

**Purified Mouse Anti-human NR2F6  
Antibody \*EM-51, monoclonal\***

Catalog number: V1031875

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	NR2F6
Clone	EM-51

**Biological Properties**

---

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

**Applications**

---

Nuclear receptor subfamily 2 group F member 6 is a 43 kDa protein that can be located in the nuclear chromatin, nucleoplasm and nucleus of cells. Sequencing of Nuclear receptor subfamily 2 group F member 6 has shown it contains a primary structural unit, the NR LBD domain. Nuclear receptor subfamily 2 group F member 6 recognizes zinc ion and sequence-specific DNA. It takes part in processes such as transcription by RNA polymerase II. Nuclear receptor subfamily 2 group F member 6 has been thought to be involved with critical functions such as DNA-binding transcription factor, nuclear receptor and steroid hormone receptor activity.