APRIL (human) monoclonal antibody (Aprily-5) (biotin conjugate)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

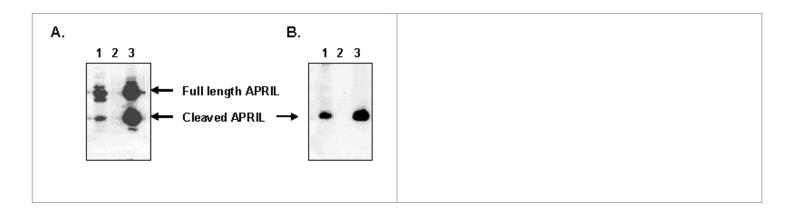
Order Online »

ALX-804-801B-C100

Biotin - 100µg

Manuals, SDS & CofA

View Online »



Handling & Storage

Short Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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

Application ELISA, IP, WB

Application Notes Excellent for Western Blot.

Clone Aprily-5

Crossreactivity Does not cross-react with mouse APRIL.

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Mouse

Immunogen Recombinant human APRIL (aa 93-233).

Isotype IgG1

Recommendation Dilutions/Conditions ELISA (detection)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

Specificity Recognizes full-length and cleaved APRIL.

UniProt ID 075888

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com