

GSDMC Polyclonal Antibody

(Catalog # A67802)

Background

cytoplasm, microtubule organizing center, mitochondrion

Description

GSDMC 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

Q9BYG8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Gasdermin-C protein (1-508AA)

Storage

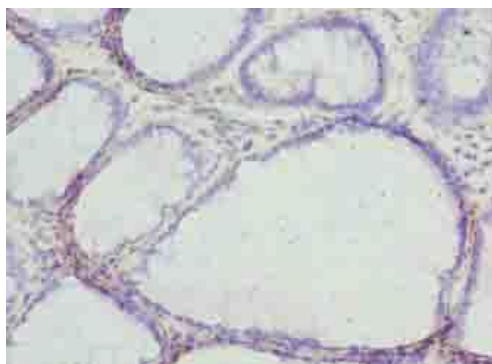
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Gasdermin-C, Melanoma-derived leucine zipper-containing extranuclear factor, GSDMC, MLZE

Application

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



Immunohistochemistry of paraffin-embedded human rectal cancer using GSDMC Polyclonal Antibody at dilution of 1:100