

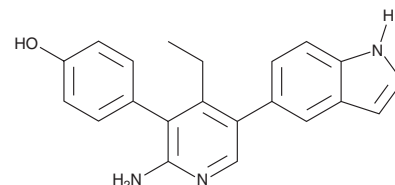
PRODUCT INFORMATION



GNE-6640

Item No. 33751

CAS Registry No.: 2009273-67-8
Formal Name: 4-[2-amino-4-ethyl-5-(1H-indazol-5-yl)-3-pyridinyl]-phenol
MF: $C_{20}H_{18}N_4O$
FW: 330.4
Purity: $\geq 98\%$
Supplied as: A 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

GNE-6640 is supplied as a solid. A stock solution may be made by dissolving the GNE-6640 in the solvent of choice, which should be purged with an inert gas. GNE-6640 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GNE-6640 in these solvents is approximately 1 mg/ml.

GNE-6640 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GNE-6640 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GNE-6640 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GNE-6640 is an inhibitor of ubiquitin-specific protease 7 (USP7; $IC_{50} = 0.75 \mu\text{M}$).¹ It is selective for USP7 over USP47 and USP5 (IC_{50} s = 20.3 and $>200 \mu\text{M}$, respectively). GNE-6640 enhances ubiquitination of the USP7 substrate MDM2 in HCT116 colon cancer cells ($IC_{50} = 0.23 \mu\text{M}$) and reduces cell viability in a panel of 108 cell lines (IC_{50} s = $<10 \mu\text{M}$ for all). GNE-6640 (10-70 μM) enhances apoptosis induced by doxorubicin (Item No. 15007) or cisplatin (Item No. 13119) in MCF-7 and U2OS cancer cells, respectively.

Reference

1. Kategaya, L., Di Lello, P., Rouge, L., et al. USP7 small-molecule inhibitors interfere with ubiquitin binding. *Nature* **550(7677)**, 534-538 (2017).

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, 11/18/2022

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