

C6orf89 Polyclonal Antibody

(Catalog # A65714)

Background

Exhibits histone deacetylase (HDAC) enhancer properties. May play a role in cell cycle progression and wound repair of bronchial epithelial cells.

Description

C6orf89 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q6UWU4

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Bombesin receptor-activated protein C6orf89 protein (120-320AA)

Storage

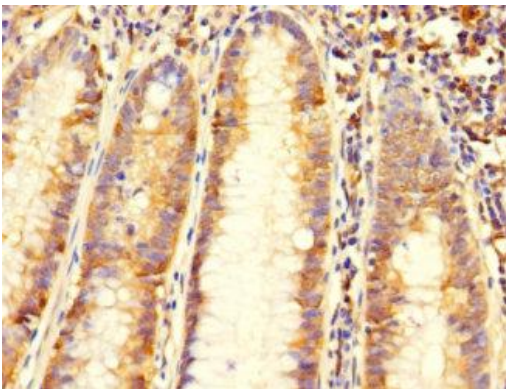
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

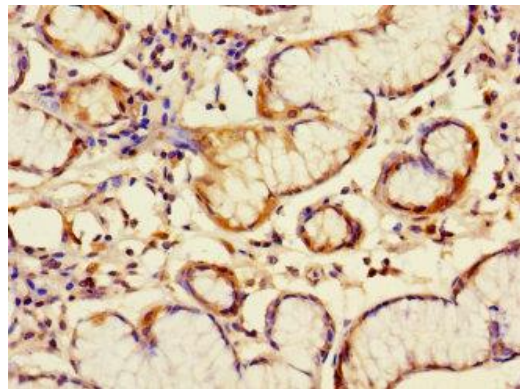
Bombesin receptor-activated protein C6orf89, Amfion, C6orf89, BRAP, UNQ177/PRO203

Application

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



Immunohistochemistry of paraffin-embedded human colon cancer using C6orf89 Polyclonal Antibody at dilution 1:100



Immunohistochemistry of paraffin-embedded human gastric cancer using C6orf89 Polyclonal Antibody at dilution 1:100