

FFAR3 Polyclonal Antibody

(Catalog # A72428)

Description

FFAR3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Mouse, Rat

Isotype

IgG

Uniprot ID

O14843

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 257-346 of human FFAR3 (NP_005295.1).

Storage

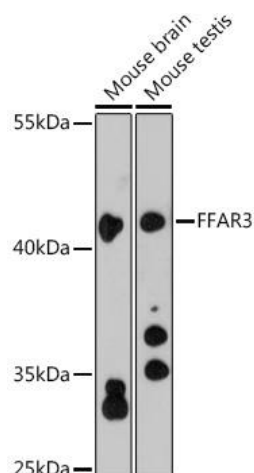
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

FFAR3; FFA3R; GPR41; GPR42; free fatty acid receptor 3

Application

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



Western blot analysis of extracts of various cell lines, using FFAR3 Polyclonal Antibody at 1:3000 dilution.

Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

Exposure time: 60s.