
RNF103 Polyclonal Antibody

(Catalog #A60622)

Background

Acts as an E2-dependent E3 ubiquitin-protein ligase, probably involved in the ER-associated protein degradation pathway.

Description

RNF103 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

O00237

Purification

>95%, Protein G purified

Immunogen

Recombinant Human E3 ubiquitin-protein ligase RNF103 protein (432-685AA)

Storage

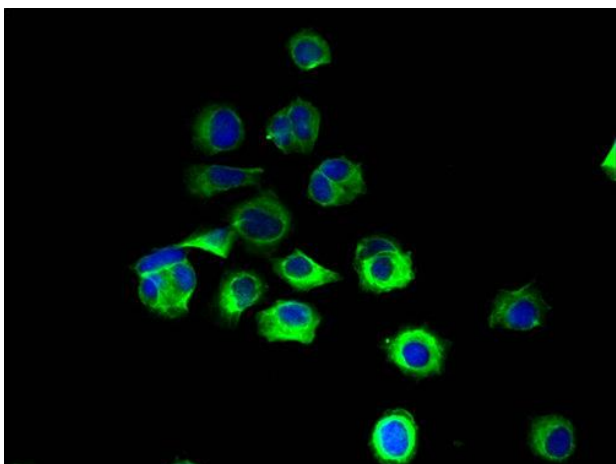
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

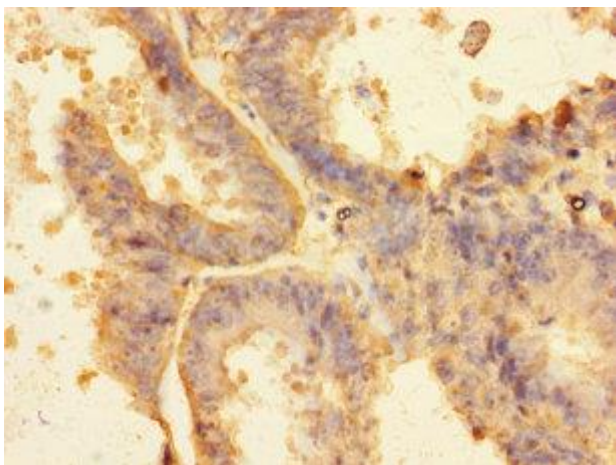
E3 ubiquitin-protein ligase RNF103 (EC:6.3.2.-), RNF103, ZFP103, KF-1, hKF-1, RING finger protein 103, Zinc finger protein 103 homolog, Zfp-103

Application

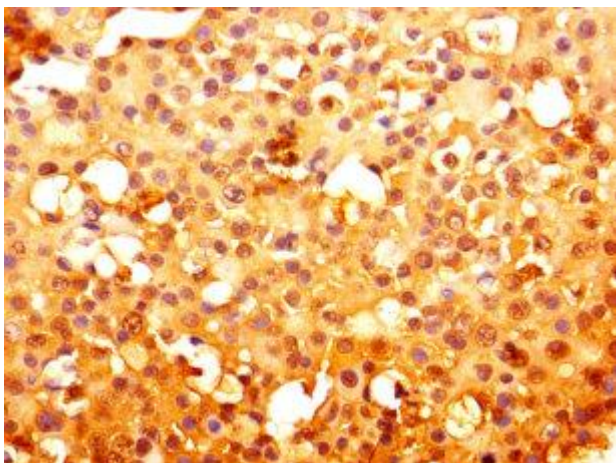
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunofluorescent analysis of MCF-7 cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



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



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