

iFluor™ 405 Anti-human CD69 Antibody
FN50Catalog number: 10690020, 10690021
Unit size: 100 tests, 500 tests**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD69 (VEA, AIM, CLEC2C)
Clone	FN50
Conjugate	iFluor™ 405

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 405 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 405
Excitation Wavelength	403 nm
Emission Wavelength	427 nm

Applications

The FN50 monoclonal antibody binds to human CD69, a 32 kD transmembrane protein frequently found on the surface of thymocytes and granulocytes. CD69 is associated with a variety of biologically interesting macromolecules/ligands. CD69 is a very popular antibody target, with over 21000 publications in the last decade. CD69 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell

types, particularly in the study of immunology and costimulatory molecules. This antibody was purified through affinity chromatography and conjugated to iFluor™ 405 (ex/em = 403/427 nm). It is compatible with the 405 nm laser and 448/45 nm bandpass filter (for example, as in the BD FACSVerser™).