

PRODUCT INFORMATION



Goat Anti-Rabbit IgG HRP

Item No. 10004301

Overview and Properties

Contents:	This vial contains 500 µl of immunoaffinity-purified IgG conjugated to HRP.
Synonym:	Horseradish Peroxidase-Conjugated Goat Anti-Rabbit IgG
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 and 50% glycerol
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

This goat anti-rabbit IgG HRP is used as the 'secondary antibody' for western blotting or ELISA where the primary antibody was generated in rabbits. This peroxidase-conjugated secondary antibody is suitable for western blotting when employing detection by commercially available chemiluminescence-based 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