

Product datasheet

Mouse beta IG H3 Matched Antibody Pair Kit ab215413

1 Image

Overview

Product name	Mouse beta IG H3 Matched Antibody Pair Kit
Detection method	Colorimetric
Assay type	ELISA set
Sensitivity	16.5 pg/ml
Range	62.5 pg/ml - 4000 pg/ml
Species reactivity	Reacts with: Mouse
Product overview	<p>Mouse beta IG H3 Matched Antibody Pair Kits include a capture and a biotinylated detector antibody pair, along with a calibrated protein standard, suitable for sandwich ELISA. The Matched Antibody Pair Kit can be used to quantify native and recombinant mouse beta IG H3.</p>

Optimization of the kit reagents to sample type, immunoassay format or instrumentation may be required. Guidelines for use of this kit in a standard 96-well microplate sandwich ELISA using HRP/TMB system of colorimetric detection is described in this assay procedure for the purposes of quantification.

Protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our [website](#). An accessory pack can be purchased which includes buffer reagents required to perform 10 x 96-well plate sandwich ELISAs ([ab210905](#)).

For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA kit [ab206987](#). Please note that while the antibody pair is the same provided in the corresponding SimpleStep ELISA Kit, due to differences in their formulation, this antibody pair cannot be used with the consumables provided with our SimpleStep ELISA Kits.

We've listened to you: due to popular demand, we will now provide our Matched Antibody Pair kits in 5x96 tests and 10x96 tests. The 2x96 tests size will be discontinued on 30th June 2020 unless inventory is depleted beforehand.

To receive an electronic copy of the Certificate of Analysis, please send an [email](#) with "CoA for matched antibody pair kit" in the subject line and the desired product number and lot number in the body of the email.

Tested applications	Suitable for: ELISA, IA
Platform	Reagents

Properties

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

Components	10 x 96 tests	5 x 96 tests
Mouse beta IG H3 Capture Antibody	1 x 100µg	1 x 50µg
Mouse beta IG H3 Detector Antibody	1 x 25µg	1 x 12.5µg
Mouse beta IG H3 Lyophilized Protein	1 x 1µg	1 vial

Function Binds to type I, II, and IV collagens. This adhesion protein may play an important role in cell-collagen interactions. In cartilage, may be involved in endochondral bone formation.

Tissue specificity Highly expressed in the corneal epithelium.

Involvement in disease Defects in TGFBI are the cause of epithelial basement membrane corneal dystrophy (EBMD) [MIM:121820]; also known as Cogan corneal dystrophy or map-dot-fingerprint type corneal dystrophy. EBMD is a bilateral anterior corneal dystrophy characterized by grayish epithelial fingerprint lines, geographic map-like lines, and dots (or microcysts) on slit-lamp examination. Pathologic studies show abnormal, redundant basement membrane and intraepithelial lacunae filled with cellular debris. Although this disorder usually is not considered to be inherited, families with autosomal dominant inheritance have been identified.

Defects in TGFBI are the cause of corneal dystrophy Groenouw type 1 (CDGG1) [MIM:121900]; also known as corneal dystrophy granular type. Inheritance is autosomal dominant. Corneal dystrophies show progressive opacification of the cornea leading to severe visual handicap.

Defects in TGFBI are the cause of corneal dystrophy lattice type 1 (CDL1) [MIM:122200]. Inheritance is autosomal dominant.

Defects in TGFBI are a cause of corneal dystrophy Thiel-Behnke type (CDTB) [MIM:602082]; also known as corneal dystrophy of Bowman layer type 2 (CDB2).

Defects in TGFBI are the cause of Reis-Buecklers corneal dystrophy (CDRB) [MIM:608470]; also known as corneal dystrophy of Bowman layer type 1 (CDB1).

Defects in TGFBI are the cause of lattice corneal dystrophy type 3A (CDL3A) [MIM:608471]. CDL3A clinically resembles to lattice corneal dystrophy type 3, but differs in that its age of onset is 70 to 90 years. It has an autosomal dominant inheritance pattern.

Defects in TGFBI are the cause of Avellino corneal dystrophy (ACD) [MIM:607541]. ACD could be considered a variant of granular dystrophy with a significant amyloidogenic tendency. Inheritance is autosomal dominant.

Sequence similarities Contains 1 EMI domain.
Contains 4 FAS1 domains.

Post-translational modifications Gamma-carboxyglutamate residues are formed by vitamin K dependent carboxylation. These residues are essential for the binding of calcium.

Cellular localization Secreted > extracellular space > extracellular matrix. May be associated both with microfibrils and with the cell surface.

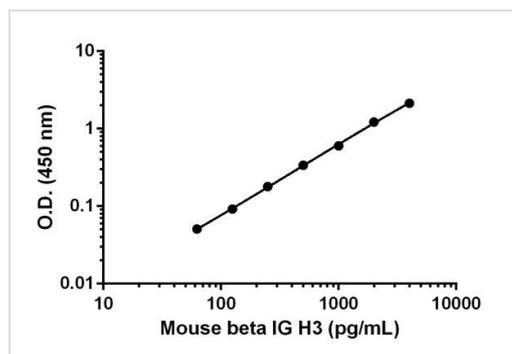
Applications

Our [Abpromise guarantee](#) covers the use of **ab215413** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
ELISA		Use at an assay dependent concentration.
IA		Use at an assay dependent concentration.

Images



Standard calibration curve. Background subtracted values are graphed.

Mouse beta IG H3 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