



Product Data Sheet

Product Name:	SARS-CoV T-cell Peptide Antigens Set	
Catalog Number:	AS-65633 (0.25 mg net)	Lot Number: See label on vial
Size:	The set contains 5 peptides packed in individual vials of 0.25 mg each in net amount.	
% Peak Area by HPLC:	≥95%	
Appearance:	Lyophilized white powder	
Solvent:	Solvent recommendation available in certificate of analysis (CofA).	
Storage:	This peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower.	
Description:	<p>The SARS-CoV T-cell peptide antigens set contains 5 peptides corresponding to SARS-CoV spike protein sequence. The sequences have been tested for HLA binding with SARS-CoV protein and includes a positive and a negative control. CV1, CV2 and CV3 sequences are also seen within the spike glycoprotein of SARS-CoV2. The peptide sequences are offered in net peptide quantities in individual vials of 0.25 mg each. The peptide epitopes can be used in ELISPOT, cytokine and CTL assays. The sequences in the set are,</p> <p>CV1- VLNDILSRL CV2- NLNESLIDL CV3- FIAGLIAIV CV4- KLPDDFMGCV (+ve control) CV5- AYRPPNAPIL (-ve control)</p>	

References:

1. Lv Y, et al. Identification of a novel conserved HLA-A* 0201-restricted epitope from the spike protein of SARS-CoV. BMC immunology. 2009 Dec 1;10(1):61.
2. Ahmed SF et al. Preliminary identification of potential vaccine targets for the COVID-19 coronavirus (SARS-CoV-2) based on SARS-CoV immunological studies. Viruses. 2020 Mar;12(3):254.

For Research Use Only