

ACTRT1 Polyclonal Antibody

(Catalog # A57994)

Background

mitophagy in response to mitochondrial depolarization, positive regulation of defense response to virus by host, xenophagy

Description

ACTRT1 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

Q8TDG2

Purification

Protein G purified

Immunogen

Recombinant Human Actin-related protein T1 protein (201-376AA)

Storage

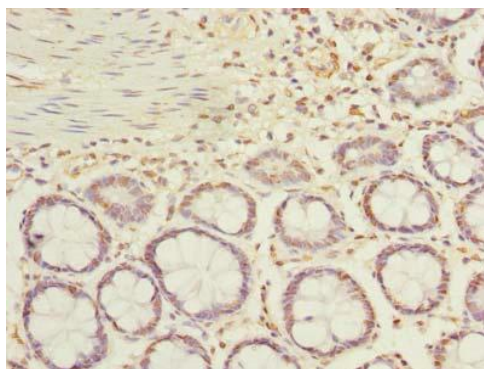
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

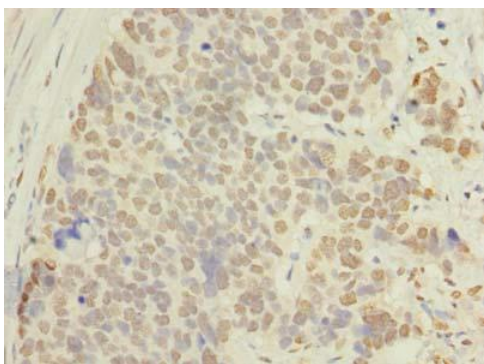
Actin-related protein T1, ARP-T1, ACTRT1, ARPT1, HSD27

Application

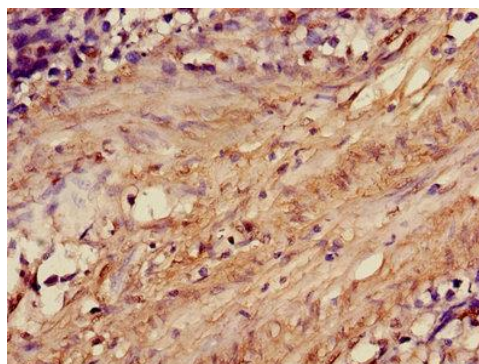
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



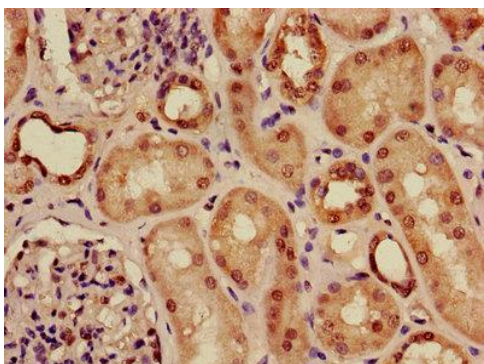
Immunohistochemistry of paraffin-embedded human rectum tissue at dilution of 1:100. Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 90s.



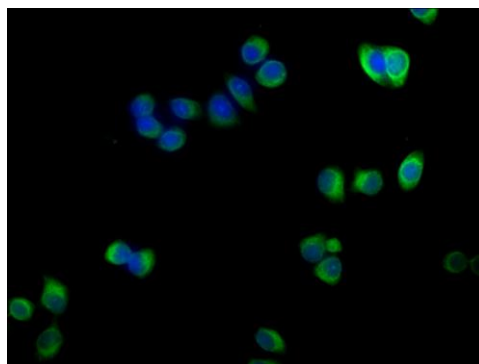
Immunohistochemistry of paraffin-embedded human gastric cancer at dilution of 1:100



Immunohistochemistry of paraffin-embedded human smooth muscle tissue at dilution of 1:100



Immunohistochemistry of paraffin-embedded human kidney tissue at dilution of 1:100



Immunofluorescence staining of PC-3 cells with ACTRT1 Polyclonal Antibody at 1:100, 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).