

Thyroxine (Total T4) ELISA Kit ab178664

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

Overview

Product name Thyroxine (Total T4) ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Sera	20			<=8.16%

Inter-assay

Sample	n	Mean	SD	CV%
Sera	10			<=8.42%

Sample type Serum, Hep Plasma, Cit plasma

Assay type Competitive

Sensitivity > 0.004 µg/ml

Range 0.044 µg/ml - 0.108 µg/ml

Recovery 97.8 %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's Thyroxine (T4) *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of T4 in serum and plasma.

A 96-well plate has been precoated with anti-Thyroxine (T4) antibodies (solid-phase). Samples and the Thyroxine (T4)-HRP conjugate are added to the wells, where Thyroxine (T4) in the sample competes with the added Thyroxine (T4)-HRP conjugate for antibody binding. 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 inversely proportional to the amount of Thyroxine (T4) in the sample and the intensity is measured at 450 nm.

Platform Microplate

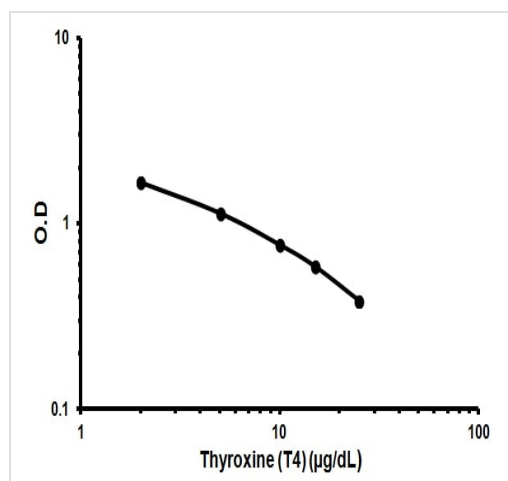
Properties

Storage instructions

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

Components	1 x 96 tests
50X Washing Solution	1 x 20ml
Anti-Thyroxine IgG Coated Microplate (12 x 8 wells)	1 unit
Conjugate Buffer	1 x 12.5ml
Cover Foil	1 unit
Stop Solution	1 x 15ml
Strip holder	1 unit
Thyroxine HRP Conjugate	1 x 1.4ml
Thyroxine Standard 0 - 0 µg/mL	1 x 1ml
Thyroxine Standard 1	1 x 1ml
Thyroxine Standard 2 - 0.050 µg/mL	1 x 1ml
Thyroxine Standard 3 - 0.100 µg/mL	1 x 1ml
Thyroxine Standard 4 - 0.150 µg/mL	1 x 1ml
Thyroxine Standard 5 - 0.250 µg/mL	1 x 1ml
TMB Substrate Solution	1 x 15ml

Images



Typical Standard Curve

Representative standard curve using ab178664.

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