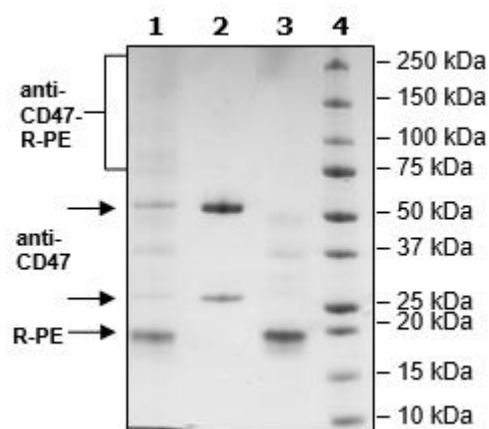


## Product Information

<b>Description:</b>	R-Phycoerythrin-labeled anti-CD47 recombinant human monoclonal recognizing the SIRP- $\alpha$ or SIRP- $\gamma$ binding region of human CD47.
<b>Label:</b>	R-Phycoerythrin (PE) is an oligomeric protein complex (270 kDa) from red algae that exhibits intensely bright red-orange fluorescence with high quantum yields. The complex consists of six heterodimers, $\alpha$ subunit (18 kDa) and $\beta$ -subunit (20 kDa), and an additional $\gamma$ -subunit (34 kDa). PE is covalently attached randomly through lysines on the target protein.
<b>Concentration:</b>	2.10 mg/ml
<b>Host Species:</b>	Human
<b>Isotype:</b>	IgG1
<b>Clonality:</b>	Monoclonal
<b>Cross Reactivity:</b>	This antibody has not been tested for cross reactivity with any other species.
<b>Formulated In:</b>	8 mM Na Phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl and 20% glycerol.
<b>Expression System:</b>	HEK293
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 12 months at -80°C. Avoid freeze/thaw cycles. Protect from light.
<b>Storage:</b>	-80°C
<b>MW:</b>	Heavy Chain: 51 kDa + PE, Light Chain: 26.5 kDa + PE
<b>Purity:</b>	≥90%
<b>Purification:</b>	Protein A affinity chromatography from HEK293 supernatants.
<b>Applications:</b>	Useful for labeling cells expressing CD47 for flow cytometry and immunofluorescence microscopy.

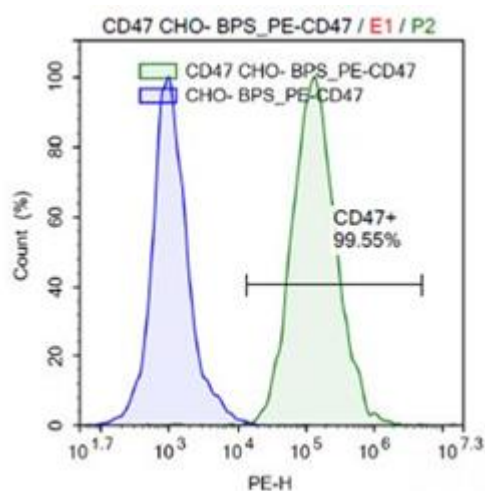
## Quality Control Data

### 4-20% SDS-Page Coomassie Staining



**Lane 1:** 8 µg anti-CD47-R-PE  
**Lane 2:** 4 µg anti-CD47  
**Lane 3:** 4 µg R-PE  
**Lane 4:** Protein Marker

### FACS Assay



Control CHO cells and CD47 expressing stable CHO cells were treated with human PE anti-CD47 antibody.