

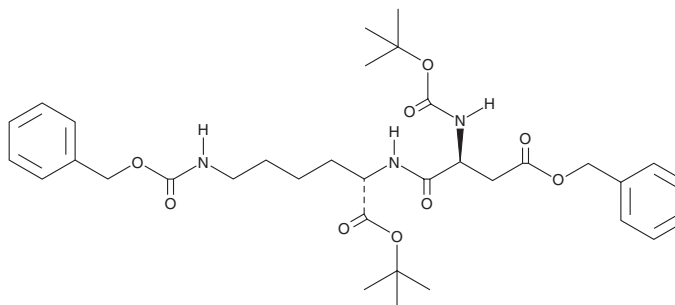
PRODUCT INFORMATION



Reversin 121

Item No. 18603

CAS Registry No.: 174630-04-7
Formal Name: 2-(1,1-dimethylethyl) 1-(phenylmethyl) ester N-[(1,1-dimethylethoxy)carbonyl]-L-aspartyl-N⁶-[(phenylmethoxy)carbonyl]-L-lysine
MF: C₃₄H₄₇N₃O₉
FW: 641.8
Purity: ≥98%
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

Reversin 121 is supplied as a crystalline solid. A stock solution may be made by dissolving the reversin 121 in the solvent of choice, which should be purged with an inert gas. Reversin 121 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of reversin 121 in these solvents is approximately 12, 20, and 25 mg/ml, respectively.

Reversin 121 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, reversin 121 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Reversin 121 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Reversin 121 is a hydrophobic peptide chemosensitizer that can reverse P-glycoprotein-mediated multidrug resistance. It binds to the P-glycoprotein multidrug transporter (MDR1) with a K_d value of 77 nM.^{1,2} Reversin 121 modulates P-glycoprotein ATPase activity in Sf9 insect cell membranes expressing human MDR1, plasma membrane vesicles from multidrug-resistant cells, and reconstituted proteoliposomes, as well as in a variety of MDR1-expressing intact tumor cells.¹

References

1. Sharom, F.J., Yu, X., Lu, P., *et al.* Interaction of the P-glycoprotein multidrug transporter (MDR1) with high affinity peptide chemosensitizers in isolated membranes, reconstituted systems, and intact cells. *Biochem. Pharmacol.* **58(4)**, 571-586 (1999).
2. Sharom, F.J., Liu, R., Romsicki, Y., *et al.* Insights into the structure and substrate interactions of the P-glycoprotein multidrug transporter from spectroscopic studies. *Biochim. Biophys. Acta.* **1461(2)**, 327-345 (1999).

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, 10/04/2022

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