

# PRODUCT INFORMATION



## (S)-CCG-1423

Item No. 20107

CAS Registry No.: 2319939-24-5

Formal Name: (S)-N-((1-((4-chlorophenyl)amino)-1-oxopropan-2-yl)oxy)-3,5-bis(trifluoromethyl)benzamide

MF:  $C_{18}H_{13}ClF_6N_2O_3$

FW: 454.8

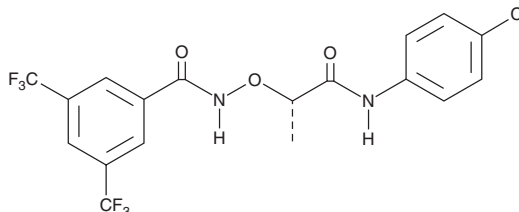
Purity:  $\geq 98\%$

UV/Vis.:  $\lambda_{max}$ : 249 nm

Supplied as: A 1 mg/ml solution in ethanol

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

Stability:  $\geq 2$  years



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

### Laboratory Procedures

(S)-CCG-1423 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of (S)-CCG-1423 in these solvents is approximately 10 mg/ml.

(S)-CCG-1423 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanol solution of (S)-CCG-1423 should be diluted with the aqueous buffer of choice. (S)-CCG-1423 has a solubility of 0.15 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

(S)-CCG-1423 is a stereoisomer of CCG-1423 (Item No. 10010350), a Rho inhibitor that blocks signaling through myocardin-related transcription factor A (MRTF-A) and serum response factor (SRF).<sup>1,2</sup> The S-isomer of CCG-1423 inhibits MRTF-A-dependent cellular events, including SRF-mediated gene expression and cell migration, more potently than the R-isomer.<sup>2</sup>

### References

1. Evelyn, C. R., Wade, S. M., Wang, Q., *et al.* CCG-1423: A small-molecule inhibitor of RhoA transcriptional signaling. *Mol. Cancer Ther.* **6**(8), 2249-2260 (2007).
2. Watanabe, B., Minami, S., Ishida, H., *et al.* Stereospecific inhibitory effects of CCG-1423 on the cellular events mediated by myocardin-related transcription factor A. *PLoS One* **10**(8), (2015).

#### 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, 11/12/2025

#### 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