

画像数 1

製品の概要

製品名	Human Interferon gamma ELISA Set (without plates)
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 5 pg/ml
検出範囲	400 pg/ml - 12.5 pg/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<p>Abcam's IFNγ Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of IFNγ in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>The Interferon gamma 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 Interferon gamma 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: < 5 pg / ml</p>
アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Reagents

製品特性

保存方法		Store at +4°C. Please refer to protocols.		
内 容	ラベル	5 x 96 tests	10 x 96 tests	15 x 96 tests
Detection biotinylated anti-IFN gamma antibody		1 x 0.55ml	2 x 0.55ml	3 x 0.55ml

内 容	ラベル	5 x 96 tests	10 x 96 tests	15 x 96 tests
IFN gamma Capture antibody		1 x 0.5ml	2 x 0.5ml	3 x 0.5ml
IFN gamma Standard (Lyophilized)	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

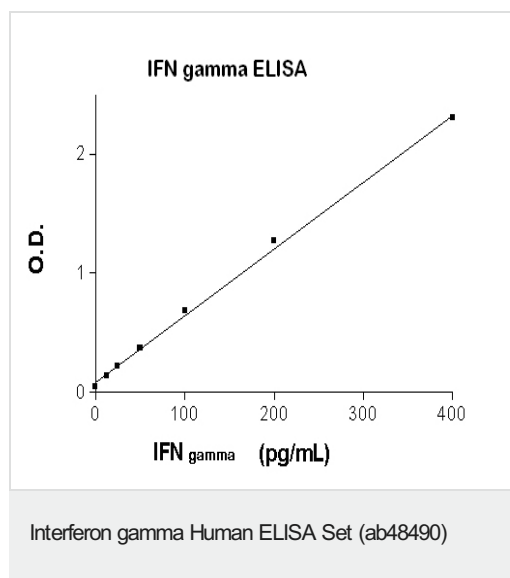
機能	Produced by lymphocytes activated by specific antigens or mitogens. IFN-gamma, in addition to having antiviral activity, has important immunoregulatory functions. It is a potent activator of macrophages, it has antiproliferative effects on transformed cells and it can potentiate the antiviral and antitumor effects of the type I interferons.
組織特異性	Released primarily from activated T lymphocytes.
関連疾患	In Caucasians, genetic variation in IFNG is associated with the risk of aplastic anemia (AA) [MIM:609135]. AA is a rare disease in which the reduction of the circulating blood cells results from damage to the stem cell pool in bone marrow. In most patients, the stem cell lesion is caused by an autoimmune attack. T-lymphocytes, activated by an endogenous or exogenous, and most often unknown antigenic stimulus, secrete cytokines, including IFN-gamma, which would in turn be able to suppress hematopoiesis.
配列類似性	Belongs to the type II (or gamma) interferon family.
翻訳後修飾	Proteolytic processing produces C-terminal heterogeneity, with proteins ending alternatively at Gly-150, Met-157 or Gly-161.
細胞内局在	Secreted.

アプリケーション

The Abpromise guarantee Abpromise保証は、次のテスト済みアプリケーションにおけるab48490の使用に適用されます
 アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



IFNgamma Standard curve ranging from 12.5 to 400 pg/mL

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