

**PE-Cy5.7 Armenian hamster Anti-human/mouse Helios Antibody \*22F6, monoclonal\***Catalog number: V103810  
Unit size: 100 tests**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
Conjugate	PE-Cy5.7

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with PE-Cy5.7 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.