

Recombinant PD-L1 / CD274 (19-239) protein

Catalog No: 81387, 81687

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.2 µg/µl

Source: Human

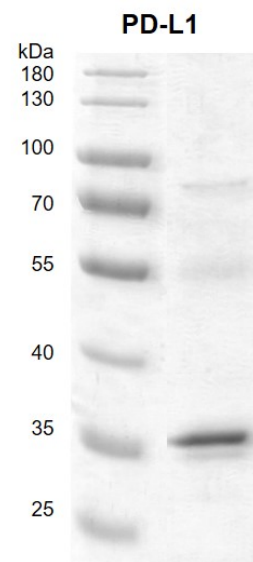
Buffer Contents: Recombinant PD-L1 / CD274 (19-239) protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

Background: PD-L1 is a ligand that binds with the receptor PD1, commonly found on T-cells, and acts to block T-cell activation. It plays a critical role in induction and maintenance of immune tolerance to self. As a ligand for the inhibitory receptor PDCD1/PD-1, PD-L1 modulates the activation threshold of T-cells and limits T-cell effector response. Through a yet unknown activating receptor, PD-L1 may costimulate T-cell subsets that predominantly produce interleukin-10 (IL10). The PDCD1-mediated inhibitory pathway is exploited by tumors to attenuate anti-tumor immunity and escape destruction by the immune system, thereby facilitating tumor survival. The interaction with PDCD1/PD-1 inhibits cytotoxic T lymphocytes (CTLs) effector function. The blockage of the PDCD1-mediated pathway results in the reversal of the exhausted T-cell phenotype and the normalization of the anti-tumor response, providing a rationale for cancer immunotherapy.

Protein Details: Recombinant PD-L1 / CD274 (19-239) protein that includes amino acids 19-239 of human PD-L1 protein (accession number NP_054862.1) was expressed in a baculovirus expression system, and contains an C-terminal FLAG tag. The molecular weight of the protein is 27.1 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of arrival.



Recombinant PD-L1 / CD274 (19-239) protein

10% SDS-PAGE Coomassie staining

MW: 27.1 kDa

Purity: >80%