

PRODUCT INFORMATION



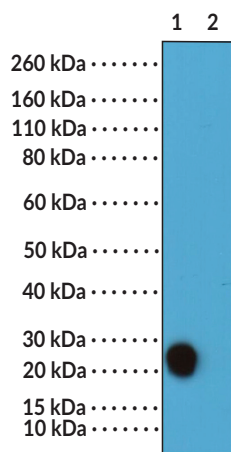
HA-tag Rabbit Monoclonal Antibody - Biotinylated (Clone RM305)

Item No. 32380

Overview and Properties

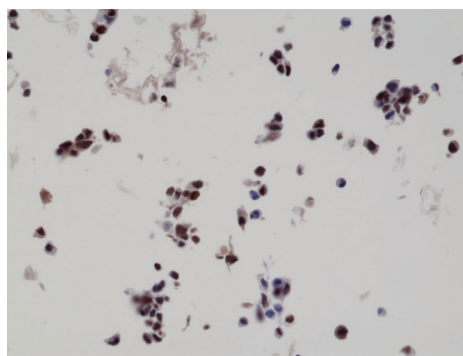
Contents:	This vial contains 50 µg of protein A-affinity purified monoclonal antibody.
Synonym:	Hemagglutinin
Immunogen:	A peptide corresponding to HA-tag
Cross Reactivity:	(+) HA-tagged proteins; (-) Endogenous mammalian or bacterial proteins
Species Reactivity:	Species Independent
Form:	Liquid
Storage:	-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:	RM305
Host:	Rabbit
Isotype:	IgG
Applications:	Flow cytometry (FC), Immunocytochemistry (ICC), Immunohistochemistry (IHC), Immunoprecipitation (IP), Western blot (WB); the recommended starting concentration is 0.1-1 µg/ml for FC, ICC, and WB, 0.01-0.5 µg/ml for IHC, and 0.5-2 µg/ml for IP. 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 Rabbit Monoclonal Antibody - Biotinylated (Clone RM305) at a concentration of 0.1 µg/ml.



Immunohistochemical staining of formalin-fixed and paraffin-embedded LNCaP cells overexpressing HA-tag Bag1 protein using HA-tag Rabbit Monoclonal Antibody - Biotinylated (Clone RM305) at a concentration of 0.01 µg/ml.

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

HA-tag Rabbit Monoclonal Antibody - Biotinylated 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.¹ 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 Rabbit Monoclonal Antibody - Biotinylated (Clone RM305) can be used for flow cytometry (FC), immunocytochemistry (ICC), immunohistochemistry (IHC), immunoprecipitation (IP), 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