

PRODUCT INFORMATION



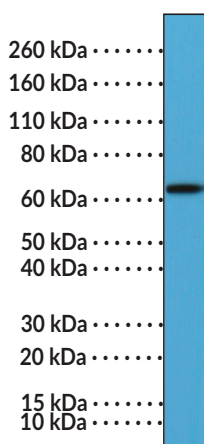
PLAP (N-Term) Rabbit Monoclonal Antibody (Clone RM317)

Item No. 32259

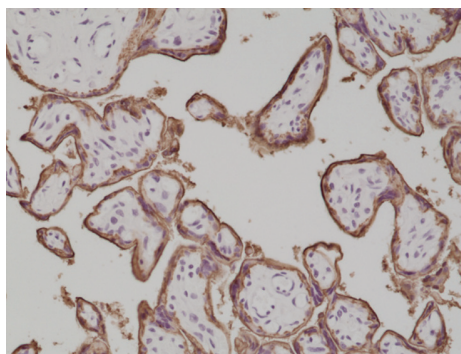
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Immunogen:	Peptide from the N-terminal region of human PLAP
Cross Reactivity:	(+) PLAP
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:	RM317
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:500-1:1,000 for IHC 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 A431 cell lysates using PLAP (N-Term) Rabbit Monoclonal Antibody (Clone RM317) at a dilution of 1:2,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human placenta tissue using PLAP (N-Term) Rabbit Monoclonal Antibody (Clone RM317) at a 1:1,000 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

Placental alkaline phosphatase (PLAP) is the placenta-specific isoenzyme of AP encoded by the *ALPP* gene in humans.^{1,2} It exists as a homodimer where each monomer is composed of an N-terminal α -helix and a crown domain that stabilize PLAP homodimers, a Zn^{2+} - and Mg^{2+} -containing active site with a catalytic serine residue, and a non-catalytic metal-binding site occupied by Ca^{2+} .^{3,4} PLAP is predominately expressed by syncytiotrophoblasts in the placenta and is tethered to the plasma membrane by a glycosylphosphatidylinositol (GPI) anchor.^{3,4} It is also present at low levels in the cervix, ovaries, and type I pneumocytes, as well as in the serum.^{1,3} Increased PLAP activity has been found in tumor tissue and cerebrospinal fluid (CSF) isolated from patients with intracranial germ cell tumors.⁵ Cayman's PLAP (N-Term) Rabbit Monoclonal Antibody (Clone RM317) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Zaher, D.M., El-Gamal, M.I., Omar, H.A., *et al.* Recent advances with alkaline phosphatase isoenzymes and their inhibitors. *Arch. Pharm. (Weinheim)* **353**(5), e2000011 (2020).
2. Wennberg, C., Kozlenkov, A., Di Mauro, S., *et al.* Structure, genomic DNA typing, and kinetic characterization of the D allozyme of placental alkaline phosphatase (PLAP/ALPP). *Hum. Mutat.* **19**(3), 258-267 (2002).
3. Millán, J.L. Alkaline phosphatases : Structure, substrate specificity and functional relatedness to other members of a large superfamily of enzymes. *Purinergic Signal.* **2**(2), 335-341 (2006).
4. Le Du, M.H., Stigbrand, T., Taussig, M.J., *et al.* Crystal structure of alkaline phosphatase from human placenta at 1.8 Å resolution. Implication for a substrate specificity. *J. Biol. Chem.* **276**(12), 9158-9165 (2001).
5. Aihara, Y., Watanabe, S., Amano, K., *et al.* Placental alkaline phosphatase levels in cerebrospinal fluid can have a decisive role in the differential diagnosis of intracranial germ cell tumors. *J. Neurosurg.* **131**(3), 687-694 (2018).