

PRODUCT INFORMATION



FcεRII/CD23 (C-Term) Rabbit Monoclonal Antibody (RM406)

Item No. 32332

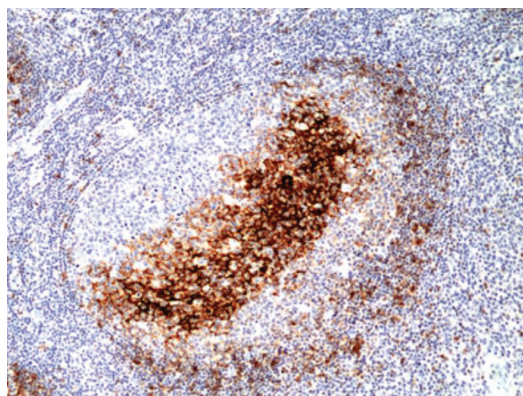
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	Fc Fragment of IgE Receptor II, FcεRII, Low-affinity Immunoglobulin ε Fc Receptor
Immunogen:	Peptide from the C-terminal region of human CD23
Cross Reactivity:	(+) CD23
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM406
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for is 1:100-1:200 for IHC and 1:500-1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



WB of Raji cell lysate using FcεRII/CD23 (C-Term) Rabbit Monoclonal Antibody (RM406) at a dilution of 1:500.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human tonsil tissue using FcεRII/CD23 (C-Term) Rabbit Monoclonal Antibody (RM406) at a dilution of 1:200.

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, 02/10/2024

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

CD23, also known as FcεRII, is a type II transmembrane glycoprotein and C-type lectin encoded by *FCER2* in humans.¹ It is composed of a short N-terminal cytoplasmic domain, a membrane-spanning region, a stalk region, and a large C-terminal globular domain. CD23 is expressed in T and B cells, polymorphonuclear leukocytes, bone marrow stromal cells, follicular dendritic cells, and intestinal epithelial cells. Alternative splicing of *FCER2* pre-mRNA produces two isoforms, CD23a and CD23b, which differ by seven amino acids in the N-terminal cytoplasmic domain and exhibit distinct functions, as well as different uptake and recycling patterns. CD23 is the low-affinity receptor for IgE and has roles in the growth and development of normal and leukemic B cells, antimicrobial immunity, and monocyte activation.^{1,2} It can also be released from cell surfaces by the metalloprotease ADAM10 as various soluble CD23 proteins (sCD23), all of which bind IgE and have cytokine-like activity.¹ Serum levels of CD23 are increased in patients with chronic lymphocytic leukemia (CLL) or hairy cell leukemia.³ Cayman's FcεRII/CD23 (C-Term) Rabbit Monoclonal Antibody (RM406) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Acharya, M., Borland, G., Edkins, A.L., *et al.* CD23/FcεRII: Molecular multi-tasking. *Clin. Exp. Immunol.* **162(1)**, 12-23 (2010).
2. Engeroff, P. and Vogel, M. The role of CD23 in the regulation of allergic responses. *Allergy* (2020).
3. Gordon, J. CD23: Novel disease marker with a split personality. *Clin. Exp. Immunol.* **86(3)**, 356-359 (1991).

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