

# PRODUCT INFORMATION



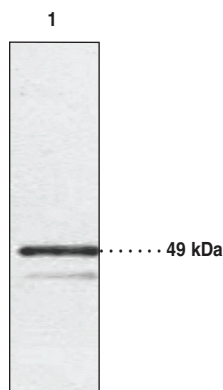
## Chemokine-Like Receptor 1 Polyclonal Antibody

Item No. 10325

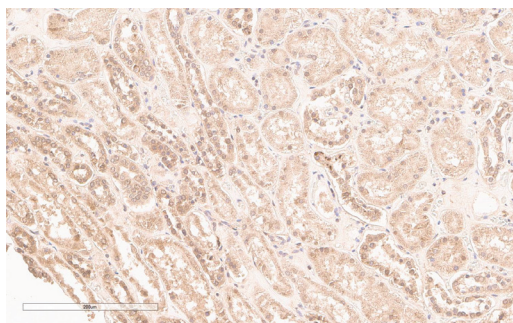
### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide affinity-purified antibody.
<b>Synonyms:</b>	CMKLR1, DEZ, GPCR ChemR23, Resolvin E1 Receptor
<b>Immunogen:</b>	Synthetic peptide from the C-terminal region of human CMKLR1
<b>Species Reactivity:</b>	(+) Human, monkey, mouse, rat
<b>Uniprot No.:</b>	Q99788
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	Flow cytometry (FC), Immunofluorescence (IF), Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1:100 for FC, IF, and WB and 1:200 for IHC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: COS-1 cell lysate (10 µg)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human kidney tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with Chemokine-like Receptor 1 Polyclonal Antibody (Item No. 10325) at a 1:200 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

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, 06/01/2020

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

# PRODUCT INFORMATION



## Description

---

Chemokine-Like Receptor 1 (CMKLR1) is a G protein-coupled receptor relevant to the cellular chemotaxis of dendritic cells and macrophages.<sup>1</sup> This receptor is also expressed in brain, liver, lung, and kidney tissues.<sup>1,2</sup> Chemerin, or TIG2, has been identified as the natural ligand for this receptor.<sup>3</sup> Resolvin E1 has also been identified as a ligand for CMKLR1, acting to dampen cellular responses to inflammation.<sup>1,4,5</sup> Chemerin is an 18 kDa protein that plays a role in immunity, inflammation, chemotaxis, and has been identified as an adipokine.<sup>6,7,2</sup> Human CMKLR1 shares 80% sequence homology with its mouse ortholog.<sup>3</sup> 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.

## 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 B<sub>4</sub> receptor BLT<sub>1</sub> 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).

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