

**iFluor™ 700 Anti-human/ rabbit/ cat/ non-human primates/ ferret CD271 Antibody  
\*NGFR5\***Catalog number: 127100J0, 127100J1  
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	iFluor™ 700

**Biological Properties**

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	iFluor™ 700
Excitation Wavelength	690 nm
Emission Wavelength	713 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 iFluor™ 700 (ex/em = 690/713 nm).