

Product Data Sheet

Product Name: Human PD-L1 inhibitor IV

Catalog Number: AS-65584 (1 mg) Lot Number: See label on vial

Sequence: H-Gly-Asn-Trp-Asp-Tyr-Asn-Ser-Gln-Arg-Ala-Gln-Leu-Tyr-Asn-Gln-OH

GNWDYNSQRAQLYNQ

Molecular Weight: 1857.0

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: 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.

Human PD-L1 inhibitor IV is a peptide-based molecule. It binds to human PD-1 and inhibits PD-1/PD-L1 binding. This peptide, hPDL-1 inhibitor IV has anchor residues (underlined), GNWDYNSQRAQLYNQ 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.

Related Products: Human PD-L1 inhibitor I, Cat# AS-65581

Human PD-L1 inhibitor II, Cat# AS-65582 Human PD-L1 inhibitor III, Cat# AS-65583 Human PD-L1 inhibitor V, Cat# AS-65585

References: 1. Bardhan K., et.al, Front Immunol **7(550)**, (2016)

2. Li Q., et.al, *Oncotarget* **7(40)**, 64967-64976 (2016) 3. D'Errico G., et.al, *Clin Transl Med* **6(1)**, 3-12 (2017)

For Research Use Only