

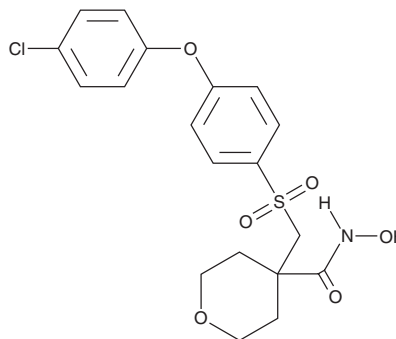
PRODUCT INFORMATION



CTS-1027

Item No. 33219

CAS Registry No.: 193022-04-7
Formal Name: 4-[[[4-(4-chlorophenoxy)phenyl]sulfonyl]methyl]tetrahydro-N-hydroxy-2H-pyran-4-carboxamide
Synonyms: Ro 113-0830, RS 130830
MF: C₁₉H₂₀ClNO₆S
FW: 425.9
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

CTS-1027 is supplied as a solid. A stock solution may be made by dissolving the CTS-1027 in the solvent of choice, which should be purged with an inert gas. CTS-1027 is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

CTS-1027 is an inhibitor of matrix metalloproteinases (MMPs).¹ It is selective for MMP-2, MMP-9, and MMP-13 (K_is= 0.22, 0.55, and 0.52 nM, respectively) over MMP-3 and MMP-1 (K_is= 9.3 and 590 nM, respectively). CTS-1027 (10 mg/kg) reduces hepatocyte apoptosis and bile infarcts and increases survival in a mouse model of bile duct ligation-induced cholestatic liver injury.²

References

1. Kim, S.-H., Pudzianowski, A.T., Leavitt, K.J., *et al.* Structure-based design of potent and selective inhibitors of collagenase-3 (MMP-13). *Bioorg. Med. Chem. Lett.* **15**(4), 1101-1106 (2005).
2. Kahraman, A., Bronk, S.F., Cazanave, S., *et al.* Matrix metalloproteinase inhibitor, CTS-1027, attenuates liver injury and fibrosis in the bile duct-ligated mouse. *Hepatol. Res.* **39**(8), 805-813 (2009).

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, 10/27/2022

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