

PRODUCT INFORMATION



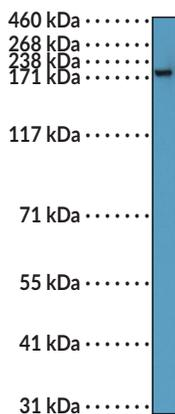
CPS1 (N-Term) Rabbit Monoclonal Antibody (RM395)

Item No. 32320

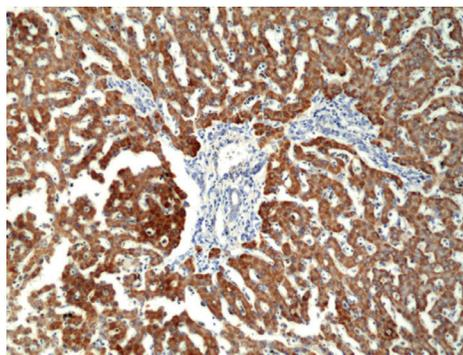
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	Carbamoyl Phosphate Synthase, mitochondrial, CPSase I
Immunogen:	Peptide from the N-terminal region of CPS1
Cross Reactivity:	(+) CPS1
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:	RM395
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200 for IHC and 1:5,000-1:10,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



WB of human T cell lysate using CPS1 (N-Term) Rabbit Monoclonal Antibody (RM395) at a dilution of 1:10,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human liver tissue using CPS1 (N-Term) Rabbit Monoclonal Antibody (RM395) 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, 02/10/2024

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

Carbamoyl phosphate synthetase I (CPS1) is a ligase involved in the urea cycle.¹ It is composed of an N-terminal domain containing an ancestral inactive glutaminase, a bicarbonate phosphorylation site, and a C-terminal domain containing a carbamate phosphorylation site and the binding site for the CPS1 allosteric activator N-acetyl-L-glutamate (NAG). It is primarily expressed in the liver and intestinal epithelial cells and is localized to the mitochondrial matrix.^{1,2} CPS1, with NAG as a cofactor, catalyzes the formation of carbamoyl phosphate from ammonia, bicarbonate, and ATP as the first, rate-limiting step in the urea cycle. Mutations in *CPS1* result in CPS1 deficiency, a disorder characterized by low levels of active CPS1, arginine, and citrulline and high levels of ammonia and glutamine, leading to neurological deficits and, potentially, death.^{3,4} Protein levels of CPS1 are increased in patient-derived rectal cancer tumors and associated with shorter disease-specific survival and metastasis-free survival.⁵ Cayman's CPS1 (N-Term) Rabbit Monoclonal Antibody (RM395) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Martínez, A.I., Pérez-Arellano, I., Pekkala, S., *et al.* Genetic, structural and biochemical basis of carbamoyl phosphate synthetase 1 deficiency. *Mol. Genet. Metab.* **101(4)**, 311-323 (2010).
2. Ryall, J.G., Nguyen, M., Bendayan, M., *et al.* Expression of nuclear genes encoding the urea cycle enzymes, carbamoyl-phosphate synthetase I and ornithine carbamoyl transferase, in rat liver and intestinal mucosa. *Eur. J. Biochem.* **152(2)**, 287-292 (1985).
3. Nitzahn, M. and Lipshutz, G.S. CPS1: Looking at an ancient enzyme in a modern light. *Mol. Genet. Metab.* **131(3)**, 289-298 (2020).
4. Pekkala, S., Martínez, A.I., Barcelona, B., *et al.* Understanding carbamoyl-phosphate synthetase I (CPS1) deficiency by using expression studies and structure-based analysis. *Hum. Mutat.* **31(7)**, 801-808 (2010).
5. Lee, Y.-Y., Li, C.-F., Lin, C.-Y., *et al.* Overexpression of CPS1 is an independent negative prognosticator in rectal cancers receiving concurrent chemoradiotherapy. *Tumour Biol.* **35(11)**, 11097-11105 (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