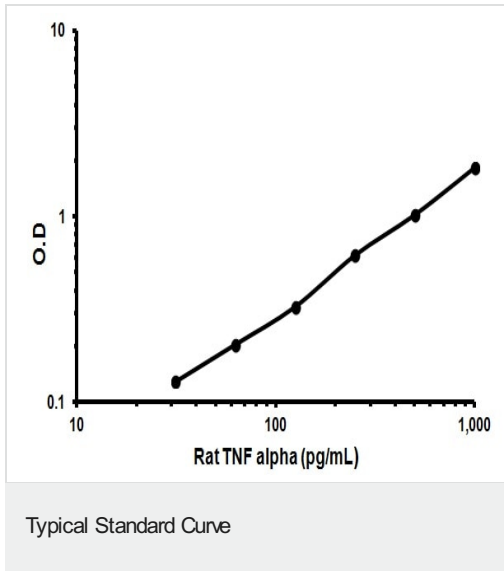
画像



Rat TNF alpha measured in culture supernatants (dilution range 1/1-1/200) from control NR8383 cells or cells stimulated for 24 hours with 50 ng x mL<sup>-1</sup> of PMA ((PMA) [ab120297](#)) or PMA with the addition of 1 ug x mL<sup>-1</sup> of LPS (Sigma) (LPS+PMA) for the last 6 hours (duplicates +/- SD).



Standard curve with background signal subtracted (duplicates; +/- SD).



Representative Standard Curve using ab46070

**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