

Purified Mouse Anti-human DAXX Antibody
DAXX-03, monoclonalCatalog number: V103555
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	DAXX
Clone	DAXX-03

Biological Properties

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

Applications

Death domain-associated protein 6 (sometimes referred to as ETS1-associated protein 1, EAP1, Fas death domain-associated protein, Daxx or hDaxx) is a protein with a molecular weight of 95 kDa, found in the nucleoplasm, nuclear body and PML body of cells. In humans, death domain-associated protein 6 has been thought to be involved with essential functions like protein homodimerization and protein kinase activator activity. Death domain-associated protein 6 is the subject of comprehensive application due to the fact that it is involved with the extrinsic apoptotic signaling pathway via death domain receptors and androgen receptor signaling pathway. It plays an important role in organismal processes, for example, cellular response to sodium arsenite, viral process and activation of JUN kinase activity, and furthermore, is thought to be essential to DNA-templated transcription, protein ubiquitination and apoptotic process. It reacts with protein kinase, p53 and protein N-terminus. It is a suppressor of DNA-templated transcription while also is a promoter of neuron death, protein phosphorylation and protein kinase activity.