

Product datasheet

Pre-coated 384 well Microplate SimpleStep ELISA® ab203359

SimpleStep ELISA

★★★★★ [1 Abreviews](#) [1 References](#) [3 Images](#)

Overview

Product name Pre-coated 384 well Microplate SimpleStep ELISA®

Product overview Abcam's 384-Well SimpleStep ELISA® Microplate can be used in place of the 96 well microplate that accompanies any of our 1x96well SimpleStep ELISA® kit from the catalogue.

To fully utilize all 384-wells, users need to combine reagents from two 1 x 96-well SimpleStep ELISA kits and it is recommended that an automated liquid handling system is used with the 384-well microplate.

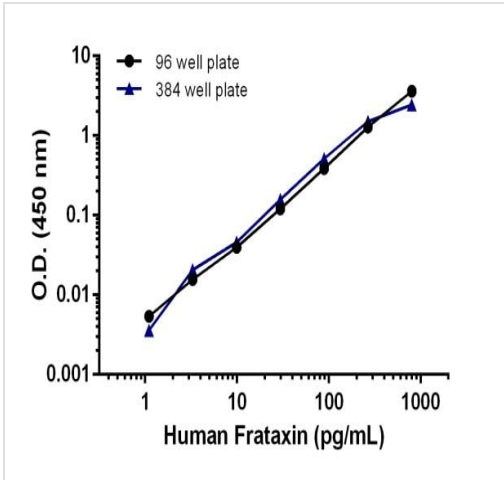
The microplate is pre-coated with the monoclonal antibody of a proprietary tag and is only suitable for use with the our SimpleStep ELISA range.

Properties

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

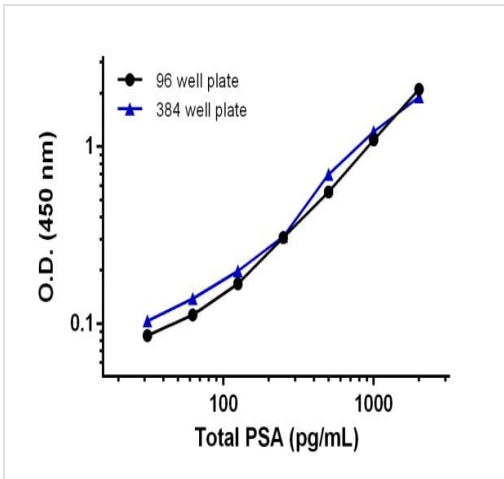
Components	1 units
Pre-Coated 384-Well Microplate	1 x 384 units

Images



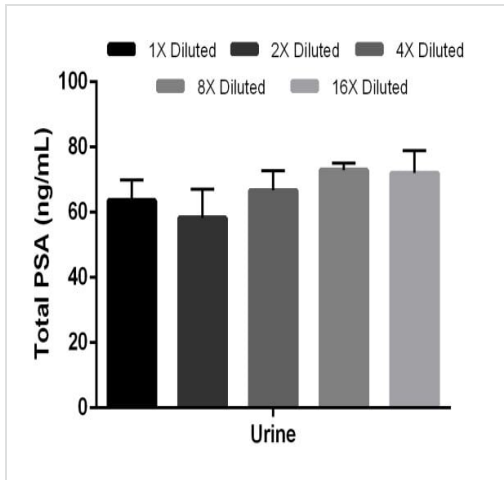
Comparison of human Frataxin protein standard curve in the 96 well microplate vs 384 well microplate format.

The protein standard from the 96 well microplate human Frataxin SimpleStep ELISA kit ([ab176112](#)) was ran on the 384 well microplate. Signal achieved on the 384 well microplate is almost identical that achieved in the 96 well format.



Comparison of the human total PSA protein standard curve in the 96 well microplate vs 384 well microplate format.

The protein standard from the 96 well microplate human total PSA SimpleStep ELISA kit ([ab188388](#)) was ran on the 384 well microplate. Signal achieved on the 384 well microplate is almost identical that achieved in the 96 well format.



Total PSA levels were measured in human urine sample, using reagents from the total PSA Human SimpleStep ELISA Kit ([ab188388](#)) on the 384 well microplate.

Linearity of dilution of human urine samples measuring total PSA ran in the 384 well microplate.

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