

**Purified Anti-human/ non-human primates  
CD184 Antibody \*12G5\***Catalog number: 11840000  
Unit size: 100 ug**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.5 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

**Antibody Properties**

Species Reactivity	Human, non-human primates
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG2a kappa
Immunogen	CD184 (CXCR4, Fusin)
Clone	12G5

**Biological Properties**

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot

**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.