

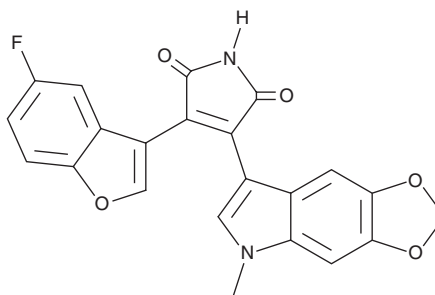
PRODUCT INFORMATION



9-ING-41

Item No. 36629

CAS Registry No.: 1034895-42-5
Formal Name: 3-(5-fluoro-3-benzofuranyl)-4-(5-methyl-5H-1,3-dioxolo[4,5-f]indol-7-yl)-1H-pyrrole-2,5-dione
Synonym: Elraglusib
MF: C₂₂H₁₃FN₂O₅
FW: 404.4
Purity: ≥98%
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

9-ING-41 is supplied as a solid. A stock solution may be made by dissolving the 9-ING-41 in the solvent of choice, which should be purged with an inert gas. 9-ING-41 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 9-ING-41 in these solvents is approximately 1 and 2 mg/ml, respectively.

Description

9-ING-41 is an inhibitor of glycogen synthase kinase 3 β (GSK3 β ; IC₅₀ = 0.71 μ M).¹ It inhibits the growth of MiaPaCa-2, BxPC-3, and HuP-T3 pancreatic cancer cells (IC₅₀s = 5, 1, and 0.6 μ M, respectively). 9-ING-41 (5 μ M) inhibits GSK3 β phosphorylation and induces apoptosis in BxPC-3 and HuP-T3 cells. *In vivo*, 9-ING-41 (40 mg/kg) reduces tumor growth in an SKOV3 mouse xenograft model.² It also reduces pleural thickening and improves lung function in a mouse model of *S. pneumoniae*-induced empyema.³

References

1. Gaisina, I.N., Gallier, F., Ougolkov, A.V., *et al.* From a natural product lead to the identification of potent and selective benzofuran-3-yl-(indol-3-yl)maleimides as glycogen synthase kinase 3 β inhibitors that suppress proliferation and survival of pancreatic cancer cells. *J. Med. Chem.* **52(7)**, 1853-1863 (2009).
2. Hilliard, T.S., Gaisina, I.N., Muehlbauer, A.G., *et al.* Glycogen synthase kinase 3 beta inhibitors induce apoptosis in ovarian cancer cells and inhibit *in vivo* tumor growth. *Anticancer Drugs* **22(10)**, 978-985 (2011).
3. Boren, J., Shryock, G., Fergis, A., *et al.* Inhibition of glycogen synthase kinase 3 β blocks mesomesenchymal transition and attenuates *Streptococcus pneumoniae*-mediated pleural injury in mice. *Am. J. Pathol.* **187(11)**, 2461-2472 (2017).

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, 10/18/2022

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