

PRODUCT INFORMATION



Toll-Like Receptor 2 Monoclonal Antibody (Clone TL2.1)

Item No. 13587

Overview

Contents:	This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
Synonym:	TLR2
Immunogen:	CHO cells transfected with human TLR2 cDNA
Cross Reactivity:	(+) Human and Canine TLR2

Properties

Storage:	4°C (≤6 months); -20°C (≥6 months);
Clone:	TL2.1
Host:	Mouse
Isotype:	IgG2a

Applications

Flow cytometry

(intracellular and cell surface): Suggested starting dilution is 1-2 µg/1x10⁶ cells

Immunoprecipitation: Suggested starting dilution is 1:10-1:500

Immunohistochemistry: Suggested starting dilution is 1-10-1:500

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, 01/24/2023

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