

**Ascites Mouse Anti-rat PLSCR1 Antibody**  
**\*13A6 [TEC-23], monoclonal\***

Catalog number: V1032030

Unit size: 0.1 ml

**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	Undiluted ascitic fluid, 15 mM sodium azide, 0.2% (w/v) BSA

**Antibody Properties**

Species Reactivity	Rat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	PLSCR1
Clone	13A6 [TEC-23]
Conjugate	Ascites

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with Ascites under optimal conditions
Application	IP, WB, FC, ICC

**Applications**

Phospholipid scramblase 1 (also known as Erythrocyte phospholipid scramblase, MmTRA1b or Ca(2+)-dependent phospholipid scramblase 1) is a transmembrane protein with a molecular weight of 35 kDa, located in the cytosol, nucleolus and collagen-containing extracellular matrix of cells. In Homo sapiens, phospholipid scramblase 1 is the subject of extensive research due to the fact that it acts in the regulation of Fc receptor mediated stimulatory signaling pathway. Phospholipid scramblase 1 plays an important role in mast cell activation. Phospholipid scramblase 1 aids in organismal processes, for instance, response to interferon- $\beta$ , defense response to virus and apoptotic process. It binds with epidermal growth factor receptor, calcium ion and S-palmitoyl cysteine. Sequencing of phospholipid scramblase 1 has shown it contains 2 types of conserved structural units: cytoplasmic and extracellular domain. Phospholipid scramblase 1 has been associated with critical functions such as phospholipid scramblase activity.