

# iFluor® 594 Anti-human CD43 Antibody

## \*HI161\*

Catalog Number: 104310C0, 104310C1

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	588 nm
Emission Wavelength	604 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® 594 (ex/em = 588/604 nm). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).