

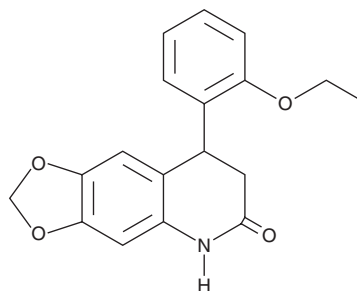
PRODUCT INFORMATION



FQI 1

Item No. 23104

CAS Registry No.: 599151-35-6
Formal Name: 8-(2-ethoxyphenyl)-7,8-dihydro-1,3-dioxolo[4,5-g]quinolin-6(5H)-one
Synonym: Factor Quinolinone Inhibitor 1
MF: C₁₈H₁₇NO₄
FW: 311.3
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 273, 302 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

FQI 1 is supplied as a crystalline solid. A stock solution may be made by dissolving the FQI 1 in the solvent of choice, which should be purged with an inert gas. FQI 1 is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

FQI 1 is an inhibitor of Late SV40 Factor (LSF) transactivation activity (IC₅₀ = 2.1 μM).¹ It inhibits growth of NIH3T3, HeLa S3, and A549 cancer cells (GI₅₀s = 3.8, 0.79, and 6.3 μM, respectively). FQI 1 decreases viability of QGY-7703 and Huh7 liver cancer cells, but not of non-cancerous Hc3716-hTERT cells and primary mouse hepatocytes, when used at a concentration of 4 μM.^{1,2} It reduces tumor growth induced by N-nitrosodiethylamine (DEN) in mice and in a QGY-7703 hepatocellular carcinoma (HCC) mouse xenograft model when administered at doses of 4 and 1 mg/kg, respectively.

References

1. Grant, T.J., Bishop, J.A., Christadore, L.M., *et al.* Antiproliferative small-molecule inhibitors of transcription factor LSF reveal oncogene addiction to LSF in hepatocellular carcinoma. *Proc. Natl. Acad. Sci. U.S.A.* **109(12)**, 4503-4508 (2012).
2. Rajasekaran, D., Siddig, A., Wiloughby, J.L., *et al.* Small molecule inhibitors of Late SV40 Factor (LSF) abrogate hepatocellular carcinoma (HCC): Evaluation using an endogenous HCC model. *Oncotarget* **6(28)**, 26266-26277 (2015).

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, 10/06/2022

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