

Mouse beta IGH3 ELISA Kit ab193717

1 Image

Overview

Product name	Mouse beta IGH3 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	3 pg/ml				
Range	3 pg/ml - 1000 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		140.1		132% - 148%
	Plasma		136		115% - 150%
	Cell culture media		143.6		137% - 149%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	Abcam's beta IGH3 mouse ELISA Kit (ab193717) is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of mouse BIGH3 in serum, plasma and cell culture supernatants.				

This assay employs an antibody specific for mouse beta IGH3 coated on a 96-well plate.

Standards and samples are pipetted into the wells and the immobilized antibody captures beta IGH3 present in the samples. The wells are washed and biotinylated anti-mouse beta IGH3 antibody 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 beta IGH3 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.

Get better reproducibility in only 90 minutes with Mouse beta IG H3 ELISA Kit ([ab206987](#)) from our SimpleStep ELISA® range.

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

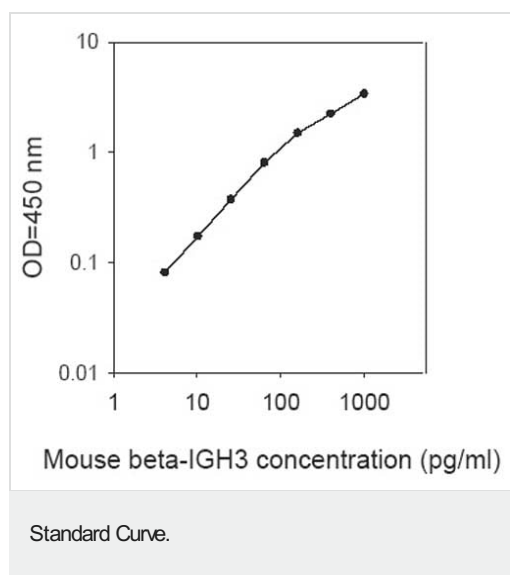
Properties

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated Mouse beta-IGH3 detection antibody	2 vials
Mouse beta-IGH3 standards (lyophilized)	2 vials
Pre-coated mouse beta-IGH3 Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Cellular localization Secreted, extracellular space, extracellular matrix.

Images



Typical Standard Curve using ab193717 BIGH3 Mouse ELISA Kit.

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