

PRODUCT INFORMATION

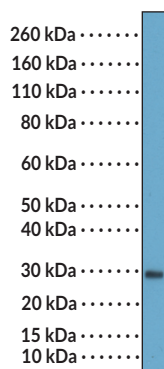
Asprosin Rabbit Monoclonal Antibody (Clone RM463)

Item No. 35811

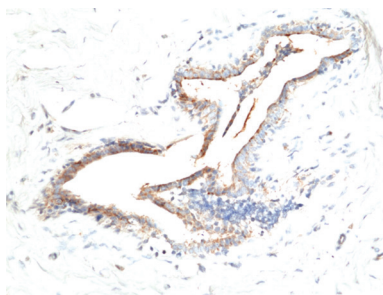
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	Cleaved FBN1, Cleaved Fibrillin-1
Immunogen:	Recombinant full-length human asprosin
Cross Reactivity:	(+) Asprosin
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:	RM463
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200 for IHC and 1:1,000-1:25,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

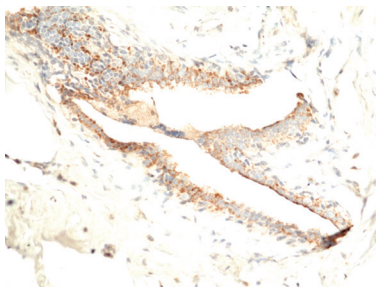
Images



WB of human plasma using Asprosin Rabbit Monoclonal Antibody (Clone RM463) at a dilution of 1:25,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast tissue using Asprosin Rabbit Monoclonal Antibody (Clone RM463) at a dilution of 1:100.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast tissue using Asprosin Rabbit Monoclonal Antibody (Clone RM463) at a dilution of 1:100.

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, 01/29/2029

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

Asprosin is a glucogenic protein hormone secreted by white adipose tissue in mammals.¹ This ~16 kDa protein is a 140-amino acid C-terminal cleavage product produced by furin cleavage of profibrillin. Asprosin increases glucose release *via* cyclic AMP (cAMP) and protein kinase A (PKA) in hepatocytes *via* interaction with a cell surface receptor, an effect that can be reversed by suramin, an inhibitor of heterotrimeric G proteins, or RP-cAMPS (Item No. 16985), an antagonist of cAMP binding to PKA. *In vivo*, elevated plasma asprosin increases hepatic glucose production but does not affect glucose uptake in peripheral organs in response to insulin. Asprosin is orexigenic, activating AgRP⁺ neurons in a cAMP-dependent manner, which leads to appetite stimulation and increased body weight.² Plasma asprosin is elevated in adult humans and mice with type 2 diabetes mellitus and other genetic conditions that confer insulin resistance.^{1,3} Cayman's Asprosin Rabbit Monoclonal Antibody (Clone RM463) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Romere, C., Duerschmid, C., Bournat, J., *et al.* Asprosin, a fasting-induced glucogenic protein hormone. *Cell* **165**(3), 566-579 (2016).
2. Duerschmid, C., He, Y., Wang, C., *et al.* Asprosin is a centrally acting orexigenic hormone. *Nat. Med.* **23**(12), 1444-1453 (2017).
3. Zhang, L., Chen, C., Zhou, N., *et al.* Circulating asprosin concentrations are increased in type 2 diabetes mellitus and independently associated with fasting glucose and triglyceride. *Clin. Chim. Acta.* **489**, 183-188 (2017).

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