
OMA1 Polyclonal Antibody

(Catalog #A67094)

Background

Metalloprotease that is part of the quality control system in the inner membrane of mitochondria. Following stress conditions that induce loss of mitochondrial membrane potential, mediates cleavage of OPA1 at S1 position, leading to OPA1 inactivation and negative regulation of mitochondrial fusion. May also cleave UQC3 under these conditions. Its role in mitochondrial quality control is essential for regulating lipid metabolism as well as to maintain body temperature and energy expenditure under cold-stress conditions.

Description

OMA1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4

Specificity

Human

Isotype

IgG

Uniprot ID

Q96E52

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Metalloendopeptidase OMA1, mitochondrial protein (22-194AA)

Storage

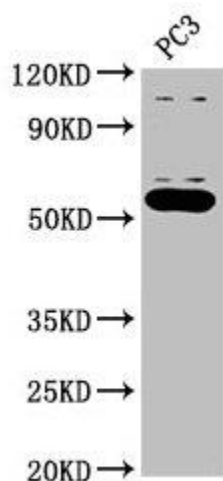
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Metalloendopeptidase OMA1, mitochondrial, Metalloprotease-related protein 1, MPRP-1, Overlapping with the m-AAA protease 1 homolog, OMA1, MPRP1

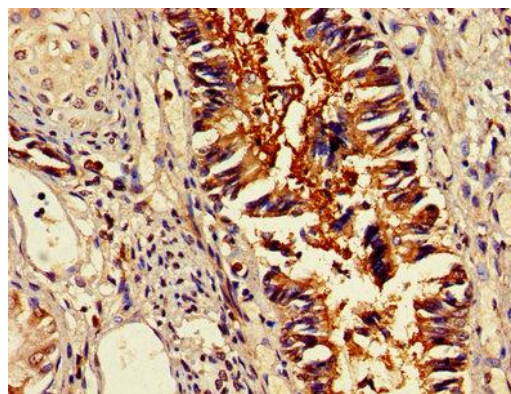
Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:200-1:500, IF:1:50-1:200

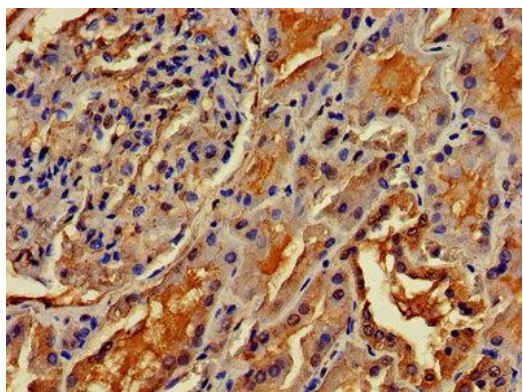


Western Blot

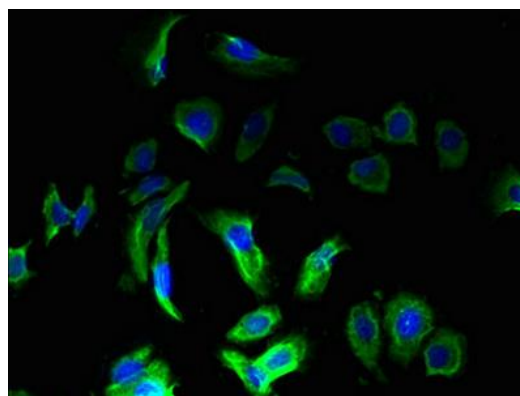
Positive WB detected in: PC-3 whole cell lysate
 All lanes: OMA1 Polyclonal Antibody at 2.7ug/ml
 Secondary
 Goat polyclonal to rabbit IgG at 1/50000 dilution
 Predicted band size: 61, 56 kDa
 Observed band size: 56 kDa



IHC image of OMA1 Polyclonal Antibody diluted at 1:400 and staining in paraffin-embedded human lung tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



IHC image of OMA1 Polyclonal Antibody diluted at 1:400 and staining in paraffin-embedded human kidney tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Immunofluorescence staining of MCF-7 cells with OMA1 Polyclonal Antibody at 1:133, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).