

**Purified Mouse Anti-human R-RAS2  
Antibody \*EM-50, monoclonal\***

Catalog number: V1032085

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	R-RAS2
Clone	EM-50

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	FC (QC TESTED), WB, ELISA, ICC

**Applications**

---

Ras-related protein R-Ras2 (sometimes called Ras-like protein TC21 or Teratocarcinoma oncogene) is a protein with a molecular weight of 23 kDa, found in the endoplasmic reticulum and extracellular exosome of cells. In humans, ras-related protein R-Ras2 reacts with GDP, GTP and S-farnesyl cysteine. it is an integral part of neuron death. Ras-related protein R-Ras2 plays an important role in organismal processes, in particular, osteoblast differentiation and ras protein signal transduction. Ras-related protein R-Ras2 has been associated with essential functions like GTPase activity. It acts to positively regulate cell migration. Mutations and abnormalities in as-related protein R-Ras2r have been associated with a number of diseases, in particular, ovarian cancer (OC) and noonan syndrome 12 (NS12).