

Product Information

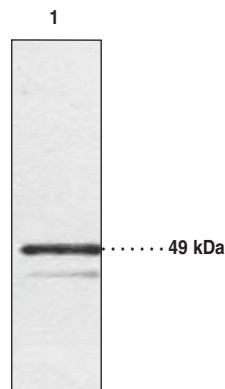


Chemokine-Like Receptor 1 Polyclonal Antibody

Item No. 10325 • Lot No. XXXX

- Contents:** This vial contains (100-500 µg of peptide affinity-purified IgG, *lot specific*) in 500 µl TBS, pH 7.4, containing 50% glycerol, 5 mg/ml BSA, and 0.02% sodium azide. This is sufficient antibody for at least 10 western blots.
- Synonyms:** CMKLR1, DEZ, GPCR ChemR23, Resolvin E1 Receptor
- Antigen:** Human CMKLR1 amino acids 358-371 (NERTSMNERETGML)
- Host:** Rabbit
- Cross Reactivity:** (+) Human, monkey, rat, and mouse CMKLR1
- Stability:** ≥1 year at -20°C
- Applications:** Flow cytometry (FC), immunocytochemistry (ICC), and western blot (WB). Recommended starting concentrations: FC - 10.0 µg/ml, ICC - 10-20 µg/ml, and WB - *lot specific:lot specific*. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.
- Concentration:** Varies by lot, from 0.2-1.0 mg/ml (100-500 µg/vial). Always 100 ml final working volume for western blotting.

Chemokine-Like Receptor 1 (CMKLR1) is a G protein-coupled receptor relevant to the cellular chemotaxis of dendritic cells and macrophages.¹ This receptor is also expressed in brain, liver, lung, and kidney tissues.^{1,2} Chemerin, or TIG2, has been identified as the natural ligand for this receptor.³ Resolvin E1 has also been identified as a ligand for CMKLR1; acting to dampen cellular responses to inflammation.^{1,4,5} Chemerin is an 18 kDa protein that plays a role in immunity, inflammation, chemotaxis, and has been identified as an adipokine.^{6,7,2} Human CMKLR1 shares 80% sequence homology with its mouse ortholog.³ The human receptor exists in two isoforms, differing only by two amino acids due to a premature stop codon. Isoform one has 363 amino acids and an expected molecular weight of 42.3 kDa. Cayman's CMKLR1 polyclonal antibody recognizes the C-terminal region of the protein and cell permeabilization is recommended prior to whole cell staining.



Lane 1: COS-1 cell lysate (10 µg)

Laboratory Procedures

Immunofluorescent staining of cultured cells

1. Prepare cell samples according to your standard protocols.
2. Wash briefly with PBS, pH 7.4, to remove excess serum (if cell culture).
3. Fix the cells with 3.5% formaldehyde in PBS, pH 7.4, for 10 minutes.
4. Wash the cells with PBS containing 2% BSA two times, 10 minutes each.
5. Incubate the cells with the antibody (recommended starting concentration of 10 µg/ml; the optimal working condition should be determined by titration) for 20-60 minutes at room temperature.
6. Wash the cells with PBS/BSA buffer three times and then incubate with a secondary antibody at a dilution suggested by the manufacturer.
7. Wash the cells three times more and examine the staining by flow cytometry or by fluorescent microscopy with an appropriate filter. Store the plate at 4°C in the dark for later analysis if desired.

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 02/03/2012

Product Information



References

1. Arita, M., Bianchini, F., Aliberti, J., *et al.* Stereochemical assignment, anti-inflammatory properties, and receptor for the omega-3 lipid mediator resolvin E1. *J. Exp. Med.* **201**(5), 713-722 (2005).
2. Takahashi, M., Takahashi, Y., Takahashi, K., *et al.* Chemerin enhances insulin signaling and potentiates insulin-stimulated glucose uptake in 3T3-L1 adipocytes. *FEBS Lett.* **582**, 573-578 (2008).
3. Wittamer, V., Grégoire, F., Robberecht, P., *et al.* The C-terminal nonapeptide of mature chemerin activates the chemerin receptor with low nanomolar potency. *Journal of Biological Chemistry* **279**(11), 9956-9962 (2004).
4. Arita, M., Ohira, T., Sun, Y.-P., *et al.* Resolvin E1 selectively interacts with leukotriene B₄ receptor BLT₁ and ChemR23 to regulate inflammation. *J. Immunol.* **178**, 3912-3917 (2007).
5. Cash, J.L., Hart, R., Russ, A., *et al.* Synthetic chemerin-derived peptides suppress inflammation through ChemR23. *JEM* **205**(4), 767-775 (2008).
6. Roh, S.-G., Song, S.-H., Choi, K.-C., *et al.* Chemerin-a new adipokine that modulates adipogenesis *via* its own receptor. *Biochem. Biophys. Res. Commun.* **362**, 1013-1018 (2007).
7. Campbell, E.L., Louis, N.A., Tomassetti, S.E., *et al.* Resolvin E1 promotes mucosal surface clearance of neutrophils: a new paradigm for inflammatory resolution. *FASEB J.* **21**, 3162-3170 (2007).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10325

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com