

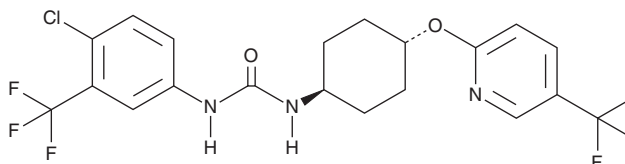
PRODUCT INFORMATION



UC2288

Item No. 37380

CAS Registry No.: 1394011-91-6
Formal Name: N-[4-chloro-3-(trifluoromethyl)phenyl]-N'-[trans-4-[[5-(trifluoromethyl)-2-pyridinyl]oxy]cyclohexyl]-urea
MF: C₂₀H₁₈ClF₆N₃O₂
FW: 481.8
Purity: ≥98%
UV/Vis.: λ_{max}: 253 nm
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

UC2288 is supplied as a solid. A stock solution may be made by dissolving the UC2288 in the solvent of choice, which should be purged with an inert gas. UC2288 soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of UC2288 in these solvents is approximately 25 mg/ml.

Description

UC2288 is an inhibitor of cyclin-dependent kinase inhibitor (p21^{WAF1}), a negative regulator of cell cycle progression.¹ It reduces the protein levels and mRNA expression of p21^{WAF1} independently of Raf and p53, respectively, when used at a concentration of 10 μM. UC2288 (8 μM) decreases the proliferation of, and phosphorylation of EGFR and extracellular-signal-related kinases 1/2 (ERK1/2) in, CNE-2 and 5-8F nasopharyngeal cell lines.² *In vivo*, UC2288, in combination with the telomerase inhibitor imetelstat, reduces tumor growth and induces apoptosis in an HCT116 colorectal cancer mouse xenograft model when administered at a dose of 15 mg/kg three times per week.³

References

1. Wettersten, H.I., Hwang, S.H., Li, C., *et al.* A novel p21 attenuator which is structurally related to sorafenib. *Cancer Biol. Ther.* **14**(3), 278-285 (2013).
2. Liang, R. and Zhu, X. UC2288 induces cell apoptosis of nasopharyngeal carcinoma cells via inhibiting EGFR/ERK pathway. *J. Cancer* **12**(4), 988-995 (2021).
3. Gupta, R., Dong, Y., Solomon, P.D., *et al.* Synergistic tumor suppression by combined inhibition of telomerase and CDKN1A. *Proc. Natl. Acad. Sci. USA* **111**(30), E3062-E3071 (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, 02/20/2023

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