

Human Lysozyme ELISA Kit (LZM) ab108877

[1 References](#) [画像数 1](#)

製品の概要

製品名 Human Lysozyme ELISA Kit (LZM)

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Overall				4.5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall				8.7%

サンプルの種類 Milk

アッセイタイプ Competitive

検出感度 = 0.09 µg/ml

検出範囲 0.093 µg/ml - 6 µg/ml

添加回収試験 = 98 %

全工程の試験時間 3h 00m

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's Lysozyme (LZM) Human *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Human Lysozyme in milk.

A Lysozyme specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated Lysozyme is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex 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 inversely proportional to the amount of Lysozyme captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

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
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	1 x 30ml
Biotinylated Human Lysozyme (Lyophilized)	1 vial
Chromogen Substrate	1 x 7ml
Lysozyme Microplate (12 x 8 well strips)	1 unit
Lysozyme Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

機能

Lysozymes have primarily a bacteriolytic function; those in tissues and body fluids are associated with the monocyte-macrophage system and enhance the activity of immunoagents.

関連疾患

Amyloidosis 8

配列類似性

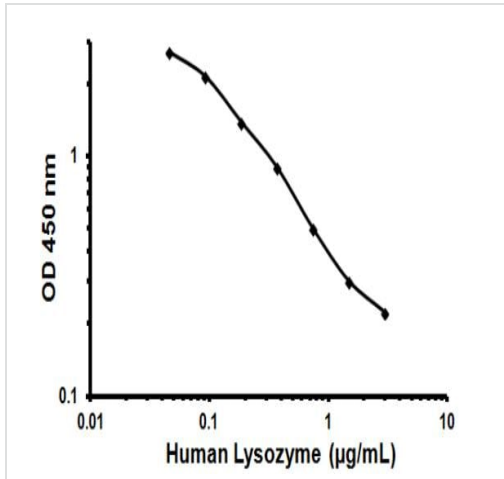
Belongs to the glycosyl hydrolase 22 family.

細胞内局在

Secreted.

画像

Representative Standard Curve using ab108877.



Typical Standard Curve

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