

---

# LMO1 Polyclonal Antibody

(Catalog # A71281)

---

## Background

This locus encodes a transcriptional regulator that contains two cysteine-rich LIM domains but lacks a DNA-binding domain. LIM domains may play a role in protein interactions; thus the encoded protein may regulate transcription by competitively binding to specific DNA-binding transcription factors. Alterations at this locus have been associated with acute lymphoblastic T-cell leukemia. Chromosomal rearrangements have been observed between this locus and at least two loci, the delta subunit of the T-cell antigen receptor gene and the LIM domain binding 1 gene. Alternatively spliced transcript variants have been described.

## Description

LMO1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

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

## Specificity

Human

## Isotype

IgG

## Uniprot ID

P25800

## Purification

Affinity Purification

## Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-156 of human LMO1 (NP\_002306.1).

## Storage

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

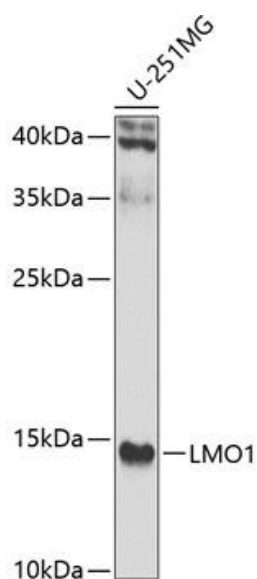
## Alternative Names

LMO1; RBTN1; RHOM1; TTG1; rhombotin-1

## Application

WB; Recommended dilution: WB: 1:500-1:2000

---



Western blot analysis of extracts of U-251MG cells, using LMO1 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: 90s.