

Human SOST ELISA Kit ab155440

1 References 1 Image

Overview

Product name	Human SOST 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	< 15 pg/ml				
Range	0.017 ng/ml - 10 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		123.3		110% - 136%
	Plasma		125.7		118% - 134%
	Cell culture media		96.04		86% - 105%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Abcam's SOST Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human SOST in serum, plasma and cell culture supernatants.				

the immobilized antibody. The wells are washed and biotinylated anti-Human SOST 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 SOST bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples.

Platform Microplate

Properties

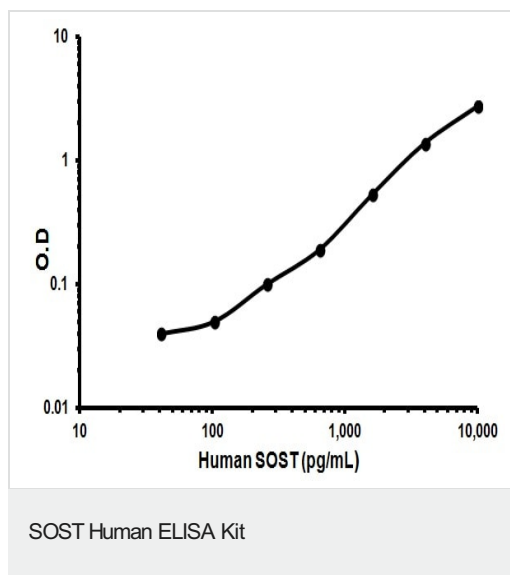
Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Biotinylated anti-Human SOST	2 vials
Recombinant Human SOST Standard (lyophilized)	2 vials
SOST Microplate (12 x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Relevance Negative regulator of bone growth that acts through inhibition of Wnt signaling and bone formation.

Cellular localization Secreted; extracellular space; extracellular matrix

Images



Representative standard curve using ab155440

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