

USP30 Polyclonal Antibody

(Catalog # A61694)

Background

Deubiquitinating enzyme tethered to the mitochondrial outer membrane that acts as a key inhibitor of mitophagy by counteracting the action of parkin (PARK2): hydrolyzes ubiquitin attached by parkin on target proteins, such as RHOT1/MIRO1 and TOMM20, thereby blocking parkin's ability to drive mitophagy (PubMed:18287522, PubMed:24896179, PubMed:25527291, PubMed:25621951). Preferentially cleaves 'Lys-6'- and 'Lys-11'-linked polyubiquitin chains, 2 types of linkage that participate to mitophagic signaling (PubMed:25621951). Does not cleave efficiently polyubiquitin phosphorylated at 'Ser-65' (PubMed:25527291). Acts as negative regulator of mitochondrial fusion by mediating deubiquitination of MFN1 and MFN2.

Description

USP30 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q70CQ3

Purification

Protein G purified

Immunogen

Recombinant Human Ubiquitin carboxyl-terminal hydrolase 30 protein (101-400AA)

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

Ubiquitin carboxyl-terminal hydrolase 30 (EC:3.4.19.121 Publication, USP30, Deubiquitinating enzyme 30, Ubiquitin thioesterase 30, Ubiquitin-specific-processing protease 30, Ub-specific protease 30

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200



Western blot

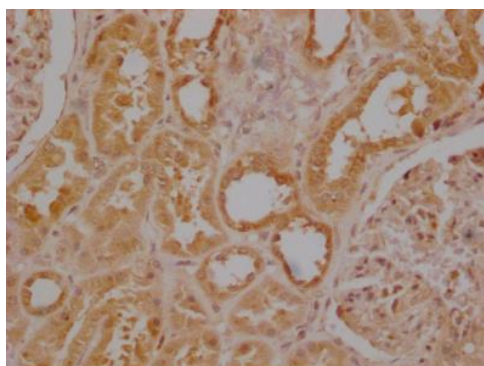
All lanes: USP30 antibody at 7ug/ml + Mouse liver tissue

Secondary

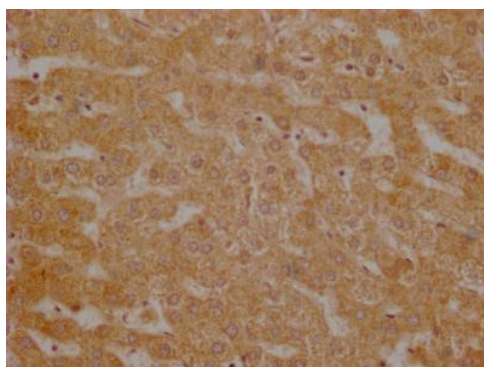
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 59 kDa

Observed band size: 59 kDa



Immunohistochemistry of paraffin-embedded human colon cancer using USP30 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human breast cancer using USP30 Antibody at dilution of 1:100