

## iFluor® 710 Anti-human CD43 Antibody \*HI161\*

Catalog Number: 104310K0, 104310K1

Unit Size: 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	717 nm
Emission Wavelength	739 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 iFluor® 710 (ex/em = 717/739 nm).