

GFRAL Polyclonal Antibody

(Catalog #A65526)

Description

GFRAL Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q6UXV0

Purification

Protein G purified

Immunogen

Recombinant Human GDNF family receptor alpha-like protein (19-351AA)

Storage

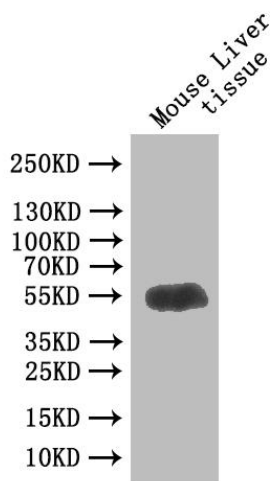
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

GDNF family receptor alpha-like, GFRAL, C6orf144, UNQ9356/PRO34128

Application

WB; Recommended dilution: WB 1:500 - 1:2000



Western Blot

Positive WB detected in: Mouse Liver tissue lysate

All lanes: GFRAL Polyclonal Antibody at 1:1000

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 45 kDa

Observed band size: 55 kDa