

Product datasheet

Human Thyroxine ELISA Kit (fT4) ab108686

[2 References](#) [1 Image](#)

Overview

Product name	Human Thyroxine ELISA Kit (fT4)
Detection method	Colorimetric
Sample type	Serum, Hep Plasma, Cit plasma
Assay type	Competitive
Sensitivity	0.5 pg/ml
Recovery	97.8 %
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	Abcam's Thyroxine (fT4) in vitro competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of free Thyroxine in Human serum and plasma.

A 96-well plate has been precoated with anti-Thyroxine antibodies. Samples and the Thyroxine-HRP conjugate are added to the wells, where Thyroxine in the sample competes with the added Thyroxine-HRP 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 in the sample and the intensity is measured at 450 nm.

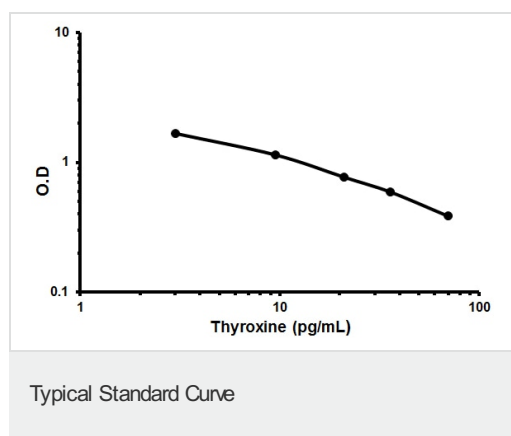
Platform	Microplate
-----------------	------------

Properties

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

Components	1 x 96 tests
1X Thyroxine-HRP Conjugate	1 x 12ml
50X Wash Solution	1 x 20ml
Anti-Thyroxine IgG Coated Microplate (12 x 8 wells)	1 unit
Cover foils	1 unit
Stop Solution	1 x 15ml
Strip holder	1 unit
Thyroxine Standard 0 – 0.0 pg/mL	1 x 1ml
Thyroxine Standard 1 – 3.0 pg/mL	1 x 1ml
Thyroxine Standard 2 – 9.5 pg/mL	1 x 1ml
Thyroxine Standard 3 – 21.0 pg/mL	1 x 1ml
Thyroxine Standard 4 – 36.0 pg/mL	1 x 1ml
Thyroxine Standard 5 – 70.0 pg/mL	1 x 1ml
TMB Substrate Solution	1 x 15ml

Images



Representative Standard Curve using ab108686.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Promise 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