

LIN28A antibody (pAb)

Catalog Nos: 61191, 61192

RRID: AB_2793546

Isotype: IgG

Application(s): ICC, IF, WB

Reactivity: Human, Mouse

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 30 kDa

Background: LIN28 is an evolutionarily conserved RNA-binding protein that serves as a marker of undifferentiated human embryonic stem cells. LIN28 can facilitate the reprogramming of human somatic cells to induced pluripotent stem cells (iPSCs).

Immunogen: This LIN28A antibody was raised against a peptide derived from the C-terminal region of human LIN28A.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

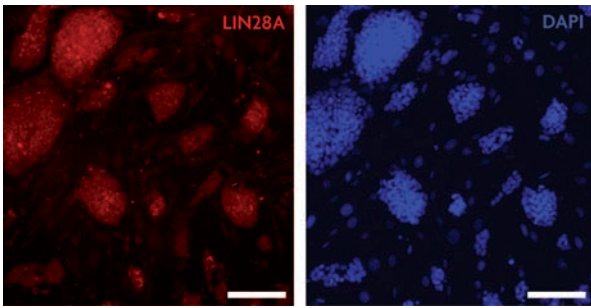
Applications Validated by Active Motif:

ICC/IF: 1:200 dilution

WB: 1:500 - 1:1,000 dilution

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



LIN28A antibody (pAb) tested by Immunocytochemistry

Mouse embryonic stem cells (mESCs) grown on mouse embryonic fibroblast feeder cells (MEFs) were fixed with 4% paraformaldehyde for 10 minutes at room temperature. Cells were then permeabilized and blocked by incubating with Blocking Solution containing 5% serum/0.1% Triton X-100 in D-PBS for 2 hours at room temperature. Cells were then incubated with LIN28A antibody at 1:200 dilution overnight at 4°C (left panel). Nuclei were stained with DAPI (right panel). Images show that LIN28A antibody specifically stains mESC colonies and does not stain MEFs. Absence of LIN28A staining in a subset of cells within the colonies suggests differentiation.

LIN28A antibody (pAb) tested by Western blot.

Cytosolic extract (40 µg) of mouse ES cells probed with LIN28A antibody at a dilution of 1:500.

