

**iFluor™ 810 Anti-human/ non-human
primates CD184 Antibody *12G5***Catalog number: 11840000, 11840001
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	0.1 mg/mL
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	iFluor™ 810

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 810 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 810
Excitation Wavelength	811 nm
Emission Wavelength	822 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 iFluor™ 810 (ex/em = 811/822 nm).