

PRODUCT INFORMATION



Aminopeptidase N/CD13 (C-Term) Rabbit Monoclonal Antibody (RM403) Item No. 3232

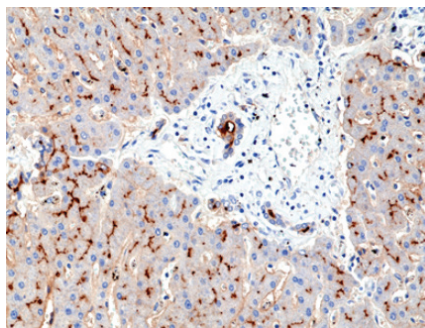
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	Alanine Aminopeptidase, APN, ANPEP, gp150, Microsomal Aminopeptidase, Particle-bound Aminopeptidase
Immunogen:	Peptide from the C-terminal region of human aminopeptidase N
Cross Reactivity:	(+) Aminopeptidase N
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:	RM403
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:100-1:200 and 1:1,000-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



WB of HepG2 cells using Aminopeptidase N/CD13 (C-Term) Rabbit Monoclonal Antibody (RM403) at a 1:2,000 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human liver tissue using Aminopeptidase N/CD13 (C-Term) Rabbit Monoclonal Antibody (RM403) 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, 02/21/2024

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

Aminopeptidase N, also known as cluster of differentiation 13 (CD13), is a zinc-dependent and membrane-bound ectopeptidase and type II membrane glycoprotein encoded by *ANPEP* in humans.^{1,2} It is composed of a short N-terminal cytosolic domain, a membrane-spanning domain, and five extracellular domains, including a C-terminal catalytic domain.^{1,3} It is ubiquitously expressed on cell membranes where it functions as a dimer to degrade peptides and proteins containing N-terminal neutral amino acids.¹ Plasma levels of soluble aminopeptidase N are increased and positively correlated with tumor progression in patients with non-small cell lung cancer (NSCLC).² *ANPEP* is overexpressed in psoriatic skin lesions, synovial fibroblasts isolated from patients with rheumatoid arthritis, and lymphocytes and peripheral blood mononuclear cells (PBMCs) isolated from patients with multiple sclerosis during acute exacerbation and chronic progression.¹ Cayman's Aminopeptidase N/CD13 (C-Term) Rabbit Monoclonal Antibody (RM403) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Lu, C., Amin, M.A., and Fox, D.A. CD13/Aminopeptidase N Is a potential therapeutic target for inflammatory disorders. *J. Immunol.* **204**(1), 3-11 (2020).
2. Wickström, K., Larsson, R., Nygren, P., *et al.* Aminopeptidase N (CD13) as a target for cancer chemotherapy. *Cancer Sci.* **102**(3), 501-508 (2011).
3. Sjöström, H., Norén, O., and Olsen, J. Structure and function of aminopeptidase. *Adv. Exp. Med. Biol.* **477**, 25-34 (2000).