

# iFluor® 610 Anti-human CD43 Antibody

## \*HI161\*

Catalog Number: 104310D0, 104310D1

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 | 610 nm |
| Emission Wavelength   | 628 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® 610 (ex/em = 610/628 nm). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).