## APRIL (human) monoclonal antibody (Aprily-2)

This antibody is covered by our Worry-Free Guarantee.

Citations: 3

View Online »

**Ordering Information** 

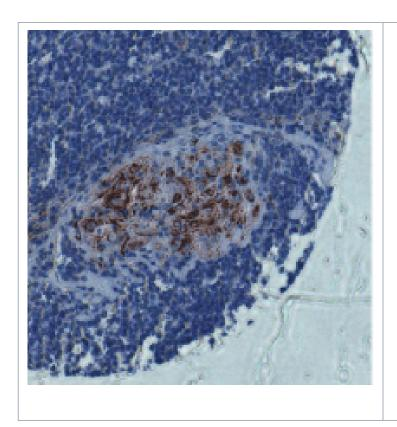
Order Online »

ALX-804-844-C100

100µg

Manuals, SDS & CofA

**View Online »** 



## **Handling & Storage**

**Use/Stability** Stable for at least 1 year after receipt when stored as recommended.

**Handling** Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name A-Proliferation-inducing ligand, TNFSF 13, CD256

**Application** IHC (PS)

**Application Notes** Immunohistochemistry (paraffin sections, 1:500).

Optimal conditions must be determined individually for

each application.

Clone Aprily-2

**Formulation** Liquid. In PBS containing 0.02% sodium azide.

**Host** Mouse

**Immunogen** Recombinant human APRIL (aa 93-233).

lgG1

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections

(1:500))Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be

determined individually for each application.

Source Purified from concentrated hybridoma tissue culture

supernatant.

Species Reactivity Human

O75888

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

