

# PRODUCT INFORMATION



## CD4 Monoclonal Antibody (Clone RIV7)

Item No. 10004599

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µg of protein A-purified IgG.
<b>Immunogen:</b>	Human peripheral lymphocytes
<b>Cross Reactivity:</b>	(+) CD4
<b>Species Reactivity:</b>	(+) Human
<b>Uniprot No.:</b>	P01730
<b>Form:</b>	Liquid
<b>Storage:</b>	4°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS, pH 7.2, with 0.02% sodium azide
<b>Clone:</b>	RIV7
<b>Host:</b>	Mouse
<b>Isotype:</b>	IgG <sub>2a</sub>
<b>Applications:</b>	Flow Cytometry (FC), Immunohistochemistry (IHC; frozen), and Immunofluorescence (IF); the recommended starting dilution for FC and IF is 10 µg/ml and 10 µg/ml (frozen) for IHC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Description

CD4 is a type I transmembrane glycoprotein that functions as a T cell receptor (TCR) co-receptor.<sup>1</sup> It exists as a single polypeptide chain composed of four extracellular immunoglobulin-like (Ig-like) domains that interact with MHC class II molecules, a transmembrane domain, and a cytoplasmic tail that associates with the tyrosine kinase LCK and mediates signal transduction to the TCR, which is essential for T cell activation.<sup>2</sup> It is expressed on the surface of, and used as a marker for, T cells, and its expression is used to characterize the development stage of thymocytes. Upon binding to antigen-displaying MHC class II molecules expressed by antigen-presenting cells (APCs), naïve CD4<sup>+</sup> T cells differentiate and proliferate in a cytokine-dependent manner into a variety of T helper (Th) cell subsets, including Th1, Th2, and Th17 cells, which enhance and direct innate and adaptive immune cell responses to numerous pathogens and have additional roles in cancer, asthma and allergy, and autoimmunity.<sup>3,4</sup> CD4 is also the receptor for HIV attachment and entry into cells, resulting in depletion of CD4<sup>+</sup> cells in patients infected with HIV.<sup>5,6</sup> Cayman's CD4 Monoclonal Antibody (Clone RIV7) can be used for flow cytometry (FC), immunocytochemistry (ICC), and immunohistochemistry (IHC; frozen) applications.

### References

1. Wittlich, M., Koenig, B.W., Hoffmann, S., *et al.* Structural characterization of the transmembrane and cytoplasmic domains of human CD4. *Biochim. Biophys. Acta* **1768**(12), 2949-2960 (2007).
2. Mak, T.W. and Saunders, M.E. The T cell receptor: Structure of its proteins and genes. *The immune response: Basic and clinical principles*. Picknett, T. and Lebedeva, V., editors, 1<sup>st</sup> edition, Elsevier Academic Press (2006).
3. Nguyen, Q.P., Deng, T.Z., Witherden, D.A., *et al.* Origins of CD4<sup>+</sup> circulating and tissue-resident memory T-cells. *Immunology* **157**(1), 3-12 (2019).
4. Zhu, J. and Paul, W.E. CD4 T cells: Fates, functions, and faults. *Blood* **112**(5), 1557-1569 (2008).
5. Wilen, C.B., Tilton, J.C., and Doms, R.W. HIV: Cell binding and entry. *Cold Spring Harb. Perspect. Med.* **2**(8), a006866 (2012).
6. Vijayan, K.K.V., Karthigeyan, K.P., Tripathi, S.P., *et al.* Pathophysiology of CD4<sup>+</sup> T-cell depletion in HIV-1 and HIV-2 infections. *Front. Immunol.* **8**, 580 (2017).

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, 08/15/2022

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