

---

# ARL11 Polyclonal Antibody

(Catalog # A71907)

---

## Background

This gene encodes a tumor suppressor related to the ADP-ribosylation factor (ARF) family of proteins. The encoded protein may play a role in apoptosis in a caspase-dependent manner. Polymorphisms in this gene have been associated with some familial cancers.

## Description

ARL11 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

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

## Specificity

Rat

## Isotype

IgG

## Uniprot ID

Q969Q4

## Purification

Affinity Purification

## Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-196 of human ARL11 (NP\_612459.1).

## Storage

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

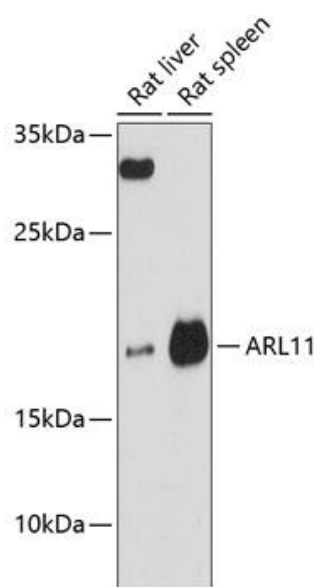
## Alternative Names

ARL11; ARLTS1; ADP ribosylation factor like GTPase 11

## Application

WB, IHC; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:100

---



Western blot analysis of extracts of various cell lines, using ARL11 Polyclonal Antibody at 1:3000 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.