

**Purified Mouse Anti-human NTAL Antibody**  
**\*NAP-03, monoclonal, Cross Adsorbed\***

Catalog number: V1031895

Unit size: 0.1 mg

**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

**Antibody Properties**

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	NTAL
Clone	NAP-03

**Biological Properties**

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse serum and then conjugated with under optimal conditions
Application	IP, WB, IHC(P)

**Applications**

Linker for activation of T-cells family member 2 (alternatively called Non-T-cell activation linker, Williams-Beuren syndrome chromosomal region 15 protein, Williams-Beuren syndrome chromosomal region 5 protein, Membrane-associated adapter molecule or Linker for activation of B-cells) is a transmembrane protein with a molecular weight of 27 kDa, expressed in the membrane raft, extracellular exosome and mast cell granule of cells. Sequencing of Williams-Beuren syndrome chromosomal region 5 protein has demonstrated it contains 2 types of conserved structural units: cytoplasmic and extracellular domain. Williams-Beuren syndrome chromosomal region 5 protein recognizes SH2 domain and S-palmitoyl cysteine. It takes part in organismal processes, for instance, mast cell degranulation, intracellular signal transduction and adaptive immune response. Williams-Beuren syndrome chromosomal region 5 protein is the subject of intensive examination in part because of the fact that it is involved with the Fc-ε receptor signaling pathway and B cell receptor signaling pathway.