

FITC Anti-human CD71 Antibody *OKT-9*Catalog number: 107131I0, 107131I1
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, κ
Immunogen	CD71 (T9, Transferrin receptor, TFRC)
Clone	OKT-9
Conjugate	FITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	FITC
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

The OKT-9 monoclonal antibody binds with human CD71, a 95 kD glycoprotein commonly expressed on the surface of reticulocytes, proliferating cells, reticulocytes and erythroid precursors. In many organisms, CD71 acts to positively regulate B cell proliferation, is a positive regulator of bone resorption and is an enhancer of T cell proliferation. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands like Transferrin. CD71 is a fairly uncommon antibody target, with a little more than 4300 publications in the last decade. Even still, CD71 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to FITC (ex/em = 491/516 nm). It is

compatible with the 488 nm laser and 530/30 nm bandpass filter (for example, as in the BD FACSAria™ II).