

PRODUCT INFORMATION



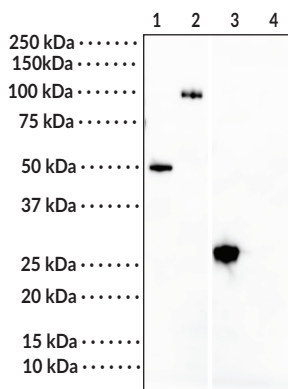
His-tag Recombinant Monoclonal Antibody

Item No. 35730

Overview and Properties

Contents:	This vial contains 100 µg of 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 0.02% sodium azide
Host:	HEK293 Cells
Isotype:	IgG2a
Applications:	ELISA and Western blot (WB); the recommended starting dilution for ELISA is 1:5,000 and 1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: CdnP (Internal His-tag) (Item No. 22809) (100 ng)
Lane 2: ENPP1 (C-terminal His-tag) (Item No. 33397) (100 ng)
Lane 3: eIF4e (N-terminal His-tag) (Item No. 25150) (50 ng)
Lane 4: eIF4e (GST-tag) (Item No. 15564) (50 ng)

Representative gel image shown; actual purity may vary between each batch.

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 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 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 6x-His tags fused to either the N- or C-terminus 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* **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