

PRODUCT INFORMATION



Toll-Like Receptor 3 Polyclonal Antibody (Clone 40C1285.6)

Item No. 13588

Overview

Contents: This vial contains 100 µg mg of protein G-purified IgG in 0.2 ml PBS containing 0.05% BSA and 0.05% sodium azide.
Synonym: TLR3
Immunogen: Synthetic peptide from human TLR3 amino acids 55-70 (VLNLTHNQLRRLPAAN)
Cross Reactivity: (+) Human and Canine TLR3

Properties

Storage: 4°C (≤6 months); -20°C (≥6 months);
Clone: 40C1285.6
Host: Mouse
Isotype: IgG_{1κ}

Applications

Western blot: Suggested starting dilution is 1-3 µg/ml
Flow cytometry
(intracellular and cell surface): Suggested starting dilution is 2-4 µg/1x10⁶ cells
Immunoprecipitation: Suggested starting dilution is 1:10 - 1:500
Immunohistochemistry: Suggested starting dilution is 2 µg/1x10⁶

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, 07/31/2015

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