

CERTIFICATE of ANALYSIS



Toll-Like Receptor 9 Monoclonal Antibody (Clone 26C593.2)

Item No. 13593 • Batch No. 0430406 • Expires 25JAN2024

Overview

Formulation:	100 µg of protein G-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide
Storage/stability:	-20°C/1yr
Applications:	WB, FC (intracellular and cell surface), IHC (frozen and paraffin-embedded sections), and ICC
Species reactivity:	(+)Human TLR9(+)Canine TLR9(+)Equine TLR9(+)Mouse TLR9(+)Rat TLR9(+)Rhesus monkey TLR9

Reviewed and approved by: Timothy Schrot

Approved date: 25JAN2023



WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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