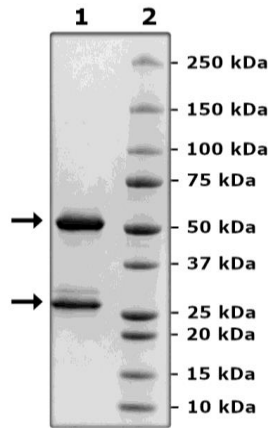


Product Information

| | |
|------------------------------|--|
| Description: | Recombinant monoclonal anti-CD25 (Basiliximab) which recognizes interleukin-2 receptor alpha subunit (IL-2RA). This affinity purified antibody has been tested for specific binding to purified human IL2RA protein (BPS Bioscience #101203) in a colorimetric ELISA binding assay. |
| Species | Human |
| Concentration: | 3.66 mg/ml |
| Expression System: | HEK293 |
| Purity: | ≥90% |
| Format: | Aqueous buffer solution. |
| Formulated In: | 8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol |
| MW: | Heavy Chain: 50 kDa; Light Chain: 23 kDa + glycans |
| Glycosylation: | This antibody runs at a higher MW by SDS-PAGE due to glycosylation. |
| Stability: | At least 12 months at -80°C. |
| Storage: | -80°C |
| Instructions for Use: | Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles. |
| Assay Conditions: | Functional validation: The antibody was validated by measuring Anti-IL-2RA binding to IL2RA antigen in ELISA assay. The IL2RA protein (BPS Bioscience #101203) was coated onto a 96-well plate overnight at 4°C (50 µl/well at a concentration of 2 µg/ml in PBS). The plate was washed 3 times with Immuno Buffer 1 (BPS Bioscience #79311) and blocked using 100 µl of Blocking Buffer 2 (BPS Bioscience #79728) for 1 hour at room temperature. After removing the blocking buffer, 50 µl/well of purified Anti-IL-2RA Antibody (BPS Bioscience #101593), serially diluted in Blocking Buffer 2, was added for 30 minutes at room temperature. After 3 more washes, the plate was incubated with Goat Anti-Human IgG Fc (HRP) (Abcam #ab97225), washed, and incubated with the Colorimetric HRP substrate. The reaction was stopped, and absorbance was read at 450 nm. The Blank value was subtracted from all values. |
| Applications: | Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling. |

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



IL2RA:Anti-IL-2RA (CD25) Activity

