

**APC Anti-human CD43 Antibody \*HI161\***Catalog Number: 104311C0, 104311C1, 104311C2  
Unit Size: 25 tests, 100 tests, 500 tests**Product Details**

---

Storage Conditions	N/A
Expiration Date	12 months upon receiving

**Chemical Properties**

---

Appearance	undefined
Molecular Weight	N/A
Soluble In	N/A

**Spectral Properties**

---

Excitation Wavelength	651 nm
Emission Wavelength	660 nm

**Applications**

---

The HI161 monoclonal antibody reacts with human CD43, a 95 - 135 kD transmembrane protein commonly located on the surface of plasma cells, thymocytes, neutrophils, myelomas and T cells. In some organisms, CD43 plays a role in the downregulation of cell adhesion, promotes tumor necrosis factor biosynthetic process and acts to negatively regulate T cell proliferation. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as EZR. CD43 is a fairly uncommon antibody target, with a little more than 5000 publications in the last decade. Even still, CD43 has been widely used in immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to APC (ex/em = 651/660 nm). It is compatible with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).