

**TRITC Anti-human CD203c Antibody**  
**\*NP4D6\***Catalog number: 120301H0, 120301H1  
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	TRITC

**Biological Properties**

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

**Spectral Properties**

Conjugate	TRITC
Excitation Wavelength	544 nm
Emission Wavelength	570 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 TRITC (ex/em = 544/570 nm). It is compatible with the 532 nm laser and 575/36 nm bandpass filter (for example, as in the Thermo Fisher Attune NxT).