

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

Uniprot

P11260

Purification

Affinity purified

Storage

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

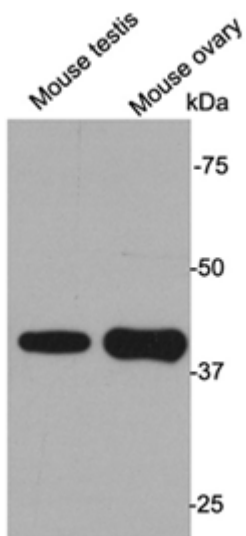
Alternative Names

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

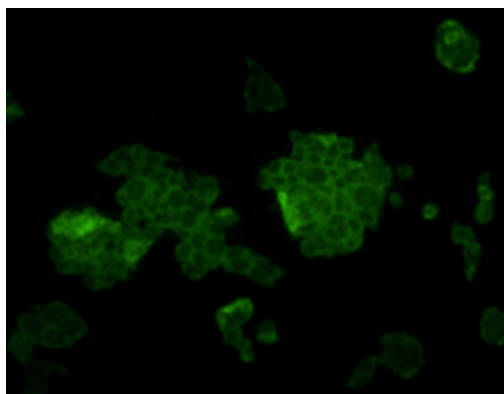
Application

WB, ICC; ; Recommended Dilution: WB: (1:1,000), ICC: (1:50-1:100)

** The BSA contained in this antibody is of negligible amounts, having undergone rigorous purification processes that include stringent heat and acid treatments. This comprehensive purification ensures the complete absence of viable microorganisms. As all BSA-containing products are strictly intended for research purposes and not for diagnostic or therapeutic use, they are exempt from certification by an authoritative body and are not subject to certification authority oversight.*



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.