

## Product datasheet

# Human Anti-Thyroglobulin ELISA Kit ab178631

1 References 1 Image

### Overview

**Product name** Human Anti-Thyroglobulin ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Pos. Serum	20			2.1%
Pos. Serum	20			1.2%

Inter-assay

Sample	n	Mean	SD	CV%
Pos. Serum	12	59.7U/ml		5.7%
Pos. Serum	12	28.9U/ml		8.5%
Neg. Serum	12	7.3U/ml		3.9%

**Sample type** Serum, Hep Plasma, Cit plasma

**Assay type** Indirect

**Sensitivity** < 0.5 U/ml

**Range** 2 U/ml - 128 U/ml

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview**

Abcam's anti-Thyroglobulin *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative determination of autoantibodies to thyroglobulin in Human serum or plasma.

A 96-well plate has been precoated with Human thyroglobulin to bind cognate antibodies. Controls or test samples are added to the wells and incubated. Following washing, a horseradish peroxidase (HRP) labelled anti-Human IgG conjugate is added to the wells, which binds to the immobilized thyroglobulin-specific antibodies. TMB is then catalyzed by the HRP to produce a blue color product that changes to yellow after adding an acidic stop solution. The intensity of

yellow coloration is directly proportional to the amount of thyroglobulin sample captured in plate.

**Platform** Microplate

## Properties

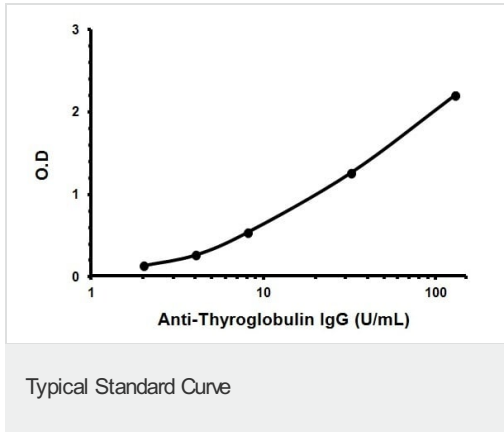
---

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

Components	Identifier	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
Cover Foil		1 unit
IgG Sample Diluent		1 x 100ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
Thyroglobulin Anti-IgG HRP Conjugate		1 x 20ml
Thyroglobulin Coated Microplate (12 x 8 wells)		1 unit
Thyroglobulin Negative Control		1 x 2ml
Thyroglobulin Positive Control		1 x 2ml
Thyroglobulin Standard A - 0 U/mL		1 x 2ml
Thyroglobulin Standard B - 2 U/mL		1 x 2ml
Thyroglobulin Standard C - 4 U/mL		1 x 2ml
Thyroglobulin Standard D - 8 U/mL		1 x 2ml
Thyroglobulin Standard E - 32 U/mL		1 x 2ml
Thyroglobulin Standard F - 128 U/mL		1 x 2ml
TMB Substrate Solution	Yellow cap	1 x 15ml

## Images

---



Representative standard curve using ab178631

**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