

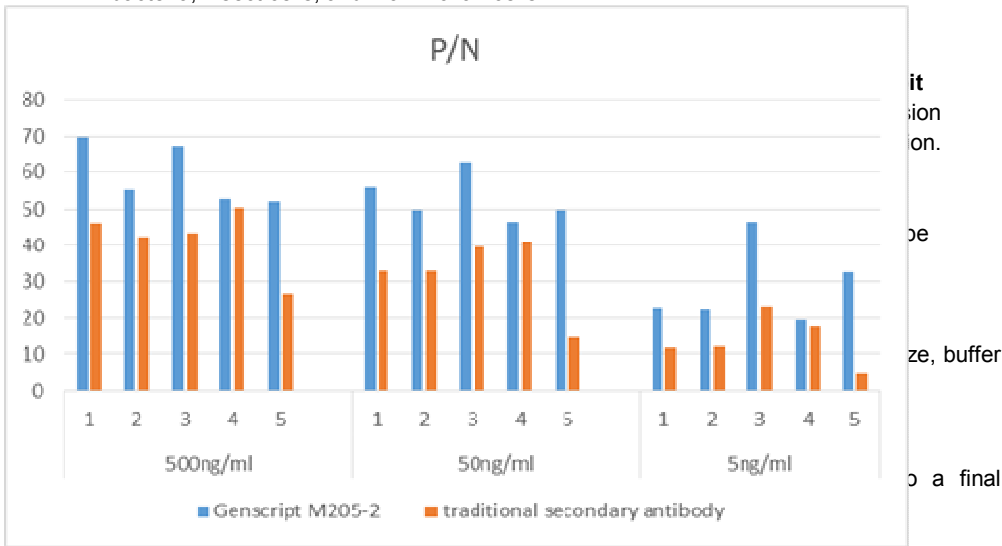
MonoRab™ Anti-His Tag (C-term) Antibody (25B6E11), mAb, Rabbit

Cat. No.: A01857-40
Host: Rabbit
Size: 40 µg
Immunogen: 6x His synthetic peptide coupled to KLH
Ig Subclass: IgG
Clone: 25B6E11
Purification: Protein A affinity column
Conjugation: Unconjugated
Version: 01/03/2017

PRODUCT INFORMATION

Description

The rabbit immune system generates antibody diversity and optimizes affinity. GenScript utilizes MonoRab™ technology to generate high affinity and specific monoclonal rabbit antibodies. **GenScript MonoRab™ Anti-His Tag (C-term) Antibody (25B6E11), mAb, Rabbit** is specific to six histidine tags placed at C-terminal and internal regions of fusion proteins. This antibody can greatly improve the effectiveness of several different kinds of immunoassays, helping researchers identify, detect, and purify polyhistidine fusion proteins in bacteria, insect cells, and mammalian cells.



Storage

GenScript MonoRab™ Anti-His Tag (C-term) Antibody (25B6E11), mAb, Rabbit 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.

