

## AAR2 Polyclonal Antibody

(Catalog # A57922)

### Background

spliceosomal tri-snRNP complex assembly

### Description

AAR2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q9Y312

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Protein AAR2 homolog protein (2-250AA)

### Storage

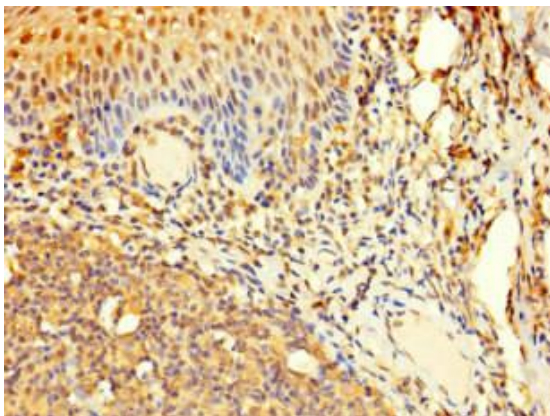
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Protein AAR2 homolog, AAR2, C20orf4, CGI-23, PRO0225, AAR2 splicing factor homolog

### Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human tonsil tissue using AAR2 Polyclonal Antibody at dilution of 1:100