

PRODUCT INFORMATION



740 Y-P (trifluoroacetate salt)

Item No. 22598

Formal Name: L-arginyl-L-glutamyl-L-isoleucyl-L-lysyl-L-isoleucyl-L-tryptophyl-L-phenylalanyl-L-glutamyl-L-asparagyl-L-arginyl-L-arginyl-L-methionyl-L-lysyl-L-tryptophyl-L-lysyl-L-lysyl-L-seryl-L- α -aspartylglycylglycyl-O-phosphono-L-tyrosyl-L-methionyl-L- α -aspartyl-L-methionyl-L-serine, trifluoroacetate salt

Arg—Gln—Ile—Lys—Ile—Trp—Phe—Gln—Asn—Arg—

Arg—Met—Lys—Trp—Lys—Lys—Ser—Asp—Gly—Gly—

pTyr—Met—Asp—Met—Ser

• XCF₃COOH

Peptide Sequence: RQIKIWFQNRRMKWKKSDGG-[pY]-MDMS

MF: C₁₄₁H₂₂₂N₄₃O₃₉PS₃ • XCF₃COOH

FW: 3,270.7

Purity: ≥85%

UV/Vis.: λ_{\max} : 210, 221, 281 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

740 Y-P (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 740 Y-P (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. 740 Y-P (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 20 mg/ml.

740 Y-P (trifluoroacetate salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 740 Y-P (trifluoroacetate salt) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 740 Y-P (trifluoroacetate salt) has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

740 Y-P is a synthetic phosphopeptide activator of PI3K.^{1,2} It contains amino acids from the Antennapedia protein, which acts as a cellular internalization sequence, followed by a sequence from PDGFR that is recognized by the SH2 domain of PI3K.¹ 740 Y-P (15 μ M) increases phosphorylation of PI3K, mammalian target of rapamycin (mTOR), and glycogen synthase kinase 3 β (GSK3 β) in HepG2 hepatocellular carcinoma cells.² It prevents glucose-oxygen deprivation-induced decreases in viability of primary rat hippocampal cells in an *in vitro* model of ischemia when used at a concentration of 1 μ M.³ 740 Y-P (1 μ M), in combination with the thrombopoietin receptor (TPOR) agonist butyzamide, hematopoietic stem cell (HSC) self-renewal agonist UM171, and polymer PCL-PVAc-PEG, has been used to expand isolated human umbilical cord blood HSCs *ex vivo* for up to 30 days.⁴

References

1. Derossi, D., Williams, E.J., Green, P.J., *et al.* Stimulation of mitogenesis by a cell-permeable PI 3-kinase binding peptide. *Biochem. Biophys. Res. Commun.* **251**(1), 148-152 (1998).
2. Qin, B., Zeng, Z., Xu, J., *et al.* Emodin inhibits invasion and migration of hepatocellular carcinoma cells via regulating autophagy-mediated degradation of snail and β -catenin. *BMC Cancer* **22**(1), 671 (2022).
3. Turovsky, E.A., Turovskaya, M.V., Gaidin, S.G., *et al.* Cytokine IL-10, activators of PI3-kinase, agonists of α -2 adrenoreceptor and antioxidants prevent ischemia-induced cell death in rat hippocampal cultures. *Arch. Biochem. Biophys.* **615**, 35-43 (2017).
4. Sakurai, M., Ishitsuka, K., Ito, R., *et al.* Chemically defined cytokine-free expansion of human haematopoietic stem cells. *Nature* **615**(7950), 127-133 (2023).

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, 11/15/2023

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