

# PRODUCT INFORMATION



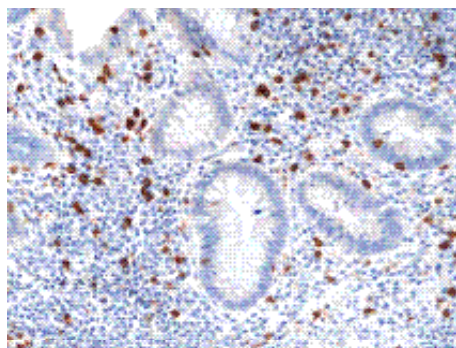
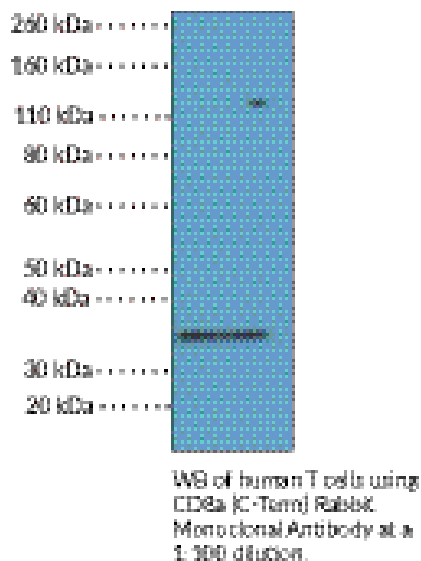
## CD8a (C-Term) Rabbit Monoclonal Antibody

Item No. 32319

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
<b>Synonyms:</b>	T Cell Surface Glycoprotein CD8 α-chain, T Cell Surface Glycoprotein CD8 Alpha Chain
<b>Immunogen:</b>	Peptide from the C-terminal region of human CD8
<b>Cross Reactivity:</b>	(+) CD8a
<b>Species Reactivity:</b>	(+) Human
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Clone:</b>	RM397
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC and WB is 1:100-1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Immunohistochemical staining of formalin-fixed and paraffin-embedded human appendix tissue using CD8a (C-Term) Rabbit Monoclonal Antibody at a 1:100 dilution.

**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, 03/18/2021

**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

---

The CD8  $\alpha$ -chain (CD8a) is a transmembrane glycoprotein that exists as a homodimer, CD8aa, or heterodimer with the CD8  $\beta$ -chain (CD8b), CD8a $\beta$ , to form the T cell receptor (TCR) co-receptor CD8.<sup>1,2</sup> CD8a is composed of an extracellular immunoglobulin variable-like (IgV) domain, which interacts with MHC class I molecules, a transmembrane domain, and a cytoplasmic tail that associates with the tyrosine kinase p56<sup>lck</sup> and mediates signal transduction.<sup>1-3</sup> CD8a $\beta$  is expressed by cytotoxic T lymphocytes (CTLs), where it acts as a TCR co-receptor by enhancing T cell signaling and activation upon binding to antigen-displaying MHC class I molecules.<sup>2,4</sup> CD8aa is expressed by CTLs and intestinal epithelial lymphocytes, as well as  $\gamma\delta$  T, natural killer (NK), and dendritic cells, where it acts as a positive or negative regulator of cell activation in a cell-dependent manner. *Cd8a*<sup>-/-</sup> mice infected with *M. tuberculosis* exhibit reduced survival.<sup>5</sup> Cayman's CD8a (C-Term) Rabbit Monoclonal Antibody can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

## References

---

1. Laugel, B., Cole, D.K., Clement, M., *et al.* The multiple roles of the CD8 coreceptor in T cell biology: Opportunities for the selective modulation of self-reactive cytotoxic T cells. *J. Leukoc. Biol.* **90(6)**, 1089-1099 (2011).
2. Cheroutre, H. and Lambolez, F. Doubting the TCR coreceptor function of CD8aa. *Immunity* **28(2)**, 149-159 (2008).
3. Irie, H.Y., Ravichandran, K.S., and Burakoff, S.J. CD8 $\beta$  chain influences CD8a chain-associated Lck kinase activity. *J. Exp. Med.* **181(4)**, 1267-1273 (1995).
4. Geng, J. and Raghavan, M. CD8aa homodimers function as a coreceptor for KIR3DL1. *Proc. Nat. Acad. Sci. U.S.A.* **115(36)**, 17951-17956 (2019).
5. Sousa, A.O., Mazzaccaro, R.J., Russell, R.G., *et al.* Relative contributions of distinct MHC class I-dependent cell populations in protection to tuberculosis infection in mice. *Proc. Nat. Acad. Sci. U.S.A.* **97(8)**, 4204-4208 (2000).

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