

PRODUCT INFORMATION



Goat Anti-Mouse IgG HRP

Item No. 10004302

Overview and Properties

Contents:	This vial contains 500 µl of immunoaffinity-purified IgG conjugated to HRP.
Synonym:	Horseradish Peroxidase-Conjugated Goat Anti-Mouse 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, 15 mg/ml BSA, and 50% glycerol
Host:	Goat
Applications:	ELISA and Western blot (WB); the recommended starting dilution is 1:10,000 and 1:4,000, respectively. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

This goat anti-mouse IgG HRP is used as the "secondary antibody" for Western blotting or ELISA where the primary antibody was generated in mice. This peroxidase conjugated secondary antibody is suitable for Western blotting when employing detection by commercially available chemiluminescence substrates. Additionally this secondary antibody conjugate can be used to develop ELISAs with substrates like ABTS (2,2'-azino-bis-[3-ethylbenzothiazoline-6-sulfonic acid]-diammonium salt), OPD (O-phenylenediamine), or TMB (3,3',5,5'-tetramethylbenzidine).

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, 10/10/2023

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