

Purified Mouse Anti-human/non-human primates p21Waf1 Antibody *WA-1, monoclonal*Catalog number: V1031940
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, non-human primates
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	p21Waf1
Clone	WA-1

Biological Properties

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

Applications

Cyclin-dependent kinase inhibitor 1 (sometimes referred to as MDA-6, Melanoma differentiation-associated protein 6, p21 or CDK-interacting protein 1) is a protein with a molecular weight of 18 kDa, found in the PCNA-p21 complex, nucleoplasm and cytosol of cells. In humans, CDK-interacting protein 1 has been found to be involved in cyclin-dependent protein serine/threonine kinase activity, cell cycle G1/S phase transition and transcription by RNA polymerase II. It is the subject of extensive research stemming from the fact that it is a member of the intrinsic apoptotic signaling pathway, cytokine-mediated signaling pathway and intrinsic apoptotic signaling pathway in response to DNA damage by p53 class mediator. It negatively regulates cardiac muscle tissue regeneration, phosphorylation and cell population proliferation. On the other hand, it also enhances programmed cell death, B cell proliferation and fibroblast proliferation. CDK-interacting protein 1 binds to protein-containing complex, protein kinase and metal ion. It plays an important role in organismal processes, namely, stress-induced premature senescence, mitotic G2 DNA damage checkpoint and DNA damage response, signal transduction by p53 class mediator resulting in transcription of p21 class mediator.