

Anti-Human cMyBP-C Antibodies (9B11C5), mAb, Mouse

PRODUCT INFORMATION

Description

cMyBP-C is not essential for sarcomere formation during embryogenesis, but is crucial for sarcomere organization and maintenance of normal cardiac function. Absence of cMyBP-C results in severe cardiac hypertrophy, increased heart-weight-to-body-weight-ratios, enlargement of ventricles, increased myofilament Ca²⁺ sensitivity and depressed diastolic and systolic function.

GenScript Anti-Human cMyBP-C Antibody (9B11C5), mAb, Mouse is produced from the hybridoma resulting from fusion of SP2/0-Ag14 myeloma and B-lymphocytes obtained from a mouse immunized with recombinant human cMyBP-C expressed in mammalian cells.

Specificity

The antibody specifically recognizes human cMyBP-C. The antibody is recommended as a capture antibody in vitro diagnostic assay development with detection antibody (GenScript, A01851-40, Anti-Human cMyBP-C Antibody (2G2F3), mAb, Mouse)

Concentration

0.5 mg/ml, lyophilized with PBS, pH 7.4, containing 0.02% sodium azide. (It could be customized for bulk order)

Note

GenScript can customize this product per customer's request including product size, buffer components, etc.

Reconstitution

Reconstitute the lyophilized antibody with deionized water (or equivalent) to a final concentration of 0.5 mg/ml.

Storage

GenScript Anti-Human cMyBP-C Antibody (9B11C5), mAb, Mouse should be stored at -20°C. This product is stable for 1 year upon receipt, when handled and stored as instructed.

Applications

Working concentrations for specific applications should be determined by the investigator. The appropriate concentrations may be affected by secondary antibody affinity, antigen concentration, the sensitivity of the method of detection, temperature, the length of the incubations, and other factors. The suitability of this antibody for applications other than those listed below has not been determined. The following concentration ranges are recommended starting points for this product.

Western Blotting: 1:2,000

ELISA: 0.5-1 µg/ml

Other applications: user-optimized

Cat. No.: A01850-40

Host: Mouse

Size: 40 µg

Immunogen: Recombinant human cMyBP-C

Ig Subclass: IgG2b,κ

Clone: 9B11C5

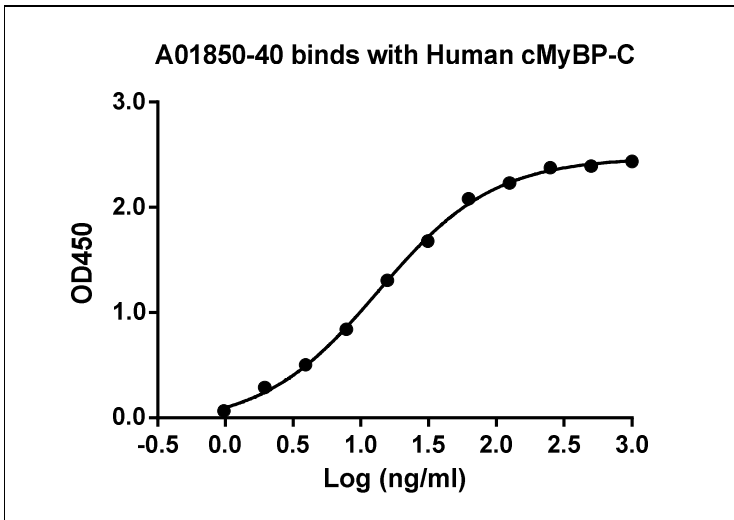
Purification: Protein A affinity column

Conjugation: Unconjugated

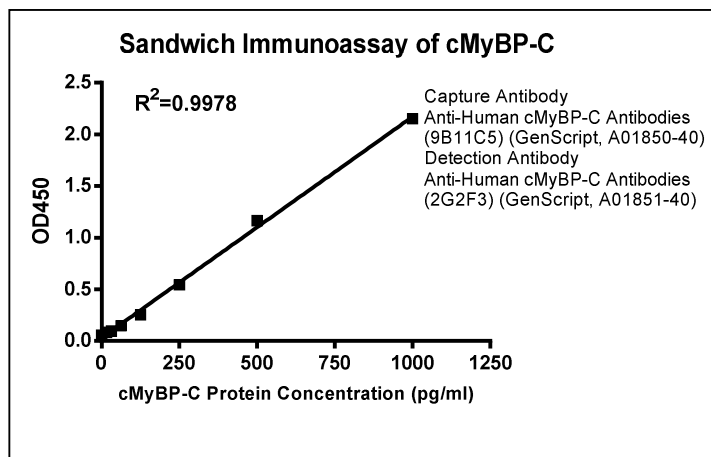
Version: 01/03/2017

** For non-clinical research use only. **

860 Centennial Ave., Piscataway, NJ 08854, USA

Example


Anti-Human cMyBP-C Antibodies (9B11C5) (GenScript, A01850-40) binds with Human cMyBP-C.
 Coating antigen: Human cMyBP-C, 1 µg/ml.
 Anti-Human cMyBP-C Antibodies (GenScript, A01850-40) dilution start from 1,000 ng/ml.
 EC_{50} = 13.4 ng/ml.



Antibody pairs analysis of cMyBP-C monoclonal antibodies by Sandwich ELISA:

General conditions:

1. Microplate was incubated with a capture antibody (GenScript, A01850-40) against cMyBP-C and blocked with PBS containing 1% BSA, followed by 3 washing cycles.
2. Incubation with cMyBP-C protein followed by 3 washing cycles.
3. Incubation with Biotin labeled detection antibody (GenScript, A01851-40) against cMyBP-C, followed by 3 washing cycles.
4. Incubation with Streptavidin-HRP, followed by 3 washing cycles.
5. Colorimetric determination of bound peroxidase activity.

** For non-clinical research use only. **

860 Centennial Ave., Piscataway, NJ 08854, USA