

PRODUCT INFORMATION



Enfuvirtide

Item No. 24097

CAS Registry No.: 159519-65-0

Formal Name: N-acetyl-L-tyrosyl-L-threonyl-L-seryl-L-leucyl-L-isoleucyl-L-histidyl-L-seryl-L-leucyl-L-isoleucyl-L- α -glutamyl-L- α -glutamyl-L-seryl-L-glutamyl-L-asparaginyl-L-glutamyl-L-glutamyl-L- α -glutamyl-L-lysyl-L-asparaginyl-L- α -glutamyl-L-glutamyl-L- α -glutamyl-L-leucyl-L-leucyl-L- α -glutamyl-L-leucyl-L- α -aspartyl-L-lysyl-L-tryptophyl-L-alanyl-L-seryl-L-leucyl-L-tryptophyl-L-asparaginyl-L-tryptophyl-L-phenylalaninamide

Ac—Tyr—Thr—Ser—Leu—Ile—His—Ser—Leu—Ile—Glu—
Glu—Ser—Gln—Asn—Gln—Gln—Glu—Lys—Asn—Glu—
Gln—Glu—Leu—Leu—Glu—Leu—Asp—Lys—Trp—Ala—
Ser—Leu—Trp—Asn—Trp—Phe—NH₂

MF: C₂₀₄H₃₀₁N₅₁O₆₄

FW: 4,491.9

Purity: $\geq 95\%$

UV/Vis.: λ_{\max} : 279 nm

Supplied as: A crystalline 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

Enfuvirtide is supplied as a crystalline solid. A stock solution may be made by dissolving the enfuvirtide in the solvent of choice, which should be purged with an inert gas. Enfuvirtide is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of enfuvirtide in these solvents is approximately 5 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of enfuvirtide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of enfuvirtide in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Enfuvirtide is a biomimetic peptide inhibitor of HIV-1 fusion with CD4⁺ cells.¹ It inhibits HIV-1 infectivity of HeLa cells stably expressing CD4 by the HXB2 strain (IC₅₀ = 692 pM) and by clinical isolates (IC₉₀₅ = 6.1-61 nM) in a single-cycle infectivity assay. It also inhibits genomic integration of HIV-1 into human intraepithelial vaginal cells and peripheral blood mononuclear cells (PBMCs; IC₅₀₅ = 51.2 and 13.58 μ M, respectively).² Enfuvirtide binds to a recombinant molecular mimic of HIV-1 glycoprotein gp41 that contains three N-terminal heptad and two C-terminal heptad repeat regions (K_d = 32 nM).³ It also binds to recombinant formyl peptide receptors (FPR) expressed in rat basophilic leukemia cells (IC₅₀ = 5 nM) and attracts and activates human peripheral blood phagocytes, but not T lymphocytes, *in vitro* when used at a concentration of 100 nM.⁴ Formulations containing enfuvirtide have been used in combination therapy for the treatment of HIV-1/AIDS.

References

1. Kariya, T., Wille, L.J., and Dage, R.C. *J. Cardiovasc. Pharmacol.* **4(3)**, 509-514 (1982).
2. Sebhat, I.K., Martin, W.J., Ye, Z., et al. *J. Med. Chem.* **45(21)**, 4589-4593 (2002).
3. Martin, W.J., McGowan, E., Cashen, D.E., et al. *Eur. J. Pharmacol.* **454(1)**, 71-79 (2002).
4. Cepoi, D., Phillips, T., Cismowski, M., et al. *Brain Res.* **1000(1-2)**, 64-71 (2004).

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, 11/03/2022

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