
IPO9 Polyclonal Antibody

(Catalog #A73416)

Background

Enables nuclear import signal receptor activity. Involved in protein import into the nucleus. Located in cytoplasm

Description

IPO9 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thimerosal, 50% glycerol, pH7.3.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q96P70

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 827-1041 of human IPO9 (NP_060555.2)

Storage

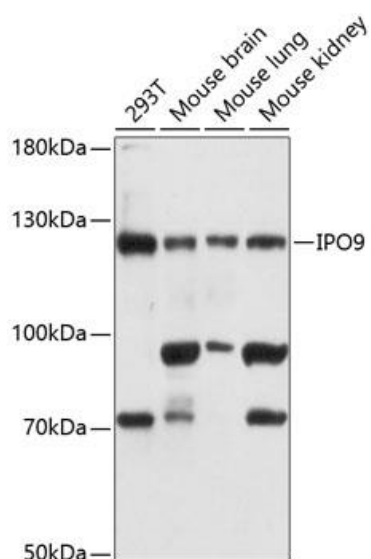
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

IPO9; Imp9

Application

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



Western blot analysis of various lysates using the IPO9 Polyclonal Antibody at 1:3000 dilution.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25µg per lane.
Blocking buffer: 3% nonfat dry milk in TBST.