

LHX2 Polyclonal Antibody

(Catalog # A72966)

Background

This gene encodes a protein belonging to a large protein family, members of which carry the LIM domain, a unique cysteine-rich zinc-binding domain. The encoded protein may function as a transcriptional regulator. The protein can recapitulate or rescue phenotypes in *Drosophila* caused by a related protein, suggesting conservation of function during evolution.

Description

LHX2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P50458

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 337-406 of human LHX2 (NP_004780.3).

Storage

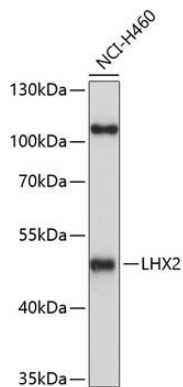
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

LHX2; LH2; hLhx2; LIM homeobox 2

Application

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



Western blot analysis of extracts of H460 cells, using LHX2 Polyclonal Antibody at 1:1000 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: 90s.