
MuERVL-Gag Polyclonal Antibody

(Catalog # A-2801)

Background

When the zygotic genome is first transcribed, a large number of retrotransposons are expressed, including the endogenous retroviruses (ERVs), LINE-1 elements, and the non-autonomous SINE elements. MuERV-L/MERVL retrovirus-like particles are transiently de-repressed and produce 3% of the transcribed mRNAs at the two-cell (2C) stage. Right after the 2C stage, the expression of MERVL-retroelement is silenced. It has been found that this regulated pattern of MERVL expression overlapped with more than one hundred 2C-specific genes that have co-opted regulatory elements from these foreign retroviruses to begin their transcription.

Description

Rabbit polyclonal antibody to MuERVL-Gag.

Formulation

1x PBS, pH 7.3; 50% glycerol; 0.03% ProClin 300

Specificity

Mouse

Isotype

IgG

Uniprot

V9H130

Purification

Affinity purified

Storage

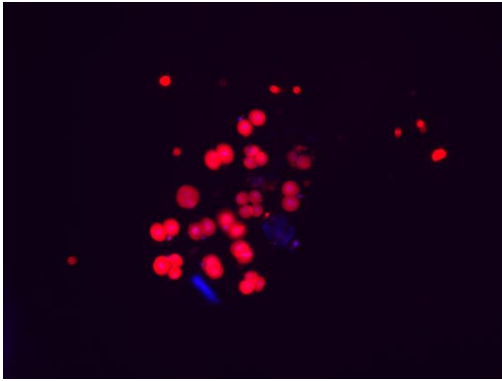
Store at 4°C after thawing. Aliquot store at -20°C or -80°C. Avoid repeated freeze / thaw cycles.

Alternative Names

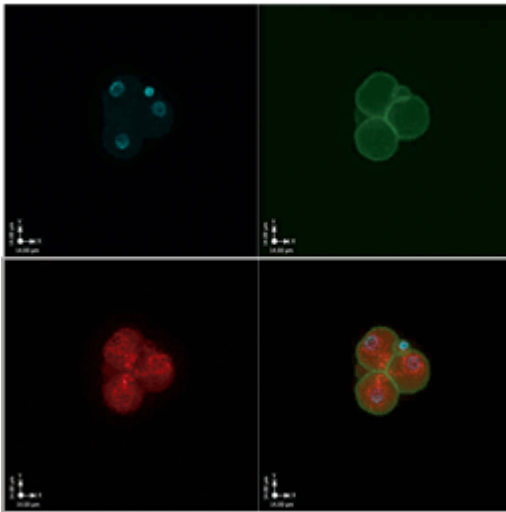
Murine Endogenous Retrovirus-Gag

Application

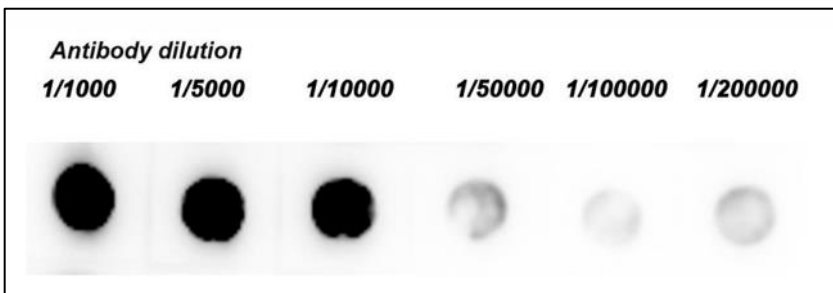
DB, IF, ICC; Recommended dilutions: DB (1:500-1:1000), IF (1:100 - 1:500), ICC (1:100-1:200)



Epifluorescence IF image using MuERVL-Gag Polyclonal Antibody in mouse preimplantation embryos from the 1 cell - 4 cell stages.



ICC staining MuERVL-Gag in mouse 3-cell stage embryo. The cells were incubated overnight at 4°C with MuERVL-Gag Polyclonal Antibody at a 1/100 dilution (shown in red) and phalloidin (shown in green). Nuclear DNA was labeled with DAPI (shown in blue).



Dot blot testing at various concentrations against 100 ng of MuERVL-Gag antigen peptide.