

Purified Armenian hamster Anti-human/mouse Helios Antibody *22F6, monoclonal*Catalog number: V103800
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	Armenian hamster
Immunogen	Helios
Clone	22F6

Biological Properties

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

Applications

Zinc finger protein helios is a 60 kDa protein that can be expressed in the nucleoplasm of cells. In humans, zinc finger protein helios has been thought to be involved with critical functions such as DNA-binding transcription factor activity, and additionally, enhances transcription by RNA polymerase II. Zinc finger protein helios binds with transcription regulatory region sequence-specific DNA, metal ion and identical protein.