

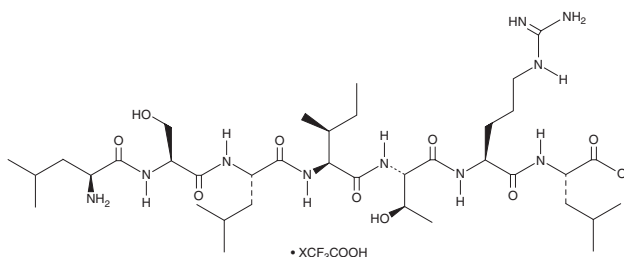
PRODUCT INFORMATION



S7 (trifluoroacetate salt)

Item No. 37467

Formal Name: L-leucyl-L-seryl-L-leucyl-L-threonyl-L-arginyl-L-leucine, trifluoroacetate salt
Synonym: Leu-Ser-Leu-Iso-Thr-Arg-Leu-OH
Peptide Sequence: LSLITRL-OH
MF: C₃₇H₇₀N₁₀O₁₀ • XCF₃COOH
FW: 815.0
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

S7 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the S7 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. S7 (trifluoroacetate salt) is soluble in 1% formic acid/water.

Description

S7 is a peptide antagonist of the IL-6 receptor.^{1,2} It inhibits IL-6 binding to the IL-6 receptor in a concentration-dependent manner.¹ S7 (50 μM) inhibits IL-6-induced increases in VEGF levels in C-33 A cervical cancer cells and RPMI-8226 B cell lymphocytes. It reduces tumor volume in an IL-6-overexpressing C-33 A cervical cancer mouse xenograft model when administered at a dose of 50 mg/kg every two days. When coupled to cysteine and the surface of lipid nanoparticles (LNPs) encapsulating doxorubicin, S7 increases LNP glioma targeting and increases survival in a U251 glioblastoma mouse xenograft model.²

References

1. Su, J.-L., Lai, K.-P., Chen, C.-A., *et al.* A novel peptide specifically binding to interleukin-6 receptor (gp80) inhibits angiogenesis and tumor growth. *Cancer Res.* **65(11)**, 4827-4835 (2005).
2. Shi, W., Cui, X., Shi, J., *et al.* Overcoming the blood-brain barrier for glioma-targeted therapy based on an interleukin-6 receptor-mediated micelle system. *RSC Adv.* **7**, 27162-27169 (2017).

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, 10/10/2023

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