

Product datasheet

Giardia lamblia antigen ELISA Kit ab178643

Overview

Product name Giardia lamblia antigen ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	12			6.86%
2	12			6.41%
3	12			6.36%
4	12			5.04%

Inter-assay

Sample	n	Mean	SD	CV%
1	11			4.77%
2	11			4.64%
3	11			7.77%
4	11			10.9%

Sample type Other biological fluids

Assay type Indirect

Sensitivity = 97.5 %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Giardia lamblia in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of Giardia lamblia antigens in faeces.

A 96-well plate has been precoated with polyclonal anti-Giardia lamblia antibodies. Controls or

test samples are added to the wells and incubated; Giardia antigens present in the stool supernatant are captured. Following washing, a horseradish peroxidase (HRP) labelled anti-Giardia antibody is added to the wells, which binds to the immobilized antibodies. TMB is then catalyzed by the HRP to produce a blue color product that changes to yellow after adding an acidic stop solution. The density of yellow coloration is directly proportional to the amount of Giardia lamblia antigen sample captured in plate.

Tested applications

Suitable for: Indirect ELISA

Platform

Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
10X Wash Solution	1 x 100ml
Anti-Giardia lamblia Coated Microplate (12 x 8 well strips)	1 unit
Anti-Giardia lamblia Conjugate	1 x 12ml
Cover Foil	1 unit
Giardia lamblia Antigen Negative Control	1 x 2ml
Giardia lamblia Antigen Positive Control	1 x 2ml
Sample Diluent	1 x 100ml
Stop Solution	1 x 15ml
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

Applications

Our [Abpromise guarantee](#) covers the use of **ab178643** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Indirect ELISA		Use at an assay dependent concentration.

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