

**Purified Mouse Anti-human/cow Ki-67
Antibody *Ki-67, monoclonal***Catalog number: V1031580
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, cow
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Ki-67
Clone	Ki-67

Biological Properties

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

Applications

Proliferation marker protein Ki-67 is a protein with a molecular weight of 359 kDa, found in the membrane, chromosome and nucleus of cells. In humans, proliferation marker protein Ki-67 is an integral part of organismal processes, namely, cell cycle and cell population proliferation. It binds with DNA, RNA and ATP. Sequencing of proliferation marker protein Ki-67 has supported it contains 2 types of conserved structural units: PP1-binding and FHA domain. It plays an important role in chromatin organization, chromosome segregation and mitotic nuclear division.