

FGFR1OP2 Polyclonal Antibody

(Catalog # A62590)

Background

May be involved in wound healing pathway.

Description

FGFR1OP2 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

Q9NVK5

Purification

>95%, Protein G purified

Immunogen

Recombinant Human FGFR1 oncogene partner 2 protein (1-172AA)

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

FGFR1 oncogene partner 2, FGFR1OP2, HSPC123

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: FGFR1OP2 Polyclonal Antibody at 6ug/ml

Lane 1: Mouse gonadal tissue

Lane 2: Human placenta tissue

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 30, 25, 21 kDa

Observed band size: 29 kDa