

Technical Data Sheet

Product Name	<i>Cytomegalovirus (CMV) Control Peptide Pool</i>																																			
Size	1.25 mg																																			
Catalog #	AS-62339																																			
Description	<p>These 5 Cytomegalovirus (CMV) peptides, at 0.25 mg each (total of 1.25 mg/vial) have been used in the stimulation of IFNγ release from CD8+ T cells in individuals with defined HLA types, these epitopes are useful in applications such as ELISPOT, intracellular cytokine and CTL assays. All peptides are provided as net weight based on peptide content.</p> <table border="1"> <thead> <tr> <th>HLA Allele</th> <th>Virus</th> <th>Peptide Sequence</th> <th>MW</th> <th>Solvent</th> </tr> </thead> <tbody> <tr> <td>A0201</td> <td>HCMV pp65</td> <td>NLVPMVATV</td> <td>944.2</td> <td>DMSO</td> </tr> <tr> <td>B18</td> <td>HCMV</td> <td>SDEEEAIVAYTL</td> <td>1340.4</td> <td>Water</td> </tr> <tr> <td>B35</td> <td>CMV pp65</td> <td>IPSINVHHY</td> <td>1080.2</td> <td>Water</td> </tr> <tr> <td>B44</td> <td>HCMV</td> <td>EFFWDANDIY</td> <td>1320.4</td> <td>Water</td> </tr> <tr> <td>B0702</td> <td>HCMV</td> <td>TPRVTGGGAM</td> <td>947.1</td> <td>Water</td> </tr> <tr> <td>All of the above</td> <td>Pooled Peptides</td> <td>All of the above</td> <td>N/A</td> <td>DMSO</td> </tr> </tbody> </table>	HLA Allele	Virus	Peptide Sequence	MW	Solvent	A0201	HCMV pp65	NLVPMVATV	944.2	DMSO	B18	HCMV	SDEEEAIVAYTL	1340.4	Water	B35	CMV pp65	IPSINVHHY	1080.2	Water	B44	HCMV	EFFWDANDIY	1320.4	Water	B0702	HCMV	TPRVTGGGAM	947.1	Water	All of the above	Pooled Peptides	All of the above	N/A	DMSO
HLA Allele	Virus	Peptide Sequence	MW	Solvent																																
A0201	HCMV pp65	NLVPMVATV	944.2	DMSO																																
B18	HCMV	SDEEEAIVAYTL	1340.4	Water																																
B35	CMV pp65	IPSINVHHY	1080.2	Water																																
B44	HCMV	EFFWDANDIY	1320.4	Water																																
B0702	HCMV	TPRVTGGGAM	947.1	Water																																
All of the above	Pooled Peptides	All of the above	N/A	DMSO																																
Storage	-20°C																																			