



Rat anti-Mouse CD274 (PD-L1) conjugated to R-PE

Product Number	D5-126
Lot#	RPE056-19-001
Concentration	100ug/mL
Amount	100ug
Host	Rat
Isotype	IgG2b, κ
Clone	10F.9G2
Store at	+4°C

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at 2-8°C in the dark. Phycobiliproteins are sensitive to freeze-thaw cycles: after reconstitution, store at 2-8°C in the dark - do not freeze.

Handling

Avoid exposure to heat and light. Prior to use, reconstitute to 1 ml with distilled deionized water, vortex and allow it to sit on ice for 20 minutes.

Buffer

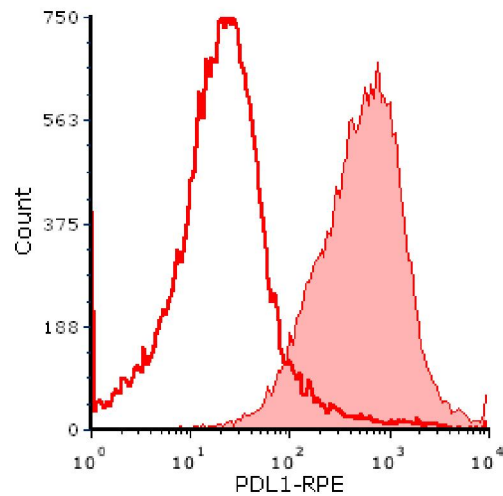
Upon reconstitution, the product is in 100mM sodium phosphate (pH 7.4), 50 mM sucrose, 100 mM sodium chloride, 0.1% BSA, and 2 mM sodium azide. The concentration of the conjugate is 100 μ g/ml.

Stability

The material is stable for up to one year if stored at 2-8°C in the dark. If further dilution of the conjugate is required, use diluted material within one week.

Note

For research use only, not for diagnostic or therapeutic use.



CD57BL/6 mouse splenocytes stained with R-PE conjugated anti-mouse CD274 Mab (filled histogram) or with matched isotype control (open histogram). Dilution = 5ug/test