

PRODUCT INFORMATION



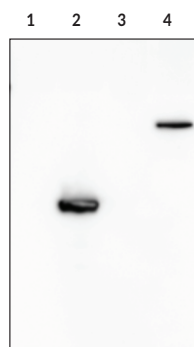
His-Tag Recombinant Monoclonal Antibody - HRP Conjugated

Item No. 35511

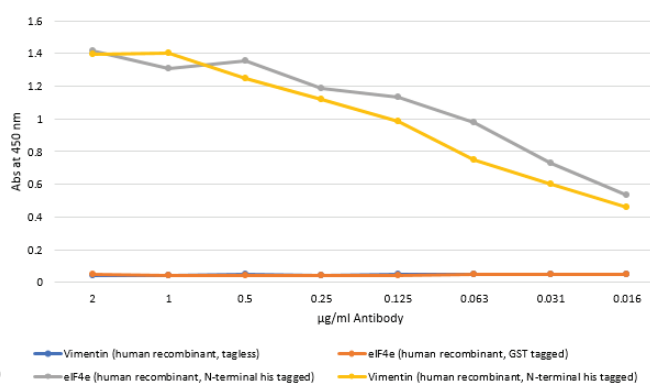
Overview and Properties

Contents:	This vial contains 100 µg of HRP-conjugated protein G-purified monoclonal antibody.
Synonyms:	Hexa-His Tag, Hexahistidine Tag, 6x-His Tag
Preparation:	This antibody, which contains a mouse IgG2a isoform, was produced as a recombinant protein expressed in HEK293-derived cells using an engineered construct based on a monoclonal anti-His antibody
Cross Reactivity:	(+) His-tagged proteins; (-) GST-tagged proteins
Species Reactivity:	Species Independent
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.05% Proclin 300
Host:	HEK293 Cells
Isotype:	IgG2a
Applications:	ELISA and Western blot (WB); the recommended starting dilution is 1:5,000 for ELISA and 1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: eIF4E (human recombinant; GST-tagged) (Item No. 15564)
Lane 2: eIF4E (human recombinant; His-tagged) (Item No. 25150)
Lane 3: Vimentin (human, recombinant, tagless)
Lane 4: Vimentin (human, recombinant) (Item No. 11234)



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 Recombinant Monoclonal Antibody - HRP Conjugated is a probe for the immunochemical detection of histidine tags on recombinant proteins. Recombinant proteins are commonly labeled with affinity tags, such as six histidine residues (6x-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 Recombinant Monoclonal Antibody - HRP Conjugated was produced as a recombinant protein using an engineered construct based on a monoclonal anti-His antibody. This antibody can be used for ELISA and Western blot (WB) applications. It recognizes proteins containing internal, N-, or C-terminal 6x-His tags and is recognized by anti-mouse 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. Priestersbach, A., Kubicek, J., Schäfer, F., *et al.* Purification of His-tagged proteins. *Methods in Enzymology*. Lorsch, J.R., editor, 1st 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