

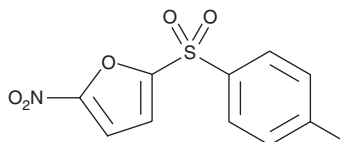
PRODUCT INFORMATION



NSC 697923

Item No. 15346

CAS Registry No.: 343351-67-7
Formal Name: 2-[(4-methylphenyl)sulfonyl]-5-nitro-furan
MF: C₁₁H₉NO₅S
FW: 267.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Protect from light



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

Laboratory Procedures

NSC 697923 is supplied as a solid. A stock solution may be made by dissolving the NSC 697923 in the solvent of choice, which should be purged with an inert gas. NSC 697923 is soluble in the organic solvent DMSO.

Description

NSC 697923 is a selective inhibitor of the E2 ubiquitin-conjugating enzyme Ubc13 when heterodimerized with Uev1A, blocking polyubiquitin chain formation at concentrations of 0.5 to 2 μ M *in vitro*.¹ It also inhibits NF- κ B activation by PMA and other agonists, presumably through its effects on Ubc13.¹ NSC 697923 blocks proliferation and induces apoptosis in diffuse large B cell lymphoma and neuroblastoma cells.^{1,2} It has antitumor efficacy *in vivo* in neuroblastoma orthotopic xenografts, when given intraperitoneally.²

References

1. Pulvino, M., Liang, Y., Oleskyn, D., *et al.* Inhibition of proliferation and survival of diffuse large B-cell lymphoma cells by a small-molecule inhibitor of the ubiquitin-conjugating enzyme Ubc13-Uev1A. *Blood* **120(8)**, 1668-1677 (2012).
2. Cheng, J., Fan, Y.-H., Xu, X., *et al.* A small-molecule inhibitor of UBE2N induces neuroblastoma cell death via activation of p53 and JNK pathways. *Cell Death Dis.* **5(e1079)**, (2014).

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, 12/15/2022

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