

iFluor™ 488 Anti-human CD203c Antibody
NP4D6Catalog number: 12030050, 12030051
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	CD203c (ENPP3)
Clone	NP4D6
Conjugate	iFluor™ 488

Biological Properties

Appearance	Orange-red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

The NP4D6 monoclonal antibody binds with human CD203c, a 150 kD transmembrane protein commonly located on the surface of mast cells and basophils. In certain organisms, CD203c is an inhibitor of mast cell proliferation, is an inhibitor of mast cell activation involved in immune response and represses inflammatory response. From a research standpoint, it is of biological interest due to its association with critical

macromolecules/ligands such as cAMP. CD203c is a relatively rare antibody target, with fewer than 500 publications in the last decade. Even still, CD203c is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology. This antibody was purified through affinity chromatography and conjugated to iFluor™ 488 (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 527/32 nm bandpass filter (for example, as in the BD FACSVerser™).