

APRIL Monoclonal Antibody [APRILY5]

(Catalog No. A-0587)

Background

The protein encoded by the gene APRIL (a proliferation-inducing ligand), also known as TNFSF13, is a member of the TNF ligand family. The TNF (tumor necrosis factors) family refers to a group of cytokines that can cause cell death (apoptosis). This protein is involved in B cell development. Experiments done *in vitro* suggested that this protein may induce apoptosis via its interaction with other TNF receptor family proteins such as TNFRSF6/FAS and TNFRSF14/HVEM. Alternative splicing causes numerous transcript variants. Brain glioblastoma multiforme and glioblastoma multiforme are two diseases associated with this protein.

Description

Mouse monoclonal antibody to APRIL, Clone APRILY5

Formulation

Liquid. In PBS containing 0.02% sodium azide.

Immunogen

Recombinant human APRIL (aa 93-233).

Specificity

Human

Isotype

IgG1

Purification

Protein A-affinity purified

Storage

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

Application

ELISA, IP, WB

Ordering Information

Products

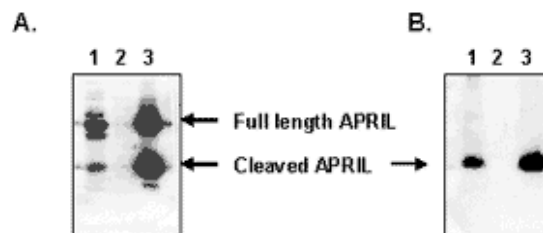
APRIL Monoclonal Antibody [APRILY5]

Size

100 µg

Cat. No.

A-0587-100



▲ **Fig. 1.** Detection of endogenous human APRIL with APRIL Monoclonal Antibody [APRILY5]. Method: Cell lysate (A) or cell supernatant (B) were run on an SDS-PAGE gel and analyzed with 2µg/ml Aprily-5 by Western blot. Dendritic cells (1), NIH3T3 (2) and NIH3T3 expressing human APRIL (3).