

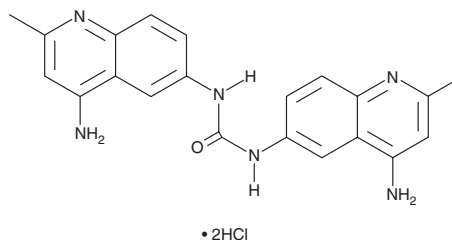
PRODUCT INFORMATION



Surfen (hydrochloride)

Item No. 26670

CAS Registry No.: 5424-37-3
Synonym: Aminoquinicarbamide
Formal Name: N,N'-bis(4-amino-2-methyl-6-quinolinyl)-urea, dihydrochloride
MF: C₂₁H₂₀N₆O • 2HCl
FW: 445.4
Purity: ≥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

Surfen (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the surfen (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Surfen (hydrochloride) is slightly soluble (0.1-1 mg/ml) in a 1:1 solution of acetonitrile:water and sparingly soluble (1-10 mg/ml) in methanol and DMSO.

Description

Surfen is an antagonist of cell surface glycosaminoglycans.^{1,2} It binds to heparin, dermatan sulfate, heparan sulfate, and chondroitin sulfate and prevents heparin-induced inhibition of Factor Xa in cell-free assays.¹ Surfen inhibits FGF2 binding to CHO cells (IC₅₀ = 5 μM) and FGF2-induced tube formation by primary mouse endothelial cells. It also inhibits neural differentiation and increases pluripotency in murine embryonic stem cells (mESCs) when used at a concentration of 5 μM.²

References

1. Schuksz, M., Fuster, M.M., Brown, J.R., *et al.* Surfen, a small molecule antagonist of heparan sulfate. *Proc. Natl. Acad. Sci. USA* **105(35)**, 13075-13080 (2008).
2. Huang, M.L., Michalak, A.L., Fisher, C.J., *et al.* Small molecule antagonist of cell surface glycosaminoglycans restricts mouse embryonic stem cells in a pluripotent state. *Stem Cells* **36(1)**, 45-54 (2018).

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, 06/26/2024

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