

PRODUCT INFORMATION



Goat Anti-Mouse IgG (H+L) Biotin

Item No. 44394

Overview and Properties

Contents:	This vial contains 500 µl of biotinylated immunoaffinity-purified IgG
Cross Reactivity:	(+) IgG
Species Reactivity:	(+) Mouse
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	0.01 M sodium phosphate, 0.25 M sodium chloride, pH 7.6, with 15 mg/ml BSA, 50% glycerol, and 0.05% sodium azide
Host:	Goat
Applications:	ELISA and Western blot (WB); the recommended starting dilutions are 1:10,000 for ELISA and 1:4,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

Goat Anti-Mouse IgG (H+L) Biotin is a polyclonal secondary antibody that can be used for Western blot (WB) and ELISA applications, where the primary antibody was generated in mice. This biotin-conjugated secondary antibody is suitable when employing detection using commercially available streptavidin- or avidin-conjugated reagents such as streptavidin-HRP (Item No. 16747), which can be quantified using various substrates, including TMB (Item Nos. 36361 | 70450).

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, 09/08/2025

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