

XFD594 Anti-human/ non-human primates CD184 Antibody

12G5

Catalog Number: 11840160, 11840161

Unit Size: 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	AF594

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF594 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
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	AF594
Excitation Wavelength	590 nm
Emission Wavelength	618 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 XFD594 (ex/em = 590/618 nm). XFD594 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 594 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).