

PRODUCT INFORMATION

Goat Anti-Rabbit IgG - Biotinylated

Item No. 44393

Overview and Properties

Contents:	This vial contains 500 µl of biotinylated immunoaffinity-purified IgG
Immunogen:	Rabbit IgG
Cross Reactivity:	(+) IgG
Species Reactivity:	(+) Rabbit; other species not tested
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 dilution is 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-Rabbit 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 rabbits. This biotin-conjugated secondary antibody is suitable for detection with 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, 09/05/2025

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