

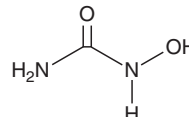
PRODUCT INFORMATION



Hydroxyurea

Item No. 23725

CAS Registry No.: 127-07-1
Formal Name: N-hydroxy-urea
Synonyms: NCI C04831, NSC 32065
MF: $\text{CH}_4\text{N}_2\text{O}_2$
FW: 76.1
Purity: $\geq 95\%$
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

Hydroxyurea is supplied as a crystalline solid. A stock solution may be made by dissolving the hydroxyurea in the solvent of choice, which should be purged with an inert gas. Hydroxyurea is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of hydroxyurea in these solvents is approximately 1 mg/ml. Hydroxyurea is also soluble in water. The solubility of hydroxyurea in water is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Hydroxyurea is an antineoplastic agent that inhibits DNA replication and cell proliferation and induces cell cycle arrest in the G_1/S phase.¹ It inhibits proliferation of NCI H460 human lung cancer cells *in vitro* ($\text{IC}_{50} = 0.56 \text{ mM}$).² Hydroxyurea reduces ribonucleoside diphosphate reductase (RNR) activity by 75% when used at a concentration of 5 mM *via* degradation of its active site tyrosyl radical.^{1,3} In a humanized mouse model of sickle cell disease (SCD), it reduces leukocyte adhesion and extravasation *via* NO production and a cGMP-dependent pathway.⁴ Hydroxyurea also inhibits HIV-1 viral replication in peripheral blood mononuclear cells (PBMCs) in a dose-dependent manner with an IC_{90} value of 0.4 mM.^{5,6} Formulations containing hydroxyurea have been used to treat cancer, SCD, and psoriasis.^{1,7,8}

References

1. Madaan, K., Kaushik, D., and Verma, T. Hydroxyurea: A key player in cancer chemotherapy. *Expert Rev. Anticancer Ther.* **12**(1), 19-29 (2012).
2. Yakisich, J.S., Azad, N., Venkatadri, R., et al. Digitoxin and its synthetic analog MonoD have potent antiproliferative effects on lung cancer cells and potentiate the effects of hydroxyurea and paclitaxel. *Oncol. Rep.* **35**(2), 878-886 (2016).
3. Krakoff, I.H., Brown, N.C., and Reichard, P. Inhibition of ribonucleoside diphosphate reductase by hydroxyurea. *Cancer Res.* **28**(8), 1559-1565 (1968).
4. Almeida, C.B., Scheiermann, C., Jang, J.E., et al. Hydroxyurea and a cGMP-amplifying agent have immediate benefits on acute vaso-occlusive events in sickle cell disease mice. *Blood* **120**(14), 2879-2888 (2012).
5. Gao, W.-Y., Johns, D.G., and Mitsuya, H. Anti-human immunodeficiency virus type 1 activity of hydroxyurea in combination with 2',3'-dideoxynucleosides. *Mol. Pharmacol.* **46**(4), 767-772 (1994).
6. Lori, F. and Lisiewicz, J. Rationale for the use of hydroxyurea as an anti-human immunodeficiency virus drug. *Clin. Infect. Dis.* **30**(Suppl 2), S193-S197 (2000).
7. Platt, O.S. Hydroxyurea for the treatment of sickle cell anemia. *N. Engl. J. Med.* **358**(13), 1362-1369 (2008).

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, 02/12/2024

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