

PRODUCT INFORMATION

(S)-Glycyl-H-1152 (hydrochloride)

Item No. 13332

CAS Registry No.: 913844-45-8

Formal Name: 2-amino-1-[(3S)-hexahydro-3-methyl-4-[(4-methyl-5-isoquinolinyl)sulfonyl]-1H-1,4-diazepin-1-yl]-ethanone, dihydrochloride

Synonym: Rho Kinase Inhibitor IV

MF: $C_{18}H_{24}N_4O_3S \cdot 2HCl$

FW: 449.4

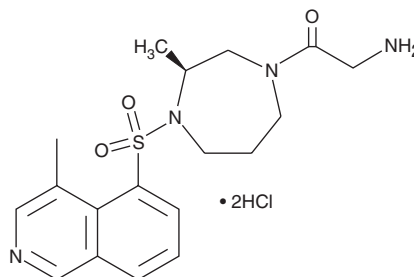
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 216, 280, 328 nm

Supplied as: A solution in methanol

Storage: $-20^{\circ}C$

Stability: $\geq XX$ years



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

Laboratory Procedures

(S)-Glycyl-H-1152 (hydrochloride) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. Solvents such as DMSO and dimethyl formamide (DMF) can be used. The solubility of (S)-glycyl-H-1152 (hydrochloride) in DMSO is approximately 10 mg/ml and approximately 5 mg/ml in DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of (S)-glycyl-H-1152 (hydrochloride) is needed, it can be prepared by evaporating the methanol and directly dissolving the neat oil in aqueous buffers. The solubility of (S)-glycyl-H-1152 (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Two Rho-associated kinases (ROCK), ROCK1 and ROCK2, act downstream of the G protein Rho to regulate cytoskeletal stability. The ROCKs play important roles in diverse cellular functions including cell adhesion and proliferation, smooth muscle contraction, and stem cell renewal.¹⁻³ Glycyl-H-1152 is a selective and potent ROCK inhibitor ($IC_{50} = 11.8$ nM for ROCK2).⁴ It is a glycyated isoquinoline compound derived from the therapeutically-important ROCK inhibitor HA-1077 (fasudil) and exhibits better specificity. Thus, it poorly inhibits Ca^{2+} /calmodulin-dependent kinase type II, protein kinase (PK) G, and Aurora A ($IC_{50} = 2.57, 2.35,$ and 3.26 μM , respectively) as well as PKA or PKC ($IC_{50} \geq 10$ μM for each).⁴ The potency of Glycyl-H-1152 is superior to that of other ROCK inhibitors, including Y-27632 ($K_i = 220$ nM) and HA-1077 ($IC_{50} = 158$ nM).^{4,5}

References

1. Olson, M.F. *Curr. Opin. Cell Biol.* **20**(2), 242-248 (2008).
2. Narumiya, S., Tanji, M., and Ishizaki, T. *Cancer Metastasis Rev.* **28**(1-2), 65-76 (2009).
3. Watanabe, K., Ueno, M., Kamiya, D., et al. *Nat. Biotechnol.* **25**(6), 681-686 (2007).
4. Tamura, M., Nakao, H., Yoshizaki, H., et al. *Biochim. Biophys. Acta* **1754**(1-2), 245-252 (2005).
5. Ishizaki, T., Uehata, M., Tamechika, I., et al. *Mol. Pharmacol.* **57**(5), 976-983 (2000).

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, 07/19/2021

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