

Product datasheet

Human Thyroxine ELISA Kit (T4) ab108661

★★★★☆ 1 Abreviews 1 Image

Overview

Product name	Human Thyroxine ELISA Kit (T4)
Detection method	Colorimetric
Sample type	Serum
Assay type	Competitive
Sensitivity	= 0.4 µg/dl
Range	0.4 µg/dl - 25 µg/dl
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	<p>ab108661, Thyroxine (T4) Human ELISA Kit is intended for the quantitative determination of Total Thyroxine (T4) concentration in Human serum.</p> <p>L-Thyroxine (T4) is a hormone that is synthesized and stored in the thyroid gland. Proteolytic cleavage of follicular thyroglobulin releases T4 into the bloodstream. Greater than 99% of Thyroxine (T4) is reversibly bound to three plasma proteins in blood - thyroxine binding globulin (TBG) binds 70%, thyroxine binding pre-albumin (TBPA) binds 20%, and albumin binds 10%. Approximately 0.03% of T4 is in the free, unbound state in blood at any one time.</p>
Platform	Microplate

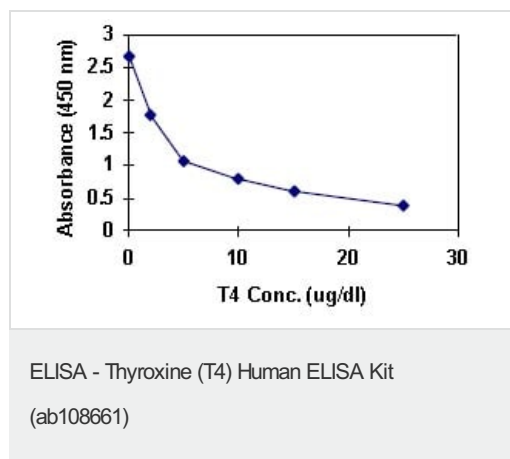
Properties

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

Components	1 x 96 tests
Enzyme Conjugate Concentrate (11x)	1 x 1.3ml
Enzyme Conjugate Diluent	1 x 13ml
Sheep anti-Thyroxine (T4) coated microtiter wells	1 x 96 units
Stop Solution	1 x 11ml

Components	1 x 96 tests
Thyroxine (T4) Reference Standard (0 µg/dl)	1 x 1ml
Thyroxine (T4) Reference Standard (10 µg/dl)	1 x 1ml
Thyroxine (T4) Reference Standard (15 µg/dl)	1 x 1ml
Thyroxine (T4) Reference Standard (2 µg/dl)	1 x 1ml
Thyroxine (T4) Reference Standard (25 µg/dl)	1 x 1ml
Thyroxine (T4) Reference Standard (5 µg/dl)	1 x 1ml
TMB Reagent	1 x 11ml

Images



Thyroxine (T4) Standard Curve

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