

iFluor™ 810 Anti-human CD71 Antibody
OKT-9Catalog number: 10713000, 10713001
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	iFluor™ 810

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 810 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 810
Excitation Wavelength	811 nm
Emission Wavelength	822 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 iFluor™ 810 (ex/em = 811/822 nm).