

# PRODUCT INFORMATION



## Human CMV pp65 (495-503)

Item No. 44036

**CAS Registry No.:** 153045-21-7  
**Formal Name:** L-asparaginyll-L-leucyl-L-valyl-L-prolyl-L-methionyl-L-valyl-L-alanyl-L-threonyl-L-valine  
**Synonyms:** Asn-Leu-Val-Pro-Met-Val-Ala-Thr-Val, Human CMV 65 kDa Phosphoprotein (495-503), Human Cytomegalovirus pp65 (495-503)

**Peptide Sequence:** NLVPMVATV-OH

**MF:** C<sub>42</sub>H<sub>74</sub>N<sub>10</sub>O<sub>12</sub>S

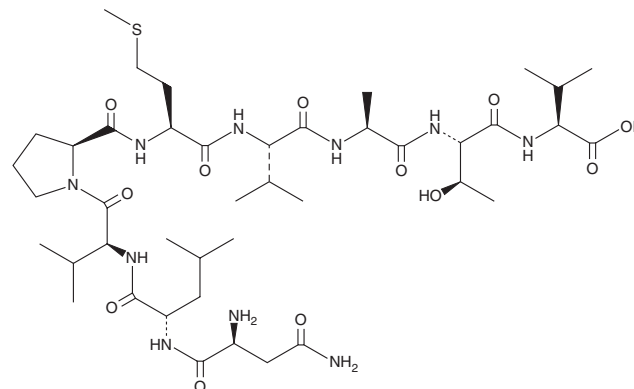
**FW:** 943.2

**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Human cytomegalovirus (CMV) pp65 (495-503) is supplied as a solid. A stock solution may be made by dissolving the human CMV pp65 (495-503) in the solvent of choice, which should be purged with an inert gas. Human CMV pp65 (495-503) is slightly soluble (0.1-1 mg/ml) in DMSO. It is also slightly soluble (0.1-1 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Human CMV pp65 (495-503) is an immunogenic T cell epitope derived from the CMV matrix protein 65 kDa phosphoprotein (pp65).<sup>1,2</sup> It binds to HLA-A\*0201 in a cell-free assay (IC<sub>50</sub> = 90 nM) and induces activation of patient-derived antigen-specific CD8<sup>+</sup> T cells.<sup>1</sup> Human CMV pp65 (495-503) has been used as an antigen to generate anti-pp65 antibodies with antiviral activity in a mouse model.<sup>2</sup>

### References

1. Gibadullin, R., Randall, C.J., Sidney, J., *et al.* Backbone modifications of HLA-A2-restricted antigens induce diverse binding and T cell activation outcomes. *J. Am. Chem. Soc.* **143(17)**, 6470-6481 (2021).
2. Chen, D., Hu, C., Hong, Y., *et al.* T cell receptor-like antibody specifically targets and eliminates cells infected with cytomegalovirus. *J. Transl. Med.* **23(1)**, 846 (2025).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/10/2025

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM