

**APC/iFluor™ 700 Anti-human CD10  
Antibody \*HI10a\***Catalog number: 101001F0, 101001F1, 101001F2  
Unit size: 25 tests, 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	CD10 (CALLA, MME, Neprilysin)
Clone	HI10a
Conjugate	APC/iFluor™ 700

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

---

Conjugate	APC/iFluor™ 700
Excitation Wavelength	685 nm
Emission Wavelength	710 nm

**Applications**

---

HI10a is an anti-human monoclonal antibody that targets the CD10 antigen. CD10 (also known as EPN, MME, CALLA or Neprilysin) is a 100 kD single-pass type II membrane protein that is expressed on the surface of cells such as NK cells and endothelial cells. CD10 has been closely linked to essential biological processes like amyloid-beta clearance, particularly amyloid-beta clearance by cellular catabolic process. Also, in some organisms, it is involved in the positive regulation of long-term synaptic potentiation and is a promoter of neurogenesis. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands like . CD10 is a fairly uncommon antibody

target, with a little more than 9600 publications in the last decade. Even still, CD10 has been widely used in 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 APC/iFluor™ 700 (ex/em = 685/710 nm).