

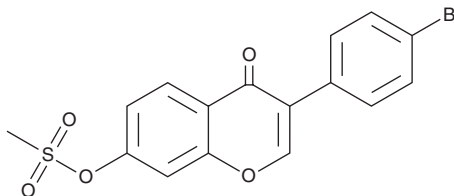
PRODUCT INFORMATION



KIN101

Item No. 37429

CAS Registry No.: 610753-87-2
Formal Name: 3-(4-bromophenyl)-7-[(methylsulfonyl)oxy]-4H-1-benzopyran-4-one
MF: $C_{16}H_{11}BrO_5S$
FW: 395.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 253 nm
Supplied as: A neat solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

KIN101 is supplied as a neat solid. A stock solution may be made by dissolving the KIN101 in the solvent of choice, which should be purged with an inert gas. KIN101 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of KIN101 in these solvents is approximately 10 and 25 mg/ml, respectively.

Description

KIN101 is an antiviral agent.¹ It inhibits the replication of hepatitis C virus 2a (HCV2a) in Huh7 cells ($EC_{50} = 0.2 \mu M$). KIN101 (10 μM) reduces HCV2a RNA levels in infected Huh7 cells and influenza A nucleoprotein levels in infected MRC-5 cells. It increases the expression of innate immune response genes, including *IRF3*, the gene encoding the transcription factor interferon regulatory factor 3 (IRF3), in Sendai virus-infected MRC-5 cells when used at a concentration of 10 μM .

Reference

1. Bedard, K.M., Wang, M.L., Proll, S.C., *et al.* Isoflavone agonists of IRF-3 dependent signaling have antiviral activity against RNA viruses. *J. Virol.* **86**(13), 7334-7344 (2012).

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, 02/22/2023

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