

Product datasheet

Human LAG3 ELISA Kit ab193707

1 Image

Overview

Product name Human LAG3 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall < 10%

Inter-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall < 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 45 pg/ml

Range 45 pg/well - 12000 pg/well

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	78.4	70% - 86%
Plasma	78.2	68% - 95%
Cell culture media	76.83	69% - 85%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's LAG3 (lymphocyte activation gene 3) Human ELISA Kit (ab193707) is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human LAG3 in serum, plasma and cell culture supernatant. (UniprotID: P18627)

This assay employs an antibody specific for human LAG3 coated on a 96-well plate. Standards and samples are pipetted into the wells and the immobilized antibody captures LAG3 present in the samples. The wells are washed and biotinylated anti-human LAG3 is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of LAG3 bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

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

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent D	1 x 15ml
5X Assay Diluent B	1 x 15ml
Biotinylated Human LAG3 detection antibody (lyophilized)	2 vials
Human LAG3 Standards (lyophilized)	2 vials
Pre-coated Human LAG3 Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

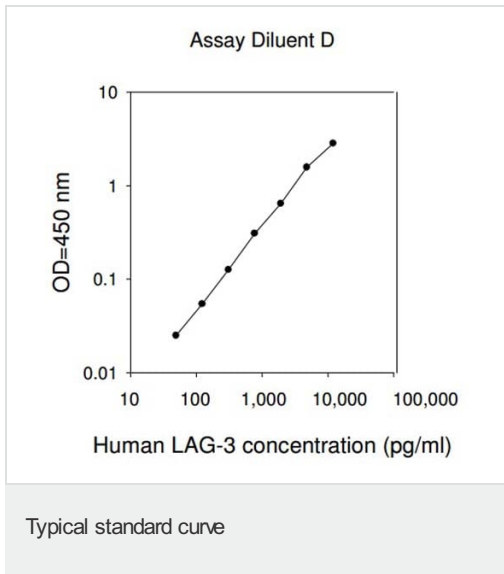
Function Involved in lymphocyte activation. Binds to HLA class-II antigens.

Tissue specificity On cell surface of activated NK and T-lymphocytes.

Sequence similarities Contains 3 Ig-like C2-type (immunoglobulin-like) domains.
Contains 1 Ig-like V-type (immunoglobulin-like) domain.

Cellular localization Membrane.

Images



Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

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.com/abpromise> or contact our technical team.

Terms and conditions

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