

PRODUCT INFORMATION



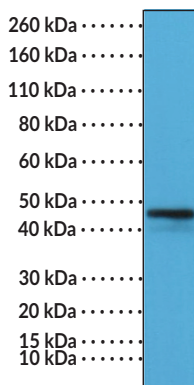
CD79a (C-Term) Rabbit Monoclonal Antibody (Clone RM297)

Item No. 32243

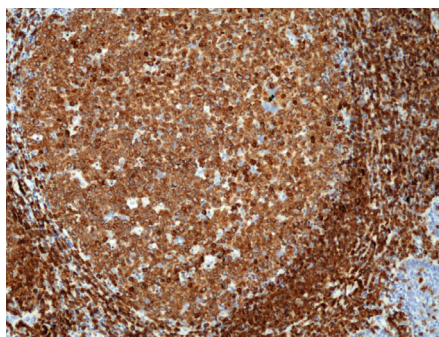
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	B Cell Antigen Receptor Complex-associated Protein α Chain, Igα, MB-1 Membrane Glycoprotein
Immunogen:	Peptide from the C-terminal region of human CD79a
Cross Reactivity:	(+) CD79a
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:	RM297
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



WB of Raji cell lysates using CD79a (C-Term) Rabbit Monoclonal Antibody (Clone RM297) at a 1:200 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human tonsil tissue using CD79a (C-Term) Rabbit Monoclonal Antibody (Clone RM297) at a 1:200 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, 01/30/2024

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

CD79a is a member of the heterodimeric signal transduction component of the B cell receptor (BCR) complex that is encoded by *CD79A/MB-1* in humans.¹ It is composed of a disulfide-linked extracellular domain, a transmembrane region, and a cytoplasmic tail containing an immunoreceptor tyrosine-based activation motif that facilitates intracellular signaling.^{1,2} CD79a is expressed exclusively in B cells and B cell progenitors, and is commonly used as a B cell marker. It is co-expressed at the cell surface with CD79b and immunoglobulins of all isotypes as the mature BCR complex.¹ Mutations in *CD79A/MB-1* have been found in patients with autosomal recessive agammaglobulinemia.³ Cayman's CD79a (C-Term) Rabbit Monoclonal Antibody (Clone RM297) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Schroeder, H.W., Jr., Imboden, J.B., and Torres, R.M. Antigen receptor genes, gene products, and co-receptors. *Clinical immunology: Principles and practice*. Fleisher, T.A., editor, 4th edition, Saunders (2013).
2. Wang, L.D. and Clark, M.R. B-cell antigen-receptor signalling in lymphocyte development. *Immunology* **110(4)**, 411-420 (2003).
3. Khalili, A., Plebani, A., Vitali, M., *et al.* Autosomal recessive agammaglobulinemia: A novel non-sense mutation in *CD79a*. *J. Clin. Immunol.* **34(2)**, 138-141 (2014).

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