

**Purified Mouse Anti-human/mouse Ikaros
Antibody *4.00E+09, monoclonal***

Catalog number: V1031550

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, mouse
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Ikaros
Clone	4.00E+09

Biological Properties

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

Applications

DNA-binding protein Ikaros is a 58 kDa protein that can be expressed in the cytoplasm, nucleus and protein-containing complex of cells. It is alternatively called Ikaros family zinc finger protein 1 and Lymphoid transcription factor LyF-1. In humans, DNA-binding protein Ikaros is an integral part of organismal processes, for example, lymphocyte differentiation, mesoderm development and chromatin organization, and furthermore, has been associated with key functions like DNA-binding transcription factor activity. DNA-binding protein Ikaros is a promoter of transcription by RNA polymerase II but also is an inhibitor of DNA-templated transcription. It reacts with transcription regulatory region sequence-specific DNA, metal ion and RNA polymerase II cis-regulatory region sequence-specific DNA. Mutations and abnormalities in DNA-binding protein Ikaros have been associated with a number of diseases, in particular, common variable immune deficiency 13 (CVID13).