

PRODUCT INFORMATION



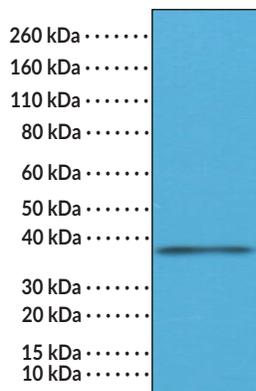
Integrin α 4/CD49D (C-Term) Rabbit Monoclonal Antibody (Clone RM268)

Item No. 32218

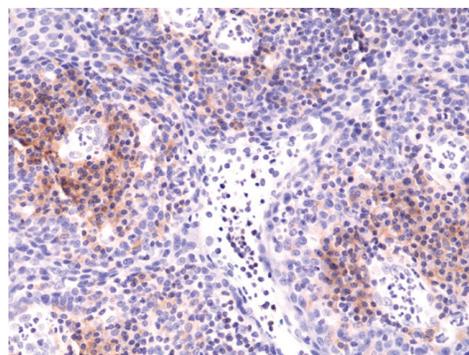
Overview and Properties

Contents:	This vial contains 100 μ l of protein A-affinity purified monoclonal antibody.
Synonyms:	CD49D, Integrin Subunit α 4, Very Late Activation Protein 4 Receptor, α 4 subunit, VLA-4 subunit α , Integrin α IV
Immunogen:	Peptide from the C-terminal region of human integrin α 4
Cross Reactivity:	(+) Integrin α 4
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	\geq 1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM268
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:1,000-1:5,000 for IHC and 1:1,000-1:2,000 for WB.

Images



WB of Jurkat cell lysates using Integrin α 4/CD49D (C-Term) Rabbit Monoclonal Antibody (Clone RM268) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human tonsil tissue using Integrin α 4/CD49D (C-Term) Rabbit Monoclonal Antibody (Clone RM268) at a dilution of 1:5,000.

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, 02/05/2024

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

Integrins are cell adhesion receptors that mediate the attachment of cells to the extracellular matrix (ECM) and facilitate cell-cell interactions.¹ The integrin family contains 18 α subunits and eight β subunits that form 24 different heterodimers grouped into familial subgroups based on subunit composition and/or their ligand binding properties. Integrin $\alpha 4$, also known as CD49D, is a subunit of the $\alpha 4\beta 1$ and $\alpha 4\beta 7$ leukocyte-specific integrins, which recognize fibronectin and vascular cell adhesion molecule 1 (VCAM-1) and have essential roles in lymphocyte migration.^{1,2} Knockout of *Itga4*, the gene encoding integrin $\alpha 4$, induces failure of chorioallantoic fusion and is embryonic lethal in mice.¹ Cayman's Integrin $\alpha 4$ /CD49D (C-Term) Rabbit Monoclonal Antibody (Clone RM268) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Barczyk, M., Carracedo, S., and Gullberg, D. Integrins. *Cell Tissue Res.* **339**(1), 269-280 (2010).
2. Gu, Z.-C., Zhang, S.-G., and Wei, J.-F. $\alpha 4\beta 7$ integrin inhibitors: A patent review. *Expert Opin. Ther. Pat.* **28**(12), 903-917 (2018).

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