

CERTIFICATE of ANALYSIS



Chemokine-Like Receptor 1 Polyclonal Antibody

Item No. 10325 • Batch No. 0803226 • Expires 03APR2028

Overview

Formulation:	500 µl of peptide affinity-purified polyclonal antibody
Concentration:	1 mg/mL
Storage/stability:	-20°C/3yr
Applications:	FC, IF, IHC, and WB
Species reactivity:	(+)Human CMKLR1(+)Monkey CMKLR1(+)Mouse CMKLR1(+)Rat CMKLR1

Reviewed and approved by: Beth Hurst

Approved date: 14APR2025

Beth Hurst

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