

PE/XFD610 Anti-human CD11c Antibody *3.9*

Catalog Number: 101131O0,
101131O1, 101131O2
Unit Size: 25 tests, 100 tests, 500 tests

Product Details

| | |
|--------------------|---|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze. |
| Expiration Date | 12 months upon receiving |
| Concentration | Lot specific (please consult certificate of analysis for given lot) |
| Formulation | Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA |

Antibody Properties

| | |
|--------------------|--|
| Species Reactivity | Human |
| Class | Primary |
| Clonality | Monoclonal |
| Host | Mouse |
| Isotype | Mouse IgG1, κ |
| Immunogen | CD11c (Integrin alpha-X, CR4, p150, ITGAX) |
| Clone | 3.9 |
| Conjugate | PE/AF610 |

Biological Properties

| | |
|-----------------------|--|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions |
| Application | Flow Cytometry (FACS) |
| Recommended Dilutions | For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. |
| | <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i> |

Spectral Properties

| | |
|-----------|----------|
| Conjugate | PE/AF610 |
|-----------|----------|

Excitation Wavelength 565 nm

Emission Wavelength 627 nm

Applications

3.9 is an anti-human monoclonal antibody that forms an immune complex with the CD11c antigen. CD11c (sometimes called CR4 or ITGAX) is a 145 - 150 kD single-pass type I membrane protein that is expressed on the surface of cells such as macrophages, B cells and dendritic cells. In some organisms, CD11c promotes angiogenesis, positively regulates myelination and is a promoter of endothelial tube morphogenesis. Moreover, it is a component of essential cellular pathways, in particular, the cytokine-mediated signaling pathway and integrin-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands like ICAM-1 and 4 and fibrinogen. CD11c is a very popular antibody target, with over 26000 publications in the last decade. CD11c has been widely used in costimulatory molecules research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE/XFD610 (ex/em = 567/627 nm). XFD610 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 610 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 615/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCytte Quanteon).