

L1ORF1p Polyclonal Antibody

(Catalog # A-2800)

Background

L1ORF1p, also known as recombinant mouse putative transposase element L1Md-A101/L1Md-A102/L1Md-A2, is a nucleic acid-binding protein. It is considered critical for retrotransposition of LINE-1 elements in the genome. L1ORF1p encodes two proteins, ORF1p and ORF2p. It may function as a nucleic acid chaperone binding its own transcript, thus preferentially mobilizing the transcript from which they are encoded.

Description

Rabbit polyclonal antibody to L1ORF1p.

Formulation

1*TBS (pH7.4), 0.5%BSA, 40%Glycerol. Preservative: 0.05% Sodium Azide.

Specificity

Mouse

Isotype

IgG

Purification

Affinity purified

Storage

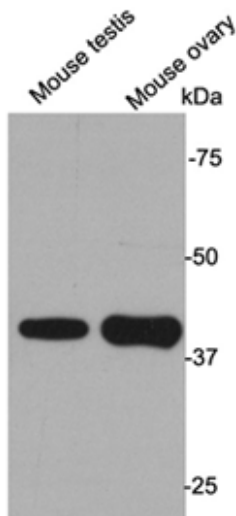
Store at 4°C after thawing. Aliquot and store at -20°C or -80°C. Avoid multiple freeze/thaw cycles.

Alternative Names

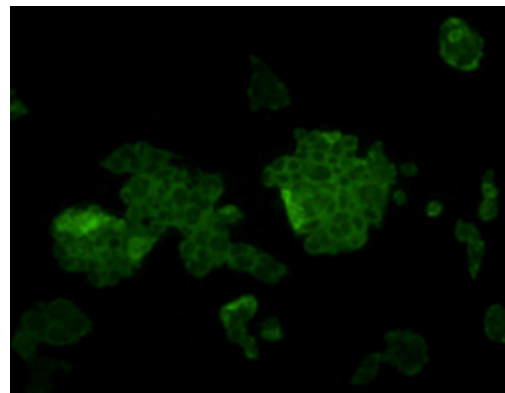
L1RE1, LORF1, Recombinant Mouse Putative transposase element L1Md-A101/L1Md-A102/L1Md-A2

Application

WB: (1:1,000), ICC: (1:50-1:100)



Western blot analysis on tissue lysates using L1ORF1p polyclonal antibody.



ICC staining L1ORF1p in F9 cells (green). Cells were fixed in paraformaldehyde, permeabilised with 0.25% Triton X100/PBS