

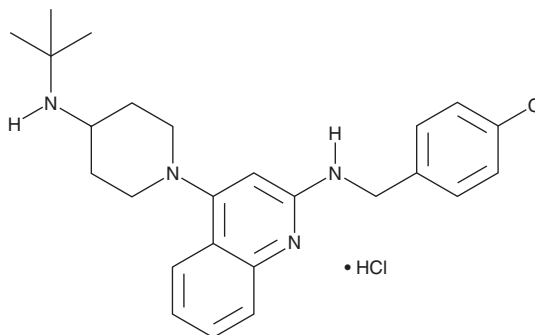
PRODUCT INFORMATION



GNS561 (hydrochloride)

Item No. 45746

CAS Registry No.: 1914148-73-4
Formal Name: N-[(4-chlorophenyl)methyl]-4-[4-[(1,1-dimethylethyl)amino]-1-piperidinyl]-2-quinolinamine, monohydrochloride
Synonym: Ezurpimtrostat
MF: $C_{25}H_{31}ClN_4 \cdot HCl$
FW: 459.5
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GNS561 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the GNS561 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. GNS561 (hydrochloride) is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

GNS561 is an inhibitor of autophagy.¹ It inhibits palmitoyl-protein thioesterase 1 (PPT1) and autophagic flux, increases lysosomal zinc(II) levels, and induces apoptosis in HepG2 hepatocellular carcinoma (HCC) cells. GNS561 reduces viability in a panel of 17 human cancer cells ($IC_{50}s = 0.22-7.27 \mu M$) and reduces alpha fetoprotein serum levels and tumor volume and weight in a patient-derived xenograft (PDX) mouse model of HCC. It also inhibits replication of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) in Vero E6 cells ($EC_{50} = 0.04 \mu M$) and reduces lung viral titers in a mouse model of SARS-CoV-2 infection.²

References

1. Brun, S., Bestion, E., Raymond, E., *et al.* GNS561, a clinical-stage PPT1 inhibitor, is efficient against hepatocellular carcinoma *via* modulation of lysosomal functions. *Autophagy* **18**(3), 678-694 (2022).
2. Bestion, E., Zandi, K., Belouzard, S., *et al.* GNS561 exhibits potent antiviral activity against SARS-CoV-2 through autophagy inhibition. *Viruses* **14**(1), 132 (2022).

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, 07/17/2026

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