

PE Anti-human CD43 Antibody *HI161*

Catalog Number: 104311L0, 104311L1, 104311L2

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 | 566 nm |
| Emission Wavelength | 574 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 PE (ex/em = 566/574 nm). It is compatible with the 561 nm laser and 582/15 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).