

Human IgD ELISA Kit ab157708

画像数 1

医薬用外毒物

医薬用外劇物

製品の概要

製品名	Human IgD ELISA Kit																				
検出方法	Colorimetric																				
再現性	Intra-Assay (同時再現性) <table><tr><th>サンプル</th><th>N</th><th>平均値</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td></td><td></td><td></td><td>< 10%</td></tr></table> Inter-Assay (日差再現性) <table><tr><th>サンプル</th><th>N</th><th>平均値</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td></td><td></td><td></td><td>< 10%</td></tr></table>	サンプル	N	平均値	SD	CV%	Overall				< 10%	サンプル	N	平均値	SD	CV%	Overall				< 10%
サンプル	N	平均値	SD	CV%																	
Overall				< 10%																	
サンプル	N	平均値	SD	CV%																	
Overall				< 10%																	
サンプルの種類	Serum, Plasma, Other biological fluids																				
アッセイタイプ	Sandwich (quantitative)																				
検出感度	4.4545 ng/ml																				
検出範囲	6.25 ng/ml - 400 ng/ml																				
添加回収試験	> 85 %																				
ステップ	Multiple steps standard assay																				
種交差性	交差種: Human																				
製品の概要	Abcam's IgD Human ELISA kit is an <i>in vitro</i> enzyme-linked immunosorbent assay (ELISA) for the																				

製品の特性

保存方法

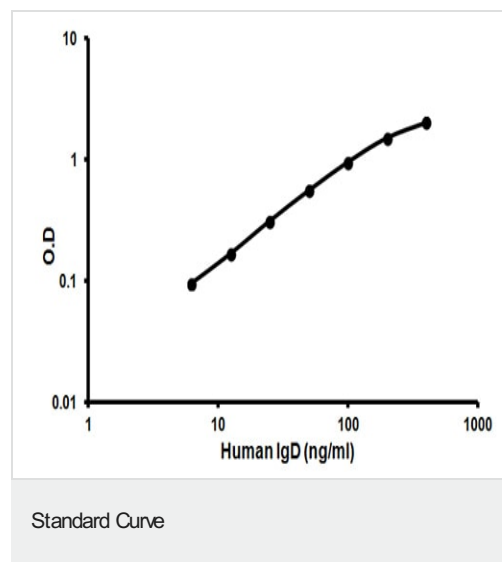
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
100X HRP-conjugated anti-Human IgD antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Chromogen Substrate Solution	1 x 12ml
Human IgD Calibrator (lyophilized)	1 vial
Human IgD ELISA Microplate	1 unit
Stop Solution	1 x 12ml

関連性

Little is known about the normal function of IgD, and few clinical signs or symptoms are associated with its absence. Individuals with low or absent levels of IgD do not appear unusually predisposed to infections. In mice, IgD may substitute for some functions of IgM when IgM is absent. IgD is expressed on peripheral B cells and is used as a pan B cell marker.

画像



Representative standard curve using ab157708 IgD Human ELISA Kit.

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