

iFluor™ 810 Anti-human CD161 Antibody
HP-3G10Catalog number: 11610000, 11610001
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
Immunogen	CD161 (NKR-P1A)
Clone	HP-3G10
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

HP-3G10 is an anti-human monoclonal antibody that recognizes the CD161 antigen. CD161 (alternatively called NKR-P1, KLRB1, NKR-P1aKLRB1a or CD161aCD161b) is a 30 kD single-pass type II membrane protein that is expressed on the surface of cells such as T cells and NK cells. CD161 is a member of essential cellular pathways, for example, the cell surface receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands such as . CD161 is a fairly uncommon antibody target, with a little more than 2000 publications in the last decade. Even still, CD161 is vital to immunology research, typically serving as a phenotypic marker for

differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 810 (ex/em = 811/822 nm).