

XFD647 Anti-mouse/ dog CD80 Antibody *16-10A1*

Catalog Number: 10800180, 10800181

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	Mouse, dog
Class	Primary
Clonality	Monoclonal
Host	Armenian
Isotype	Armenian Hamster IgG2
Immunogen	CD80 (B7-1, B7, BB1)
Clone	16-10A1
Conjugate	AF647

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF647 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 AF647

Excitation Wavelength 650 nm

Emission Wavelength 671 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 XFD647 (ex/em = 650/671 nm). XFD647 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 647 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 642 nm laser and 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).