

**Purified Rabbit Anti-mouse/rat NTAL  
Antibody \*PAb (487), monoclonal, Cross  
Adsorbed\***Catalog number: V1031915  
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 | Mouse, rat |
| Class              | Primary    |
| Clonality          | Monoclonal |
| Host               | Rabbit     |
| Immunogen          | NTAL       |
| Clone              | PAb (487)  |

**Biological Properties**

|             |                                                                                                                                    |
|-------------|------------------------------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography, cross-adsorbed against human serum and then conjugated with under optimal conditions |
| Application | WB, ICC                                                                                                                            |

**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.