

PRODUCT INFORMATION

Goat Anti-Rabbit IgG (H+L) Alexa Fluor® 647

Item No. 44395

Overview and Properties

Contents:	This vial contains 500 µl of immunoaffinity-purified IgG conjugated to Alexa Fluor® 647.
Immunogen:	Rabbit IgG
Cross Reactivity:	(+) IgG
Species Reactivity:	(+) Rabbit
Form:	Liquid
Ex./Em. Max:	650/668 nm
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	0.01M sodium phosphate, 0.25M sodium chloride, pH 7.6, with 15 mg/ml BSA, 50% glycerol, and 0.05% sodium azide
Host:	Goat
Applications:	Immunofluorescence (IF) and flow cytometry (FC); the recommended starting dilution is 1:500 for IF and FC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

Goat Anti-Rabbit IgG (H+L) Alexa Fluor® 647 is a polyclonal secondary antibody that can be used for flow cytometry (FC) and immunofluorescence (IF) applications, where the primary antibody was generated in rabbits. This antibody is composed of immunoaffinity-purified IgG conjugated to Alexa Fluor® 647, which displays excitation/emission maxima of 650/668 nm, respectively.¹

Reference

1. Suzuki, T., Matsuzaki, T., Hagiwara, H., *et al.* Recent advances in fluorescent labeling techniques for fluorescence microscopy. *Acta Histochem. Cytochem.* **40(5)**, 131-137 (2007).

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, 12/05/2025

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

PRODUCT INFORMATION



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