

PRODUCT INFORMATION



SL 0101-1

Item No. 11704

CAS Registry No.: 77307-50-7
Formal Name: 3-[[3,4-di-O-acetyl-6-deoxy- α -L-mannopyranosyl]oxy]-5,7-dihydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one

MF: C₂₅H₂₄O₁₂

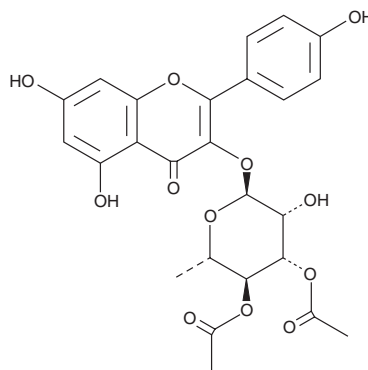
FW: 516.5

Purity: \geq 95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

SL 0101-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the SL 0101-1 in the solvent of choice, which should be purged with an inert gas. SL 0101-1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SL 0101-1 in these solvents is approximately 0.16 and 0.33 mg/ml, respectively.

Description

The p90 ribosomal S6 kinases (RSK) 1-4 are downstream members of the extracellular signal-regulated kinase (ERK)/MAPK cascade. SL 0101-1 is a kaempferol glycoside, isolated from the tropical plant *F. refracta*, that selectively inhibits RSK2 with an IC₅₀ value of 89 nM ($K_i = 1 \mu\text{M}$) without interfering with upstream activators of RSK, including ERK, MEK, EGFR, and PKC.¹ At 100 μM , SL 0101-1 inhibits the proliferation of MCF-7 breast cancer cells, arresting cells in the G₁ phase of the cell cycle.¹ SL 0101-1 has been used to characterize the intracellular signaling events associated with angiotensin II-induced tyrosine kinase activation with a concentration of 30 μM attenuating angiotensin II-induced cell proliferation.²

References

1. Smith, J.A., Poteet-Smith, C.E., Xu, Y., *et al.* Identification of the first specific inhibitor of p90 ribosomal S6 kinase (RSK) reveals an unexpected role for RSK in cancer cell proliferation. *Cancer Res.* **65**(3), 1027-1034 (2005).
2. Godeny, M.D. and Sayeski, P.P. ERK1/2 regulates ANG II-dependent cell proliferation via cytoplasmic activation of RSK2 and nuclear activation of elk1. *Am. J. Physiol. Cell Physiol.* **291**, C1308-C1317 (2006).

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, 10/13/2022

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