

**PE/iFluor™ 700 Anti-mouse/ dog CD80
Antibody *16-10A1***Catalog number: 108001X0, 108001X1, 108001X2
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	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	Mouse, dog
Class	Primary
Clonality	Monoclonal
Host	Armenian
Isotype	Armenian Hamster IgG2
Immunogen	CD80 (B7-1, B7, BB1)
Clone	16-10A1
Conjugate	PE/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 700
Excitation Wavelength	566 nm
Emission Wavelength	708 nm

Applications

The 16-10A1 monoclonal antibody reacts with mouse/ dog CD80, a 60 kD transmembrane glycoprotein commonly located on the surface of dendritic cells, macrophages, T cells, Tregs and B cells. CD80 is associated with a variety of biologically interesting macromolecules/ligands, in particular, CD152 (CTLA-4) and CD28. CD80 is a moderately popular antibody target, with over 19000 publications in the last decade. CD80 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology and neuroscience cell markers. This antibody was purified through affinity chromatography and conjugated to PE/iFluor™ 700 (ex/em = 566/708

nm). It is compatible with the 561 nm laser and 710/50 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).