

# PRODUCT INFORMATION



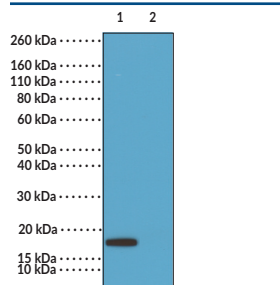
## His-tag Monoclonal Antibody (Clone RMH01)

Item No. 32187

### Overview and Properties

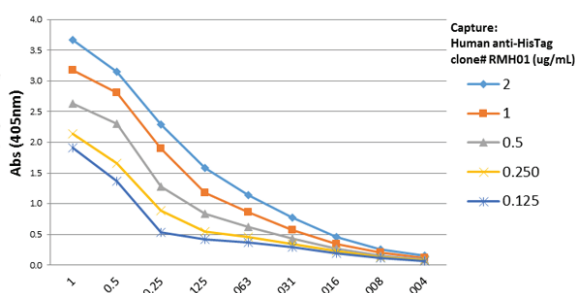
<b>Contents:</b>	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
<b>Synonyms:</b>	Deca-His Tag, Decahistidine Tag, Hexa-His Tag, Hexahistidine Tag, Poly-His Tag
<b>Immunogen:</b>	Mixture of a peptide with a 6x His tag at the N-terminus and a peptide with a 6x His tag at the C-terminus
<b>Cross Reactivity:</b>	(+) His-tagged proteins; (-) 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>Concentration:</b>	1 mg/ml
<b>Clone:</b>	RMH01
<b>Host:</b>	Chimeric Monoclonal Antibody
<b>Isotype:</b>	IgG
<b>Applications:</b>	ELISA and Western blot (WB); the recommended starting concentration is 0.1-2 µg/ml for capture and 0.01-0.5 µg/ml for detection for ELISA and 0.1-1 µg/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



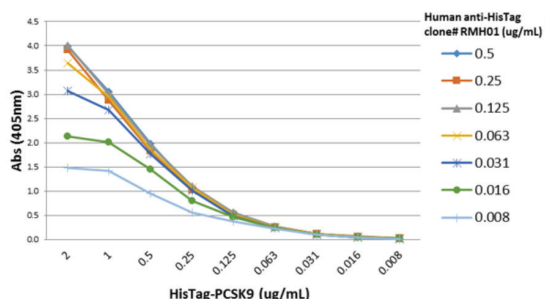
Lane 1: 293T cells transfected  
Lane 2: 293T cells untransfected

WB of 293T cells transfected with DNA construct encoding His-Tag histone H3 (G34W) protein or left untransfected using His-tag Monoclonal Antibody (Clone RMH01) at a concentration of 0.2 µg/ml, followed by a HRP-conjugated



Whole Lysate of 294T Expressing HisTag-H3K36M (mg/mL, 50uL/well)

Detection of His-Tag histone H3 (K36M) protein in whole lysate of 294T cells transfected with a DNA construct encoding His-tag histone H3 (K36M) protein. Sandwich ELISA using His-tag Monoclonal Antibody (Clone RMH01) as the capture antibody (50 µg/well) and Histone H3K36M Monoclonal Antibody (Item No. 31547) as the detection antibody, followed by an alkaline phosphatase-conjugated rabbit IgG as the secondary antibody.



A Titer ELISA using His-tag Monoclonal Antibody. The plate was coated with different amounts of purified HisTag-PCSK9 recombinant protein. A serial dilution of His-tag Monoclonal Antibody (Clone RMH01) was used as the primary antibody and an alkaline phosphatase-conjugated anti-human IgG as the secondary antibody.

**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, 08/16/2024

**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

---

His-tag Monoclonal Antibody is a probe for the immunochemical detection of histidine tags on recombinant proteins. Recombinant proteins are commonly labeled with affinity tags, such as 6-10 histidine residues (6x-10x His), to facilitate both their detection and purification.<sup>1</sup> Poly-His tags are commonly utilized because of their small size, low potential to interfere in protein folding or activity, weak immunogenicity, and high affinity for transition metal ion matrices, such as Ni<sup>2+</sup>, used in immobilized metal-affinity chromatography (IMAC) for protein purification.<sup>1,2</sup> Cayman's His-tag Monoclonal Antibody (Clone RMH01) is comprised of rabbit IgG variable domains and the human IgG1 constant domain. This antibody can be used for ELISA and Western blot (WB) applications. It recognizes proteins containing 6x or 10x His tags fused to either the N- or C-terminus and is recognized by anti-human IgG secondary antibodies.

## References

---

1. Terpe, K. Overview of tag protein fusions: From molecular and biochemical fundamentals to commercial systems. *Appl. Microbiol. Biotechnol.* **60(5)**, 523-533 (2003).
2. Spriestersbach, A., Kubicek, J., Schäfer, F., *et al.* Purification of His-tagged proteins. *Methods in Enzymology*. Lorsch, J.R., editor, 1<sup>st</sup> edition, *Academic Press* (2015).

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