
ENTPD8 Polyclonal Antibody

(Catalog # A62482)

Background

Canalicular ectonucleoside NTPDase responsible for the main hepatic NTPDase activity. Ectonucleoside NTPDases catalyze the hydrolysis of gamma- and beta-phosphate residues of nucleotides, playing a central role in concentration of extracellular nucleotides. Has activity toward ATP, ADP, UTP and UDP, but not toward AMP.

Description

ENTPD8 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q5MY95

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Ectonucleoside triphosphate diphosphohydrolase 8 protein (30-434AA)

Storage

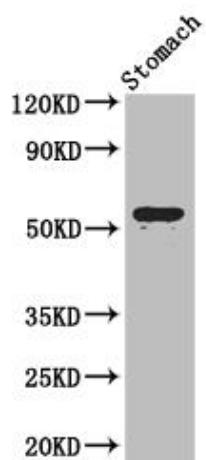
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Ectonucleoside triphosphate diphosphohydrolase 8, E-NTPDase 8, NTPDase 8, NTPDase8, ENTPD8, UNQ2492/PRO5779

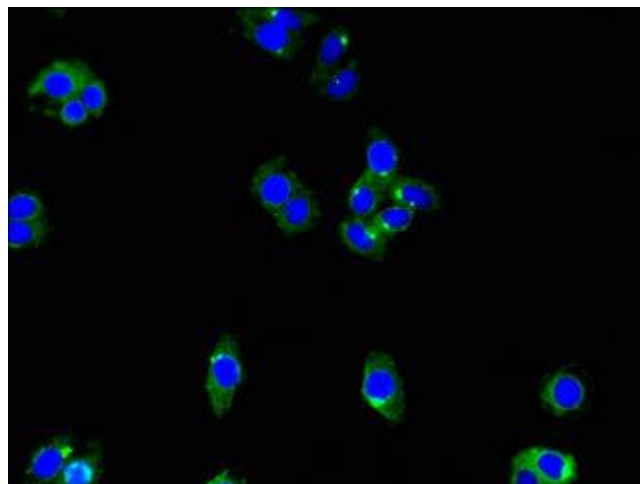
Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:2000-1:5000, IHC:1:20-1:200, IF:1:50-1:200

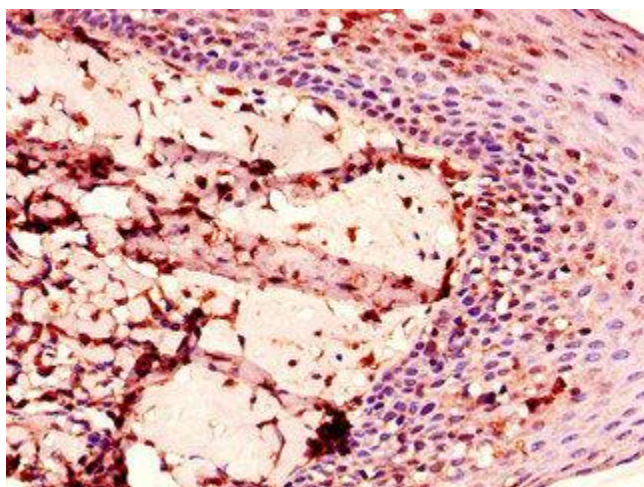


Western Blot

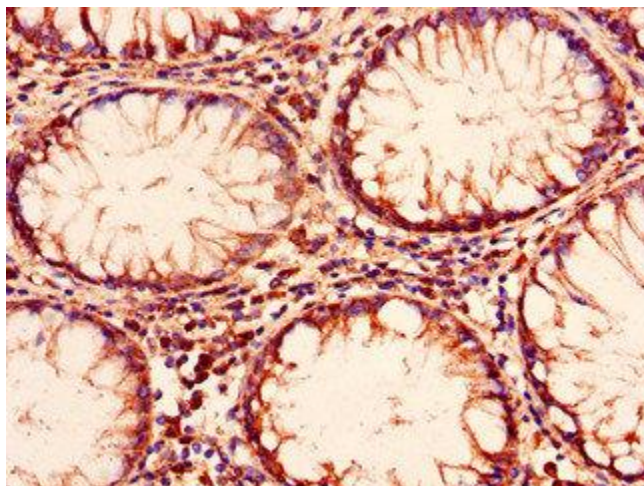
Positive WB detected in: Mouse stomach tissue
 All lanes: ENTPD8 Polyclonal Antibody at 2ug/ml
 Secondary
 Goat polyclonal to rabbit IgG at 1/50000 dilution
 Predicted band size: 54, 50 kDa
 Observed band size: 54 kDa



Immunofluorescent analysis of HepG2 cells at a dilution of 1:100
 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit
 IgG(H+L)



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



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