

# PRODUCT INFORMATION

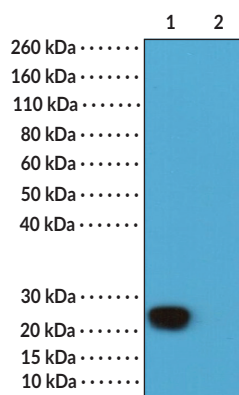


## HA-Tag Chimeric Rabbit-Human Monoclonal Antibody (Clone RMH02) Item No. 35900

### Overview and Properties

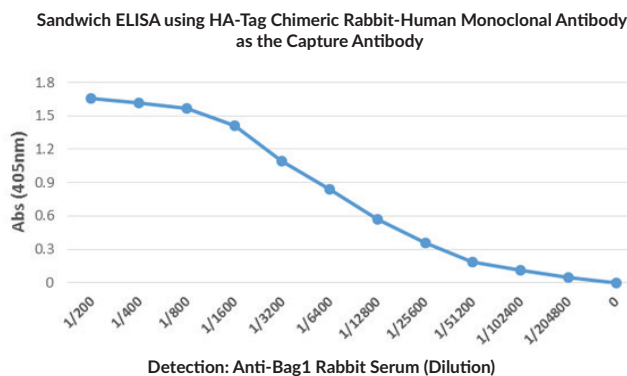
|                            |   |
|----------------------------|---|
| <b>Contents:</b>           | This vial contains 100 µg of protein A-affinity purified monoclonal antibody.   |
| <b>Synonym:</b>            | Hemagglutinin   |
| <b>Immunogen:</b>          | A peptide corresponding to HA-tag   |
| <b>Cross Reactivity:</b>   | (+) HA-tagged proteins<br>(-) Endogenous mammalian or bacterial proteins  |
| <b>Species Reactivity:</b> | (+) Species independent   |
| <b>Form:</b>               | Liquid  |
| <b>Storage:</b>            | -20°C (as supplied)   |
| <b>Stability:</b>          | ≥1 year   |
| <b>Storage Buffer:</b>     | PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide   |
| <b>Clone:</b>              | RMH02   |
| <b>Host:</b>               | This antibody is composed of a fusion of the antigen binding region of rabbit anti-HA-Tag monoclonal antibody (clone RMH02) with the constant domain of human IgG1  |
| <b>Isotype:</b>            | Rabbit Fv fused to human IgG1 CH and Cκ   |
| <b>Applications:</b>       | ELISA and Western blot (WB); the recommended starting concentrations for ELISA are 0.5-2 µg/ml (capturer) and 0.01-0.5 µg/ml (detector) and is 0.1-1 µg/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

### Images



Lane 1: 293T cells transfected  
Lane 2: 293T cells untransfected

WB of 293T cells transfected with DNA construct encoding HA-Tag KRas fusion protein or left untransfected using HA-Tag Chimeric Rabbit-Human Monoclonal Antibody (Clone RMH02) at a concentration of 0.1 µg/ml, followed by a HRP-conjugated anti-Human IgG secondary antibody.



Detection of HA-tag Bag1 protein in Rabbit Serum. Sandwich ELISA using HA-Tag Chimeric Rabbit-Human Monoclonal Antibody (Clone RMH02) (Item No. 35900) (1 µg/ml) as the capture antibody and Anti-Bag1 rabbit serum as the detection antibody, followed by an alkaline phosphatase-conjugated anti-Rabbit IgG as the secondary antibody. The plate was coated with 50 µl/well of HA-Tag Chimeric Rabbit-Human Monoclonal Antibody (Clone RMH02) (Item No. 35900).

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/30/2024

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

HA-Tag Chimeric Rabbit-Human Monoclonal Antibody is a probe for the immunochemical detection of HA tags on recombinant proteins. Recombinant proteins are commonly labeled with affinity tags, such as HA, which corresponds to amino acids 98 to 106 from human influenza hemagglutinin (HA), to facilitate both their detection and purification.<sup>1</sup> HA tags are commonly utilized because of their small size, low potential to interfere in protein folding or activity, and strong immunoreactivity. Cayman's HA-Tag Chimeric Rabbit-Human Monoclonal Antibody (Clone RMH02) is composed of human IgG1 CH and C $\kappa$  constant domains fused to the antigen-binding domain of a rabbit anti-HA-tag monoclonal antibody. It can be used for ELISA and Western blot (WB) applications. The antibody recognizes proteins containing HA tags fused to either the N- or C-terminus.

## Reference

---

1. Zhao, X., Li, G., and Liang, S. Several affinity tags commonly used in chromatographic purification. *J. Anal. Methods Chem.* 581093 (2013).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM