

PE/XFD610 Anti-human CD274 Antibody *ACH0271*

Catalog Number: 12740100,
12740101, 12740102
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	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
Class	Primary
Clonality	Monoclonal
Host	Rabbit
Isotype	IgG
Immunogen	CD274 (PD-L1, B7-H1)
Clone	ACH0271
Conjugate	PE/AF610

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions
Application	Flow Cytometry (FACS)
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	PE/AF610
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 627 nm

Applications

ACH0271 is an anti-human monoclonal antibody that is specific for the CD274 antigen. CD274 (sometimes referred to as PD-L1 or B7-H1) is a 40 kD member of the immunoglobulin receptor superfamily that is expressed on the surface of cells such as NK cells, dendritic cells, B cells and macrophages. CD274 is associated with a variety of biologically interesting macromolecules/ligands, namely, PD1. CD274 is a fairly uncommon antibody target, with a little more than 4700 publications in the last decade. Even still, CD274 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of cancer biomarkers, costimulatory molecules and immunology. This antibody was purified through affinity chromatography and conjugated to PE/XFD610 (ex/em = 567/627 nm). XFD610 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 610 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).