

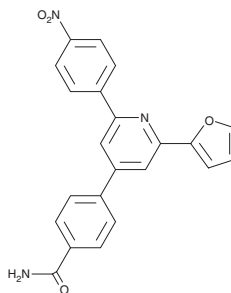
PRODUCT INFORMATION



KJ Pyr 9

Item No. 19116

CAS Registry No.: 581073-80-5
Formal Name: 4-[2-(2-furanyl)-6-(4-nitrophenyl)-4-pyridinyl]-benzamide
MF: C₂₂H₁₅N₃O₄
FW: 385.4
Purity: ≥98%
UV/Vis.: λ_{max}: 280, 345 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

KJ Pyr 9 is supplied as a crystalline solid. A stock solution may be made by dissolving the KJ Pyr 9 in the solvent of choice, which should be purged with an inert gas. KJ Pyr 9 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of KJ Pyr 9 in these solvents is approximately 20 and 25 mg/ml, respectively.

KJ Pyr 9 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, KJ Pyr 9 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. KJ Pyr 9 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

KJ Pyr 9 is a cell-permeable inhibitor of c-Myc ($K_d = 6.5$ nM) that interferes with Myc-Max complex formation in cells.^{1,2} It preferentially blocks the proliferation of c-Myc-overexpressing cells, reducing Myc-driven gene expression.¹ KJ Pyr 9 also potently blocks cell growth directed by overexpression of v-Myc, while less effectively blocking N-Myc- and L-Myc-dependent proliferation.² It is effective *in vivo*, reducing the growth of breast cancer MDA-MB-231 xenografts in mice.¹

References

1. Hart, J.A., Garner, A.L., Yu, J., *et al.* Inhibitor of MYC identified in a Kröhnke pyridine library. *Proc. Natl. Acad. Sci. USA* **111**(34), 12556-12561 (2014).
2. Raffener, P., Röck, R., Schrafl, A., *et al.* *In vivo* quantification and perturbation of Myc-Max interactions and the impact on oncogenic potential. *Oncotarget* **5**(19), 8869-8878 (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, 12/14/2022

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