

**Purified Mouse Anti-human ARAP1  
Antibody \*ARAP1-2, monoclonal\***

Catalog number: V103125

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	ARAP1
Clone	ARAP1-2

**Biological Properties**

---

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

**Applications**

---

Arf-GAP with p-GAP domain, ANK repeat and PH domain-containing protein 1 is a protein with a molecular weight of 188 kDa, located in the plasma membrane, cytosol and Golgi apparatus of cells. In Homo sapiens, Arf-GAP with p-GAP domain, ANK repeat and PH domain-containing protein 1 is the subject of extensive examination stemming from the fact that it is a component of the positive regulation of epidermal growth factor receptor signaling pathway. Arf-GAP with p-GAP domain, ANK repeat and PH domain-containing protein 1 has been associated with critical functions such as GTPase activator activity. It reacts with metal ion, phosphatidylinositol-3,4,5-trisphosphate and type 1 angiotensin receptor. Sequencing of Arf-GAP with p-GAP domain, ANK repeat and PH domain-containing protein 1 has shown it contains 8 conserved structural units: Arf-GAP, Ras-associating, p-GAP, SAM, PH 1, PH 2, PH 3, and PH 4 domain. Arf-GAP with p-GAP domain, ANK repeat and PH domain-containing protein 1 inhibits stress fiber assembly, and it also upregulates GTPase activity, filopodium assembly and receptor recycling. It is an integral part of small GTPase mediated signal transduction and cellular component movement. Arf-GAP with p-GAP domain, ANK repeat and PH domain-containing protein 1 is an integral part of organismal processes, in particular, actin filament reorganization involved in cell cycle and signal transduction.