

## PE/XFD610 Anti-human/ non-human primates CD184 Antibody \*12G5\*

Catalog Number: 118401N0,  
118401N1, 118401N2  
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, non-human primates |
| Class              | Primary                   |
| Clonality          | Monoclonal                |
| Host               | Mouse                     |
| Isotype            | Mouse IgG2a kappa         |
| Immunogen          | CD184 (CXCR4, Fusin)      |
| Clone              | 12G5                      |
| 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.<br><br><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

---

The 12G5 monoclonal antibody recognizes human/ non-human primates CD184, a 45 kD member of the Rhodopsin family commonly found on the surface of T cells, endothelial cells, hematopoietic progenitors, granulocytes and Tregs. CD184 is associated with a variety of biologically interesting macromolecules/ligands, for example, CXCL12, viral MIP-2 and SDF-1. CD184 is a relatively rare antibody target, with fewer than 1000 publications in the last decade. Even still, CD184 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of cell biology, neuroscience and neuroinflammation. 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 610/20 nm bandpass filter (for example, as in the BD FACSAria™ III).