



Product Information Sheet

Product Name: CEF Control Peptide Pool

The CEF Control Peptide Pool is a group of 32 peptides, 8-12 amino acids in length, with sequences derived from the human Cytomegalovirus, Epstein-Barr Virus and Influenza Virus. Used in the stimulation of IFN γ 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.

Size: These are pooled peptides
 cat# AS-61036-05, 0.5 mg each of the 32 peptides for a total of 16 mg/vial
 cat# AS-61036-025, 0.25 mg each, total 8 mg/vial
 cat# AS-61036-003, 0.03 mg each, total 1 mg/vial

Catalog Number: AS-61036-003

HPLC Purity: > 95%

Storage: -20°C

Peptides: (see table of 32 peptides)

HLA Allele	Virus	Peptide Purity (%)	Peptide Sequence	Solubility (1 mg/ml)	Size	Cat #s *
A1	Influenza A	>95	VSDGGPNLY	Water	1 mg	AS-28332
	Influenza A	>95	CTELKLSDY	Water	1 mg	AS-28333
A2	Influenza M	>95	GILGFVFTL	Water	1 mg	AS-28310
	Influenza A	>95	FMYSDHFHI	DMSO	1 mg	AS-28311
	EBV LMP2A	>95	CLGGLLTMV	Water	1 mg	AS-28319
	EBV BMLF1	>95	GLCTLVAML	DMSO	1 mg	AS-28320
A0201	HCMV pp65	>95	NLVPMVATV	Water	1 mg	AS-28328
AA68	Influenza NP	>95	KTGGPIYKR	Water	1 mg	n/a
A3	Influenza NP	>95	RVLSFIKGTK	Water	1 mg	n/a
	Influenza A	>95	ILRGSVAHK	Water	1 mg	AS-28334
	EBV	>95	RVRAYTYSK	Water	1 mg	n/a
	EBV	>95	RLRAEAQVK	Water	1 mg	n/a
A3, A11, A60B1	Influenza M	>95	SIIPSGPLK	Water	1 mg	n/a
A11	EBV EBNA 4NP	>95	AVFDRKSDAK	Water	1 mg	n/a



HLA Allele	Virus	Peptide Purity (%)	Peptide Sequence	Solubility	Size	Cat. # *
	EBV	>95	IVTDFSVIK	Water	1 mg	n/a
	EBV	>95	ATIGTAMYK	Water	1 mg	AS-28338
A24	EBV RTA	>95	DYCNVLNKEF	Water	1 mg	n/a
B7	Influenza NP	>95	LPFDKTTVM	Water	1 mg	n/a
	EBV	>95	RPPIFIRRL	Water	1 mg	n/a
B8	Influenza NP	>95	ELRSRYWAI	Water	1 mg	n/a
	EBV BZLF-1	>95	RAKFKQLL	Water	1 mg	n/a
	EBV EBNA 3A	>95	FLRGRAYGL	Water	1 mg	n/a
	EBV EBNA 3	>95	QAKWRLQTL	Water	1 mg	n/a
B18	HCMV	>95	SDEEEAIVAYTL	1% NH4OH	1 mg	n/a
B27	Influenza NP	>95	SRYWAIRTR	Water	1 mg	AS-28317
	Influenza M	>95	ASCMGLIY	Water	1 mg	n/a
	EBV EBNA 3C	>95	RRIYDLIEL	Water	1 mg	n/a
B35	EBV EBNA3A	>95	YPLHEQHGM	Water	1 mg	AS-28327
	CMV pp65	>95	IPSINVHHY	Water	1 mg	n/a
B44	EBV	>95	EENLLDFVRF	1% NH4OH	1 mg	n/a
	HCMV	>95	EFFWDANDIY	1% NH4OH	1 mg	n/a
B0702	HCMV	>95	TPRVTTGGAM	Water	1 mg	n/a
All of the above	Pooled Peptides	>95	All of the above	DMSO	16 mg	AS-61036-05
All of the above	Pooled Peptides	>95	All of the above	DMSO	8 mg	AS-61036-025
All of the above	Pooled Peptides	>95	All of the above	DMSO	1 mg	AS-61036-003

* n/a indicates the peptide is not sold separately, but may be available through bulk custom order

References: NIH AIDS Research & Reference Reagent Program
 Currier, JR. et al. J. Immunol. Methods 260, 157 (2002); Mwau, M. et al. AIDS Res. Hum. Retroviruses 18, 611 (2002)