



Product Data Sheet

Product Name:	SARS-CoV-2 T-cell Peptide Antigens Set	
Catalog Number:	AS-65640 (0.25 mg net)	Lot Number: See label on vial
Size:	The set contains 6 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>Identified from in silico studies, the set of 6 peptides offered are unique T-cell epitopes from the SARS-CoV-2 spike protein sequence. The peptide sequences are offered in net peptide quantities in individual vials of 0.25 mg each. The sequences in the set are,</p> <p>CV2-1 YLQPRTFLL CV2-2 GYVFASTK CV2-3 KLPDDFTGCV CV2-4 SIIAYTMSL CV2-5 NYNYLYRLFR CV2-6 GYLQPRTFLL</p>	

References:

1. 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.
2. Baruah V, & Bose S. Immunoinformatics-aided identification of T cell and B cell epitopes in the surface glycoprotein of 2019-nCoV. *Journal of medical virology*. 2020 May;92(5):495-500.
3. Grifoni A et al. A sequence homology and bioinformatic approach can predict candidate targets for immune responses to SARS-CoV-2. *Cell host & microbe*. 2020 Mar 16.
4. Coutard B et al. The spike glycoprotein of the new coronavirus 2019-nCoV contains a furin-like cleavage site absent in CoV of the same clade. *Antiviral research*. 2020 Apr 1;176:104742.
5. Kiyotani K et al. Bioinformatic prediction of potential T cell epitopes for SARS-Cov-2. *Journal of Human Genetics* (2020) 65:569–575.

For Research Use Only