

PRODUCT INFORMATION



Transportan (trifluoroacetate salt)

Item No. 37497

Formal Name: glycyl-L-tryptophyl-L-threonyl-L-leucyl-L-asparaginyl-L-seryl-L-alanylglycyl-L-tyrosyl-L-leucyl-L-leucylglycyl-L-lysyl-L-isoleucyl-L-asparaginyl-L-leucyl-L-lysyl-L-alanyl-L-leucyl-L-alanyl-L-alanyl-L-leucyl-L-alanyl-L-lysyl-L-lysyl-L-isoleucyl-L-leucinamide, trifluoroacetate salt

Peptide Sequence: GWTLSAGYLLGKINLKALAALAKKIL-NH₂

MF: C₁₃₄H₂₂₇N₃₅O₃₂ • XCF₃COOH

FW: 2,840.5

Purity: ≥90%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

H-Gly-Trp-Thr-Leu-Asn-Ser-Ala-Gly-Tyr-Leu-Leu-Gly-Lys-Ile-Asn-Leu-Lys-Ala-Leu-Ala-Ala-Leu-NH₂ • XCF₃COOH

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

Laboratory Procedures

Transportan (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the transportan (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Transportan (trifluoroacetate salt) is soluble in a 10:90 solution of acetonitrile:water. We do not recommend storing the aqueous solution for more than one day.

Description

Transportan is a cell-penetrating peptide.¹ It is composed of 12 amino acids corresponding to amino acids 1-12 of the neuropeptide galanin (Item No. 33139) linked by a lysine residue to the mast cell degranulating peptide mastoparan (Item No. 20240). Transportan has been used in the surface functionalization of liposomes for drug delivery *in vitro* and *in vivo* and for the delivery of fluorescent cargo *in vitro*.^{2,3}

References

1. Shah, V.R., Shah, Y.D., and Athalye, M.N. Novel approaches in development of cell penetrating peptides. *J. Appl. Pharm. Res.* **9(1)**, 8-24 (2021).
2. Patel, S.G., Sayers, E.J., He, L., et al. Cell-penetrating peptide sequence and modification dependent uptake and subcellular distribution of green fluorescent protein in different cell lines. *Sci. Rep.* **9(1)**, 6298 (2019).
3. Li, Y.-X., Wang, N., Hasan, M.M., et al. Co-administration of transportan peptide enhances the cellular entry of liposomes in the bystander manner both *in vitro* and *in vivo*. *Mol. Pharm.* **19(11)**, 4123-4134 (2022).

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, 10/03/2023

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