

## APRIL Monoclonal Antibody [APRILY2]

(Catalog No. A-0617)

### Background

Cytokine that binds to TNFRSF13B/TACI and to TNFRSF17/BCMA. Plays a role in the regulation of tumor cell growth. May be involved in monocyte/macrophage-mediated immunological processes.

### Description

Mouse monoclonal antibody to APRIL, Clone APRILY2

### Formulation

Liquid. In PBS containing 0.02% sodium azide.

### Immunogen

Recombinant human APRIL (aa 93-233)

### Specificity

Human

### Isotype

IgG1

### Uniprot ID

O75888

### Purification

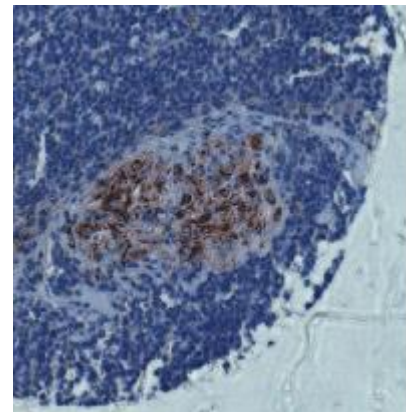
Serum

### Storage

Store at 4°C for short-term. Aliquot and store at -20°C for long term. Avoid multiple freeze/thaw cycles.

### Alternative Names

Interleukin-33, IL-1F11, NF-HEV



▲ Biopsy of tumor lesion from a patient with lymphoma (DLBCL) stained with APRIL Monoclonal Antibody [APRILY2]. The biopsy was stained by standard immunohistochemistry.

### Ordering Information

#### Products

APRIL Monoclonal Antibody [APRILY2]

#### Size

50 µg

100 µg

#### Cat. No.

A-0617-050

A-0617-100

*This product is for research purposes only. Not intended for use in diagnostic procedures.*