

Product datasheet

Human Thyroid Stimulating Hormone ELISA Kit (TSH) ab108659

[1 Abreviews](#) [1 Image](#)

Overview

Product name	Human Thyroid Stimulating Hormone ELISA Kit (TSH)
Detection method	Colorimetric
Sample type	Serum
Assay type	Sandwich (quantitative)
Sensitivity	= 0.05 U/ml
Range	0.05 µIU/ml - 10 µIU/ml
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	<p>Abcam's Thyroid Stimulating Hormone <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Thyroid Stimulating Hormone in Human serum.</p> <p>A 96-well plate has been precoated with anti-Thyroid Stimulating Hormone antibodies. Samples and standards are added to the wells, where Thyroid Stimulating Hormone in the sample and standards binds to the precoated antibody. After incubation and washing, added Anti-Thyroid Stimulating Hormone HRP conjugate binds to the antibody-Thyroid Stimulating Hormone complex. After incubation, the wells are washed to remove unbound material and TMB substrate is then added which is catalyzed by HRP to produce blue coloration. The reaction is terminated by addition of Stop Solution which stops the color development and produces a color change from blue to yellow. The intensity of signal is directly proportional to the amount of Thyroid Stimulating Hormone in the sample and the intensity is measured at 450 nm.</p>
Tested applications	Suitable for: Sandwich ELISA
Platform	Microplate

Properties

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

Components	1 x 96 tests
Anti-Thyroid Stimulating Hormone Coated Microplate (12 x 8 wells)	1 unit

Components	1 x 96 tests
Anti-Thyroid Stimulating Hormone HRP Conjugate Reagent	1 x 13ml
Stop Solution	1 x 11ml
TMB Reagent	1 x 11ml
TSH Standard 2 - 0.5 μ IU/mL (Lyophilized)	1 vial
TSH Standard 3 - 2.0 μ IU/mL (Lyophilized)	1 vial
TSH Standard 4 - 5 μ IU/mL (Lyophilized)	1 vial
TSH Standard 5 - 10 μ IU/mL (Lyophilized)	1 vial
TSH Standard 6 - 25 μ IU/mL (Lyophilized)	1 vial
TSH Standard 1 - 0 μ IU/mL (Lyophilized)	1 vial

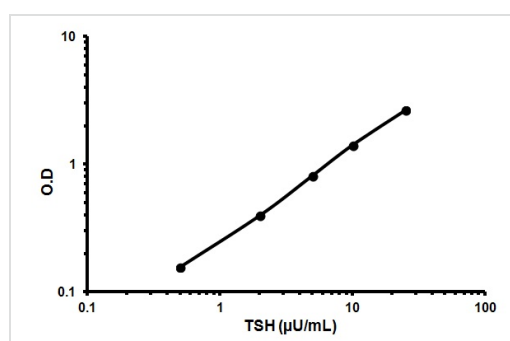
Applications

Our [Abpromise guarantee](#) covers the use of **ab108659** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



Representative standard curve using
ab108659

Typical Standard Curve

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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