

PRODUCT INFORMATION

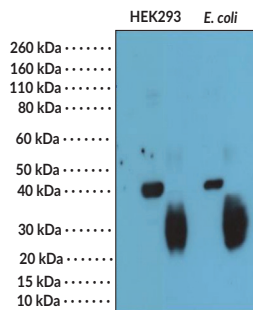
His-Tag Rabbit Monoclonal Antibody (Clone RM146X)

Item No. 42809

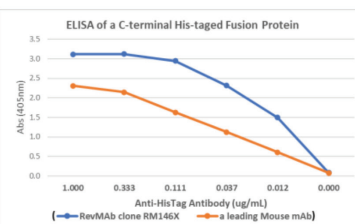
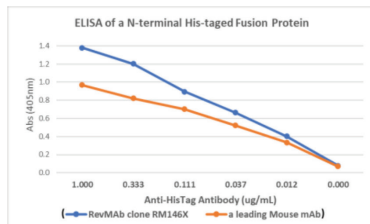
Overview and Properties

Contents:	This vial contains 100 µg of protein A-affinity purified monoclonal antibody
Synonyms:	6x-His Tag, Deca-His Tag, Decahistidine Tag, Hexa-His Tag, Hexahistidine Tag, Poly-His Tag, Polyhistidine Tag
Immunogen:	6xHis-tag coupled to KLH
Cross Reactivity:	(+) N- and C-terminal His-tagged recombinant proteins; (-) Very low non-specific reactivity with endogenous proteins in mammalian or bacterial cells compared with other His-Tag antibodies
Species Reactivity:	Species Independent
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
Concentration:	1.0 mg/ml
Clone:	RM146X
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA, Immunocytochemistry (ICC), Immunoprecipitation (IP), and Western blot (WB); the recommended concentration is 0.01-0.2 µg/ml for ELISA, 0.5-2 µg/ml for ICC and IP, and 0.01-0.1 µg/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

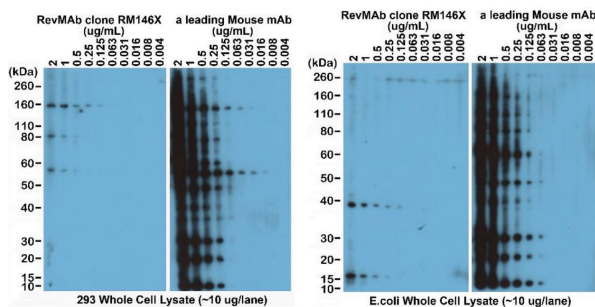
Images



WB of 293 cell or E. coli whole cell lysate with or without N-terminal His-tag fusion protein X or C-terminal His-tag fusion protein Y using His-Tag Rabbit Monoclonal Antibody (Clone RM146X).



ELISA comparison of a N-terminal His-tag fusion protein (50 µg/well) and a C-terminal His-tag fusion protein (50 µg/well), using His-Tag Rabbit Monoclonal Antibody (Clone RM146X), versus a leading anti-His-tag mouse monoclonal antibody.



Western blot of His-Tag Rabbit Monoclonal Antibody (Clone RM146X) with endogenous proteins in mammalian 293 cells or E. coli bacteria cells, compared with a leading anti-His-tag mouse monoclonal 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, 03/28/2025

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 Rabbit 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.¹ 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 Ni²⁺, used in immobilized metal-affinity chromatography (IMAC) for protein purification.^{1,2} Cayman's His-Tag Rabbit Monoclonal Antibody (Clone RM146X) can be used for ELISA, immunocytochemistry (ICC), immunoprecipitation (IP), and Western blot (WB) applications. The antibody recognizes proteins containing His-tags fused to either the N- or C-terminus.

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. Priestersbach, A., Kubicek, J., Schäfer, F., *et al.* Purification of His-tagged proteins. *Methods in Enzymology* **559**, 1-15 (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