

PRODUCT INFORMATION



Chroman 1 (hydrochloride hydrate)

Item No. 34682

Formal Name: (S)-N-(2-(2-(dimethylamino)ethoxy)-4-(1H-pyrazol-4-yl)phenyl)-6-methoxychromane-3-carboxamide, dihydrochloride hydrate

Synonym: (S)-Chroman 1

MF: C₂₄H₂₈N₄O₄ • 2HCl [XH₂O]

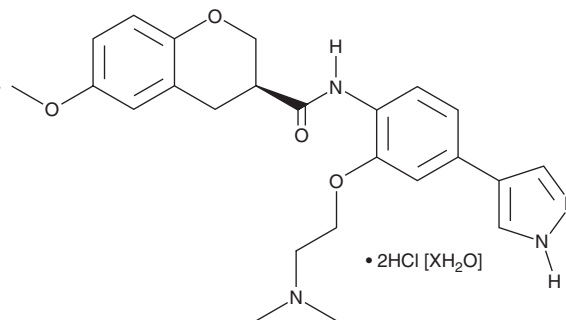
FW: 527.4

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

Chroman 1 (hydrochloride hydrate) is supplied as a solid. A stock solution may be made by dissolving the chroman 1 (hydrochloride hydrate) in the solvent of choice, which should be purged with an inert gas. Chroman 1 (hydrochloride hydrate) is soluble in the organic solvent DMSO. It is also soluble in water. The solubility of chroman 1 (hydrochloride hydrate) in DMSO and water is approximately 100 mM and 5 mM, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Chroman 1 is an inhibitor of Rho-associated kinase 2 (ROCK2; IC₅₀ = <1 nM).¹ It is selective for ROCK2 over protein kinase A (PKA) and Cdc42-binding kinase (MRCKα; IC₅₀s = 1,740 and 127 nM, respectively). Chroman 1 inhibits phosphorylation of myosin light chain (MLC) in a cell-based assay (IC₅₀ = <4 nM).

Reference

1. Chen, Y.T., Vojkovsky, T., Fang, X., *et al.* Asymmetric synthesis of potent chroman-based Rho kinase (ROCK-II) inhibitors. *Med. Chem. Commun.* **2**(1), 73-75 (2011).

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, 12/12/2022

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