
PODXL Polyclonal Antibody

(Catalog #A71496)

Background

This gene encodes a member of the sialomucin protein family. The encoded protein was originally identified as an important component of glomerular podocytes. Podocytes are highly differentiated epithelial cells with interdigitating foot processes covering the outer aspect of the glomerular basement membrane. Other biological activities of the encoded protein include: binding in a membrane protein complex with Na⁺/H⁺ exchanger regulatory factor to intracellular cytoskeletal elements, playing a role in hematopoietic cell differentiation, and being expressed in vascular endothelium cells and binding to L-selectin.

Description

PODXL Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

O00592

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 287-526 of human PODXL (NP_005388.2).

Storage

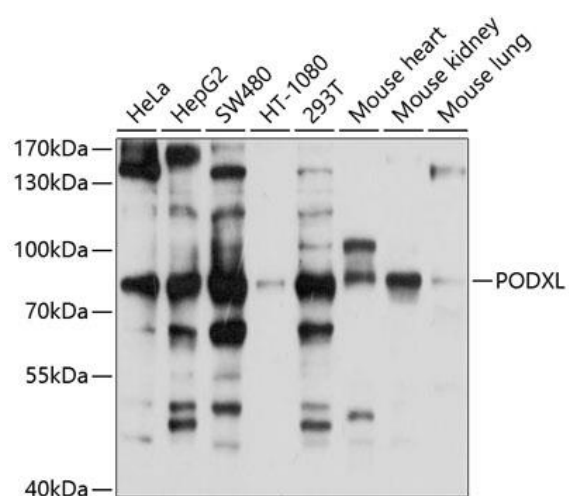
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

PODXL; Gp200; PC; PCLP; PCLP-1; podocalyxin

Application

WB; Recommended dilution: WB 1:1000 - 1:2000



Western blot analysis of extracts of various cell lines,
using PODXL 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: 3s.