

PRODUCT INFORMATION



AUNP-12

Item No. 28053

CAS Registry No.: 1353563-85-5

Formal Name: N²,N⁶-bis(L-seryl-L-asparaginyl-L-threonyl-L-seryl-L-α-glutamyl-L-seryl-L-phenylalanyl)-L-lysyl-L-phenylalanyl-L-arginyl-L-valyl-L-threonyl-L-glutamyl-L-leucyl-L-alanyl-L-prolyl-L-lysyl-L-alanyl-L-glutamyl-L-isoleucyl-L-lysyl-L-α-glutamine

Synonyms: AUR-012, SNTSEFK(SNTSESF) FRVTQLAPKAQIKE-NH₂

MF: C₁₄₂H₂₂₆N₄₀O₄₈

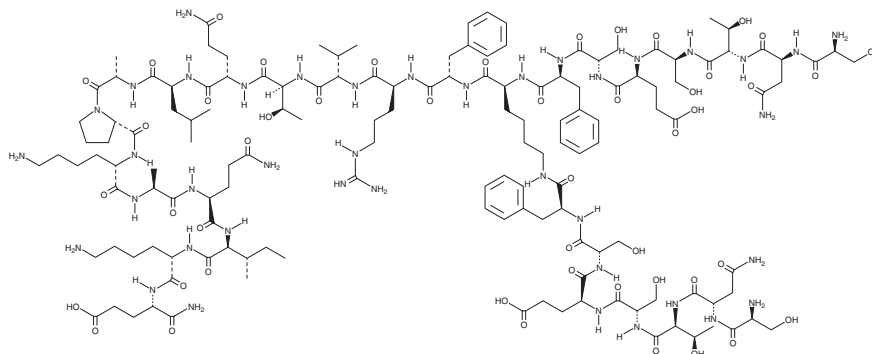
FW: 3,261.6

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

AUNP-12 is supplied as a solid. A stock solution may be made by dissolving the AUNP-12 in water. We do not recommend storing the aqueous solution for more than one day.

Description

AUNP-12 is an inhibitor of the interaction between programmed cell death 1 (PD-1) and its ligands, PD-L1 and PD-L2.¹ It inhibits binding of soluble PD-1 to HEK293 cells expressing human PD-L2 (EC₅₀ = 0.72 nM). AUNP-12 reverses PD-L1-induced inhibition of proliferation of PMA-stimulated, PD-1-expressing rat peripheral blood mononuclear cells (PBMCs; EC₅₀ = 0.41 nM). It reduces tumor growth and lung metastasis by 44 and greater than 60%, respectively, in a B16/F10 mouse melanoma model when administered at a dose of 5 mg/kg per day.

Reference

1. Sasikumar, P.G.N., Ramachandra, M., Vadlamani, S.K., *et al.* Immunosuppression modulating compounds. *Aurigen Discovery Technologies Limited*. **WO2011161699A2** (2011).

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, 12/14/2022

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