

**Purified Mouse Anti-human GRAP2
Antibody *UW40, monoclonal***

Catalog number: V103765

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	GRAP2
Clone	UW40

Biological Properties

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

Applications

GRB2-related adapter protein 2 is a 38 kDa protein that can be expressed in the nucleus, nucleoplasm and plasma membrane of cells. It is alternatively called Hematopoietic cell-associated adapter protein GrpL, Adapter protein GRID and Grf40 adapter protein. Sequencing of GRB2-related adapter protein 2 has supported it contains 3 types of conserved structural units: SH2, SH3 1 and SH3 2 domain. It is the subject of extensive examination stemming from the fact that it is a component of the Fc-ε receptor signaling pathway and T cell receptor signaling pathway. It takes part in organismal processes, for instance, T cell costimulation, cell-cell signaling and Ras protein signal transduction.