

Human Thyroglobulin ELISA Kit ab155441

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

Overview

Product name Human Thyroglobulin ELISA Kit

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
overall				< 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 0.22 ng/ml

Range 0.205 ng/ml - 50 ng/ml

Recovery Sample specific recovery

Sample type	Average %	Range
Serum	115.5	105% - 125%
Plasma	98.71	88% - 106%
Cell culture media	96.01	88% - 103%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Thyroglobulin Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of Human Thyroglobulin in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human Thyroglobulin coated on a 96-well plate. Standards and samples are pipetted into the wells and Thyroglobulin present in a sample is

bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Thyroglobulin antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Thyroglobulin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples.

Platform Microplate

Properties

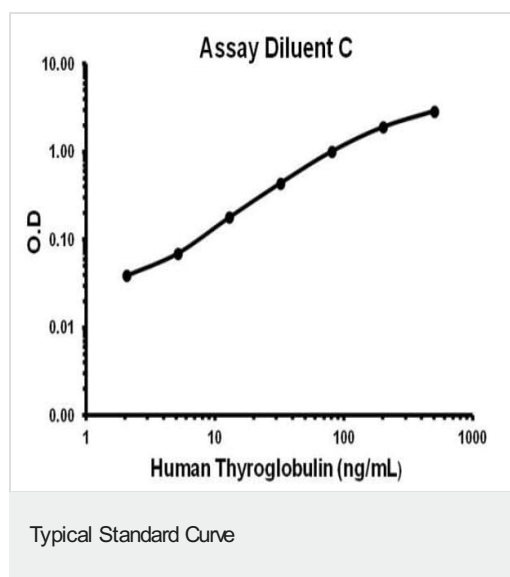
Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent C	1 x 30ml
Biotinylated anti-Human Thyroglobulin	2 vials
Recombinant Human Thyroglobulin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
Thyroglobulin Microplate (12 x 8 wells)	1 unit
TMB One-Step Substrate Reagent	1 x 12ml

Relevance Precursor of the iodinated thyroid hormones thyroxine (T4) and triiodothyronine (T3).

Cellular localization Secreted.

Images



Representative standard curve using ab155441 - Assay Diluent C.

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