
BAFF Monoclonal Antibody [BUFFY1], FITC Conjugated

(Catalog No. A-0531)

Background

The protein encoded by this gene is a cytokine that belongs to the tumor necrosis factor (TNF) ligand family. This cytokine is a ligand for receptors TNFRSF13B/TACI, TNFRSF17/BCMA, and TNFRSF13C/BAFFR. This cytokine is expressed in B cell lineage cells, and acts as a potent B cell activator. It has been also shown to play an important role in the proliferation and differentiation of B cells. Alternatively spliced transcript variants encoding distinct isoforms have been identified. Diseases associated with TNFSF13B include Sialadenitis and Hypersensitivity Reaction Type II Disease.

Description

Rat monoclonal antibody to BAFF, Clone BUFFY1 (FITC conjugated)

Formulation

Liquid. In PBS containing 0.02% sodium azide.

Immunogen

Recombinant human soluble BAFF (aa 83-285)

Specificity

Human

Isotype

IgG2a

Uniprot ID

Q9Y275

Storage

Store at 4°C. Avoid multiple freeze/thaw cycles.

Alternative Names

B cell-activating factor, BLyS, TALL-1, THANK, zTNF4, TNFSF 13B/20, CD257

Application

Flow Cytometry, ICC, WB

Ordering Information

Products

BAFF Monoclonal Antibody [BUFFY1], FITC Conjugated

Size

100 µl

Cat. No.

A-0531-100

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