

XFD555 Anti-human/ rabbit/ cat/ non-human primates/ ferret CD271 Antibody
NGFR5, XFD555 Same Structure to Alexa Fluor™ 555

Catalog number: 12710160, 12710161
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, rabbit, cat, non-human primates, ferret
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD271 (p75NTR, TNFRSF16, NGFR, Gp80-LNGFR)
Clone	NGFR5
Conjugate	AF555

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF555 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	AF555
Excitation Wavelength	553 nm
Emission Wavelength	568 nm

Applications

NGFR5 is an anti-human/ rabbit/ cat/ non-human primates/ ferret monoclonal antibody that is specific for the CD271 antigen. CD271 (sometimes referred to as TNFRSF16 or Gp80-LNGFR) is a 75 kD transmembrane protein that is found on the surface of cells like dendritic cells. CD271 is a component of important cellular pathways, namely, the neurotrophin TRK receptor signaling pathway, negative regulation of fibroblast growth factor receptor signaling pathway and positive regulation of apoptotic signaling pathway. Also, it has been associated with critical biological processes like glucose homeostasis, especially cellular glucose homeostasis. In certain organisms, CD271 plays a role in the upregulation of pri-miRNA transcription by RNA polymerase II, positively regulates apoptotic process and is a negative regulator of hair follicle development. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands like NT-3 and NT-4. CD271 is a fairly uncommon antibody target, with a little more than 1700 publications in the last decade. Even still, CD271 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of stem cells, immunology and neuroscience. This antibody was purified through affinity chromatography and conjugated to XFD555 (ex/em = 553/568 nm). XFD555 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 555 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 561 nm laser and 577/35 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).