

**Purified Mouse Anti-human LRRC32
Antibody *GARP5, monoclonal***

Catalog number: V1031690

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	LRRC32
Clone	GARP5

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB

Applications

Transforming growth factor β activator LRRC32 is a 72 kda transmembrane protein that can be located in the integral component of plasma membrane, nucleoplasm and extracellular space of cells. It is alternatively called Leucine-rich repeat-containing protein 32, Garpin and Glycoprotein A repetitions predominant. In humans, LRRC32 is the subject of intensive study stemming from the fact that it acts in the transforming growth factor β receptor signaling pathway. Sequencing of LRRC32 has shown evidence of it contains 4 types of conserved structural units: extracellular, cytoplasmic, LRRNT and LRRC32 domain. LRRC32 is an integral part of transforming growth factor β activation and transforming growth factor beta3 activation. It downregulates activated T cell proliferation and cytokine secretion, on the other hand, it also enhances gene expression. LRRC32 binds to transforming growth factor β and Growth factor. LRRC32 takes part in organismal processes, for example, secondary palate development.