

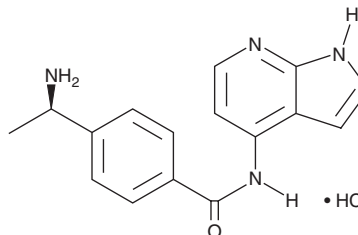
PRODUCT INFORMATION



Y-33075 (hydrochloride)

Item No. 36216

CAS Registry No.: 471843-75-1
Formal Name: 4-[(1R)-1-aminoethyl]-N-1H-pyrrolo[2,3-b]pyridin-4-ylbenzamide, monohydrochloride
Synonyms: RKI-983, SNJ-1656, Y-39983
MF: C₁₆H₁₆N₄O • HCl
FW: 316.8
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 227, 298 nm
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

Y-33075 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the Y-33075 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Y-33075 (hydrochloride) is soluble in DMSO.

Description

Y-33075 is an inhibitor of Rho-associated kinase 2 (ROCK2; IC₅₀ = 3.6 nM).¹ It is selective for ROCK2 over PKC and calcium/calmodulin-dependent protein kinase II (CaMKII; IC₅₀s = 420 and 810 nM, respectively). Topical application of Y-33075 (0.05% v/v) reduces intraocular pressure in rabbit and cynomolgus monkeys. Y-33075 (10 and 100 μM) increases axonal regeneration at the crush site in a cat model of surgically-induced optic nerve damage.² It also inhibits demyelination and reduces the incidence of relapse episodes in a mouse model of chronic relapsing experimental autoimmune encephalomyelitis (EAE).³

References

1. Tokushige, H., Inatani, M., Nemoto, S., *et al.* Effects of topical administration of Y-39983, a selective rho-associated protein kinase inhibitor, on ocular tissues in rabbits and monkeys. *Invest. Ophthalmol. Vis. Sci.* **48(7)**, 3216-3222 (2007).
2. Sagawa, H., Terasaki, H., Nakamura, M., *et al.* A novel ROCK inhibitor, Y-39983, promotes regeneration of crushed axons of retinal ganglion cells into the optic nerve of adult cats. *Exp. Neurol.* **205(1)**, 230-240 (2007).
3. Gao, C., Huang, L., Long, Y., *et al.* Y-39983, a selective Rho-kinase inhibitor, attenuates experimental autoimmune encephalomyelitis via inhibition of demyelination. *Neuroimmunomodulation* **20(6)**, 334-340 (2013).

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, 10/25/2022

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