

Product datasheet

Pig IgG ELISA Kit ab291065

1 Image

Overview

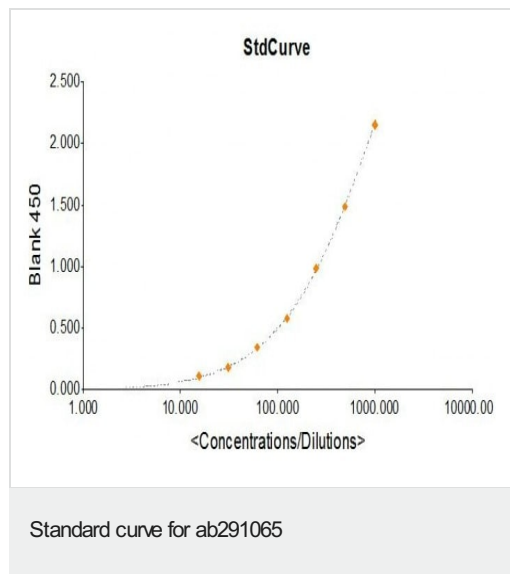
Product name	Pig IgG ELISA Kit
Detection method	Colorimetric/Fluorometric
Sample type	Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	= 7.08 ng/ml
Range	15.625 ng/ml - 1000 ng/ml
Assay time	0h 70m
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Pig
Product overview	Pig IgG ELISA Kit (ab291065) is a highly sensitive two-site enzyme linked immunoassay (ELISA) for measuring IgG in pig biological samples.
Platform	Microplate (12 x 8 well strips)

Properties

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

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Calibrator	1 vial
Chromogen Substrate Solution	1 x 12ml
ELISA Micro Plate, antibody coated	1 unit
Enzyme Conjugated Detection Antibody	1 x 150µl
Stop Solution	1 x 12ml

Images



Typical Standard Curve: These standard curves are for demonstration only. A standard curve must be run with each assay.

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