

Triiodothyronine (T3) ELISA Kit ab108685

★★★★★ [1 Abreviews](#) [4 References](#) [画像数 1](#)

製品の概要

製品名	Triiodothyronine (T3) ELISA Kit										
検出方法	Colorimetric										
再現性	Intra-Assay (同時再現性)										
	<table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td>40</td> <td></td> <td></td> <td>< 10.7%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall	40			< 10.7%
サンプル	N	平均値	SD	CV%							
Overall	40			< 10.7%							
	Inter-Assay (日差再現性)										
	<table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td>3</td> <td></td> <td></td> <td>< 9.1%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall	3			< 9.1%
サンプル	N	平均値	SD	CV%							
Overall	3			< 9.1%							
サンプルの種類	Serum, Hep Plasma, Cit plasma										
アッセイタイプ	Competitive										
検出感度	0.05 ng/ml										
添加回収試験	97.5 %										
ステップ	Multiple steps standard assay										
種交差性	交差種: Human										
製品の概要	Abcam's Triiodothyronine (T3) <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of total Triiodothyronine in serum and plasma.										

A 96-well plate has been precoated with anti-T3. Samples and the T3-HRP conjugate are added to the wells, where Triiodothyronine in the sample competes with the added T3-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 Triiodothyronine in the sample and the intensity is measured at 450 nm.

製品の特性

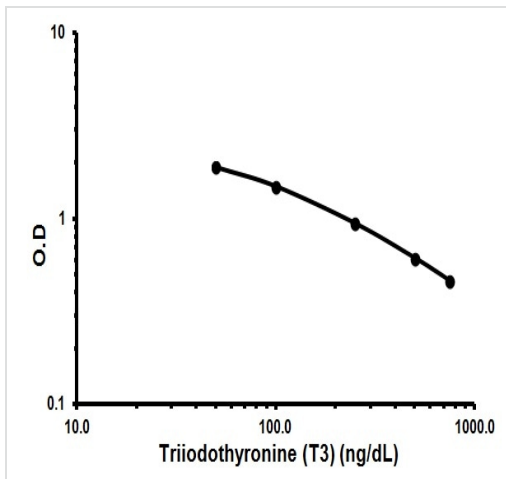
保存方法

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

内容	1 x 96 tests
11X T3-HRP Conjugate	1 x 1.4ml
50X Wash Solution	1 x 20ml
Anti-T3 IgG Coated Microplate (12 x 8 wells)	1 unit
Conjugate Buffer	1 x 12.5ml
Cover foils	1 unit
Stop Solution	1 x 15ml
Strip holder	1 unit
T3 Standard 0 - 0.0 ng/mL	1 x 1ml
T3 Standard 1 - 0.5 ng/mL	1 x 1ml
T3 Standard 2 - 1.0 ng/mL	1 x 1ml
T3 Standard 3 - 2.5 ng/mL	1 x 1ml
T3 Standard 4 - 5.0 ng/mL	1 x 1ml
T3 Standard 5 - 7.5 ng/mL	1 x 1ml
TMB Substrate Solution	1 x 15ml

画像

Representative Standard Curve using ab108685



Typical 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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors