
ZNRF3 Polyclonal Antibody, HRP Conjugated

(Catalog # A68039)

Background

E3 ubiquitin-protein ligase that acts as a negative regulator of the Wnt signaling pathway by mediating the ubiquitination and subsequent degradation of Wnt receptor complex components Frizzled and LRP6. Acts on both canonical and non-canonical Wnt signaling pathway. Acts as a tumor suppressor in the intestinal stem cell zone by inhibiting the Wnt signaling pathway, thereby restricting the size of the intestinal stem cell zone.

Description

ZNRF3 Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q9ULT6

Purification

>95%, Protein G purified

Immunogen

Recombinant Human E3 ubiquitin-protein ligase ZNRF3 protein (56-219AA)

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 ZNRF3, RING finger protein 203, Zinc/RING finger protein 3, ZNRF3, KIAA1133, RNF203

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
