

PRODUCT INFORMATION



PINK1 Polyclonal Antibody

Item No. 10006283

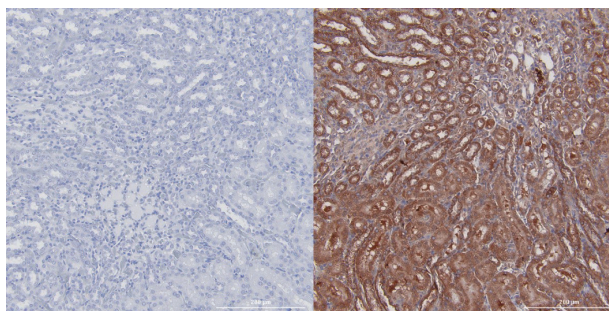
Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified antibody.
Synonyms:	BRPK, PARK6, PTEN Induced Putative Kinase 1
Immunogen:	Synthetic peptide from the C-terminal region of human PINK1
Species Reactivity:	(+) Human, mouse, and rat; other species not tested
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:100 and 1:200 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

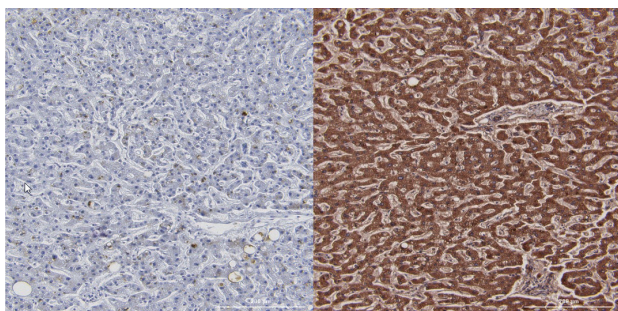
Images



Lane 1: Mouse liver 100,000 x g pellet (30 µg)



Immunohistochemical staining of formalin-fixed paraffin-embedded (FFPE) rat kidney tissue after heat induced antigen retrieval in citrate buffer, pH 6.0, after incubation with PINK1 Polyclonal Antibody (Item no. 10006283) at a 1:100 dilution (left panel, secondary alone).



Immunohistochemical staining of formalin-fixed paraffin-embedded (FFPE) human liver tissue after heat induced antigen retrieval in citrate buffer, pH 6.0, after incubation with PINK1 Polyclonal Antibody (Item no. 10006283) at a 1:100 dilution (left panel, secondary alone).

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, 12/11/2023

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

Phosphatase and tensin homolog (PTEN) dephosphorylates lipids such as phosphatidylinositol 3,4,5-triphosphate (PIP₃) and its defects contribute to a variety of human cancers. PTEN induced putative kinase 1 (PINK1) was first identified when studying the tumor-suppressive function of the PTEN signaling pathway and is thus believed to be involved in human cancer pathology.¹ The mRNA of PINK1 is expressed ubiquitously among adult tissues with most abundant expression in the heart, skeletal muscle, and testis.¹ PINK1 is located in mitochondria and its homozygous C-terminal mutation is associated with early onset of Parkinson's disease.²

Cayman's PINK1 Polyclonal Antibody recognizes primarily the full length protein at about 66 kDa in human, mouse, and rat tissues. In addition, a truncated form of the protein at about 33 kDa is also detected.³

References

1. Unoki, M. and Nakamura, Y. Growth-suppressive effects of BPOZ and EGR2, two genes involved in the PTEN signaling pathway. *Oncogene* **20**, 4457-4465 (2001).
2. Valente, E.M., Abou-Sleiman, P.M., Caputo, V., *et al.* Hereditary early-onset Parkinson's disease caused by mutations in PINK1. *Science* **304**, 1158-1160 (2004).
3. Mammalian Gene Collection (MGC) Program Team Generation and initial analysis of more than 15,000 full-length human and mouse cDNA sequences. *Proc. Natl. Acad. Sci. USA* **99(26)**, 16899-16903 (2002).

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