ELISA: 0.05-0.2 µg/ml

Western Blot: 0.1-0.2 µg/ml

Immunoprecipitation (IP): 2 µg/ml

Immunofluorescent staining: 1 µg/ml

Flow cytometry (FACS): 1 µg/ml

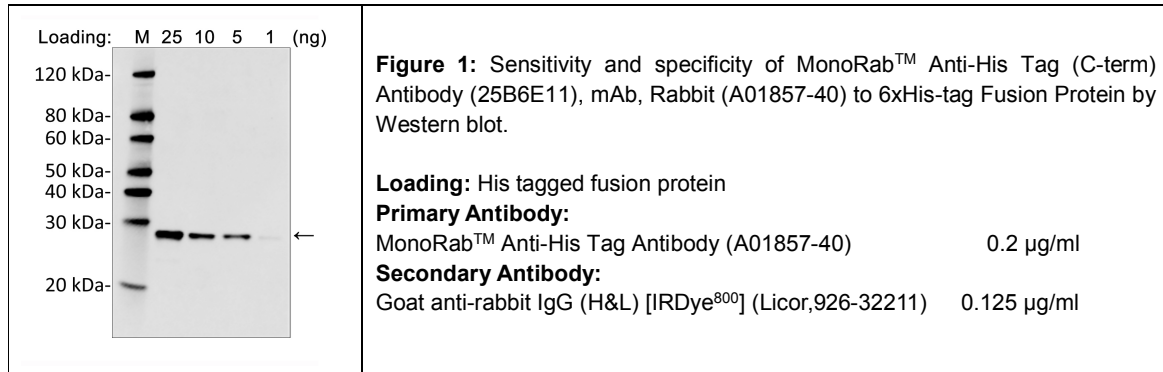
Other applications: user-optimized

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

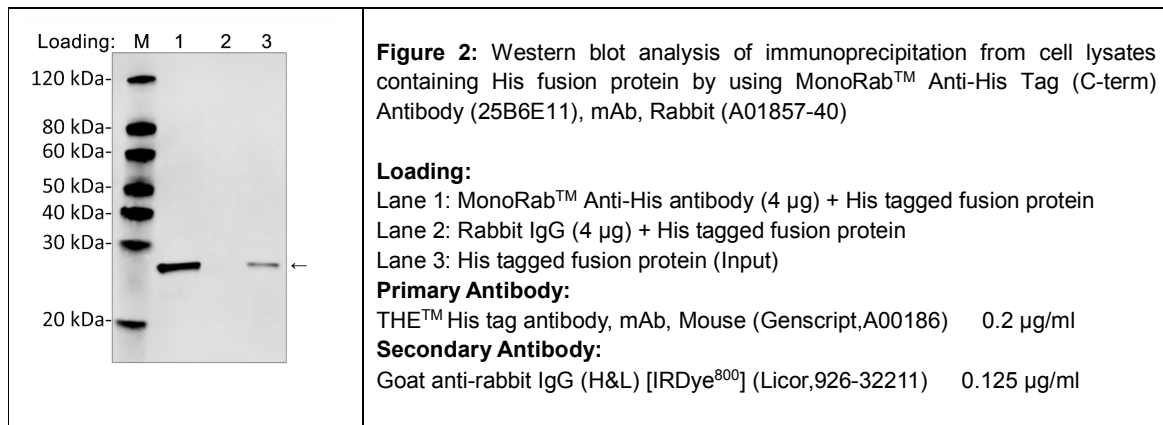
860 Centennial Ave., Piscataway, NJ 08854, USA

Example

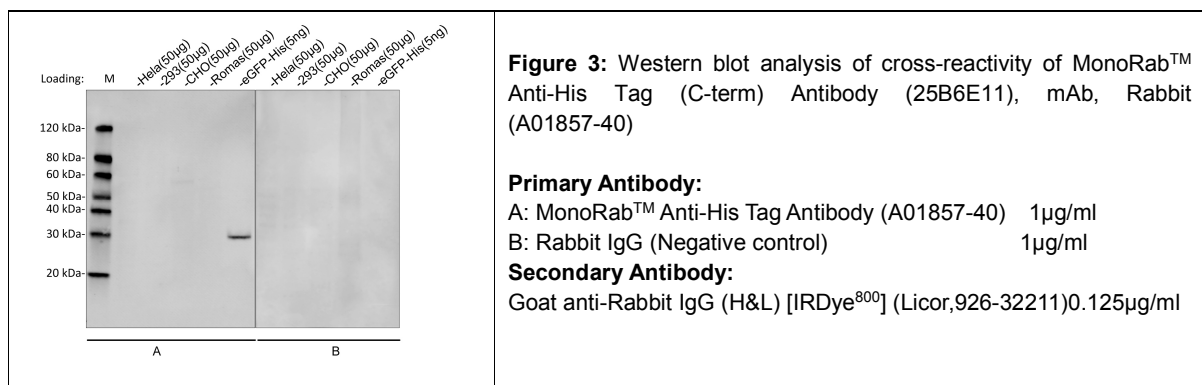
- Western Blot of 6xHis-tag Fusion Protein using MonoRab™ Anti-His Tag (C-term) Antibody (25B6E11), mAb, Rabbit (A01857-40)



- Western blot analysis of immunoprecipitation using MonoRab™ Anti-His Tag (C-term) Antibody (25B6E11), mAb, Rabbit (A01857-40)



- Western blot analysis of cross-reactivity using MonoRab™ Anti-His Tag (C-term) Antibody (25B6E11), mAb, Rabbit (A01857-40)



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

860 Centennial Ave., Piscataway, NJ 08854, USA