

Extra Standard for Interferon gamma ELISA Kit ab59555

製品の概要

製品名	Extra Standard for Interferon gamma ELISA Kit
アッセイタイプ	Sandwich
製品の概要	Extra Standard for Interferon gamma ELISA Kit (ab59555) is part of the Human IFN gamma ELISA Kit (ab46025).
	Refer to the kit protocol booklet for instructions of use. Please contact our Scientific Support team for more lot specific information.

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	1 units
IFN gamma Standard (Lyophilized)	Yellow	1 vial

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.

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