

Human Heparin ELISA Kit ab285267

1 Image

Overview

Product name	Human Heparin ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma, Tissue Homogenate
Assay type	Sandwich (quantitative)
Range	9.3 ng/ml - 300 ng/ml
Assay time	0h 90m
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	This Heparin ELISA kit ab285267 (previously called QuickDetect Heparin (Human) ELISA Kit, E4464) is a sandwich ELISA assay for measuring the quantity of human Heparin in serum, plasma and cell culture and other biological fluid. The entire process can be performed within 90 minutes with only a single incubation and a wash step respectively, resulting in fewer handling steps, fewer errors and more consistent results.
Notes	This product is manufactured by BioVision, an Abcam company and was previously called E4464 QuickDetect™ Heparin (Human) ELISA Kit. E4464-100 is the same size as the 96 test size of ab285267.
Platform	Microplate (12 x 8 well strips)

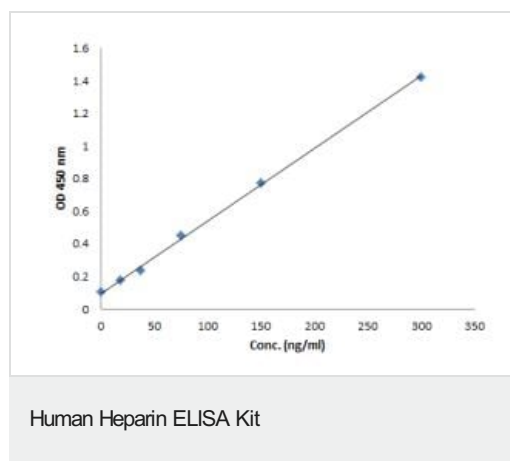
Properties

Storage instructionsStore at +4°C. Please refer to protocols.

Components	96 tests	96 tests
20X Wash buffer	1 x 25ml	1 x 25ml
Chromogen Solution A	1 x 6ml	1 x 6ml
Chromogen Solution B	1 x 6ml	1 x 6ml
HRP-Conjugate Reagent	1 x 6ml	1 x 6ml

Components	96 tests	96 tests
Micro ELISA Plate	1 unit	1 unit
Microplate sealers	2 units	2 units
Special Diluent	1 x 6ml	1 x 6ml
Standard (S0-S5)	1 x 0.6ml	1 x 0.6ml
Standard Diluent	1 x 6ml	1 x 6ml
Stop Solution	1 x 6ml	1 x 6ml

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