



Product Data Sheet

Product Name:	Human PD-L1 inhibitor I	
Catalog Number:	AS-65581 (1 mg)	Lot Number: See label on vial
Sequence:	H-Phe-Asn-Trp-Asp-Tyr-Ser-Trp-Lys-Ser-Glu-Arg-Leu-Lys-Glu-Ala-Tyr-Asp-Leu-OH FNWDYSWKSERLKEAYDL	
Molecular Weight:	2350.7	
Peptide Purity:	Peak Area by HPLC ≥95%	
Storage:	This peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at –20°C or lower.	
Description:	<p>Programed cell death protein 1 (PD-1) is a cell surface receptor. PD-1 and its ligand PD-L1 are crucial immune checkpoints to downregulate T cell activation, tolerance and immunopathology. Several types of cancer cells overexpress PD-L1 in order to escape from the PD-1/PD-L1 immunosurveillance mechanism.</p> <p>Human PD-L1 inhibitor I is a peptide-based molecule. It binds to human PD-1 and inhibits PD-1/PD-L1 binding. This peptide, PDL-1 inhibitor I has anchor residues, WDY that influence binding of hPD-L1 to hPD-1. Developing inhibitors specifically blocking the PD-1/PD-L1 pathway has become a popular approach toward cancer treatment.</p>	
Related Products:	Human PD-L1 inhibitor II, Cat# AS-65582 Human PD-L1 inhibitor III, Cat# AS-65583 Human PD-L1 inhibitor IV, Cat# AS-65584 Human PD-L1 inhibitor V, Cat# AS-65585	
References:	<ol style="list-style-type: none">1. Bardhan K., et.al, <i>Front Immunol</i> 7(550), (2016)2. Li Q., et.al, <i>Oncotarget</i> 7(40), 64967-64976 (2016)3. D'Errico G., et.al, <i>Clin Transl Med</i> 6(1), 3-12 (2017)	

For Research Use Only