

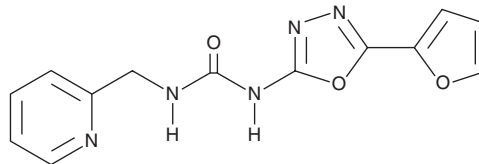
PRODUCT INFORMATION



NK 252

Item No. 19131

CAS Registry No.: 1414963-82-8
Formal Name: N-[5-(2-furanyl)-1,3,4-oxadiazol-2-yl]-N'-(2-pyridinylmethyl)-urea
MF: C₁₃H₁₁N₅O₃
FW: 285.3
Purity: ≥98%
UV/Vis.: λ_{max}: 275 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

NK 252 is supplied as a crystalline solid. A stock solution may be made by dissolving the NK 252 in the solvent of choice. NK 252 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of NK 252 in these solvents is approximately 14 and 20 mg/ml, respectively.

NK 252 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NK 252 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. NK 252 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NK 252 is a Nrf2 activator that interacts directly with the domain containing the Nrf2-binding site of Keap1.¹ At 1.36 μM, it is reported to activate the NQO1-antioxidant response element in a luciferase reporter gene assay two-fold above background.¹ NK 252 has been shown to downregulate the expression of fibrogenic genes, producing antifibrotic effects in a rat model of non-alcoholic steatohepatitis.¹

Reference

1. Shimozono, R., Asaoka, Y., Yoshizawa, Y., *et al.* Nrf2 activators attenuate the progression of nonalcoholic steatohepatitis-related fibrosis in a dietary rat model. *Mol. Pharmacol.* **84**(1), 62-70 (2013).

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, 11/29/2022

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