

Mouse Sonic Hedgehog-N ELISA Kit ab155450

1 Image

Overview

Product name Mouse Sonic Hedgehog-N 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 Lysate, Tissue Lysate

Assay type Sandwich (quantitative)

Sensitivity < 5 pg/ml

Range 5 pg/ml - 1200 pg/ml

Recovery	Sample specific recovery		
	Sample type	Average %	Range
	Cell Lysate	126.2	114% - 137%
	Tissue Lysate	130.3	114% - 150%

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Mouse

Product overview Abcam’s Shh-N Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of mouse Shh-N in cell lysates and tissue lysates.

This assay employs an antibody specific for mouse Shh-N coated on a 96-well plate. Standards and samples are pipetted into the wells and Shh-N present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse Shh-N antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells

and color develops in proportion to the amount of Shh-N bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

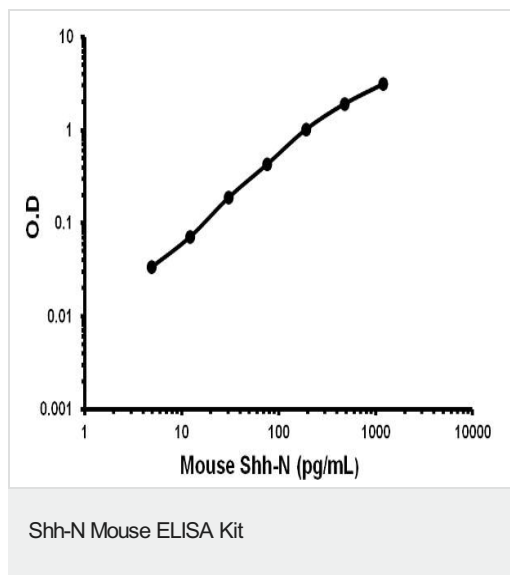
Notes	Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.
Platform	Microplate

Properties

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
2X Cell Lysis Buffer	1 x 5ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
5X Sample Diluent Buffer	1 x 10ml
Biotinylated anti-mouse Shh-N	2 vials
Recombinant Mouse Shh-N Standard (lyophilized)	2 vials
Shh-N Microplate (12 x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Images



Representative standard curve using ab155450

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