

PRODUCT INFORMATION



GMP Y-27632 (hydrochloride) Ancillary Material

Item No. 44675

CAS Registry No.: 129830-38-2
Formal Name: *trans*-4-[(1R)-1-aminoethyl]-N-4-pyridinyl-cyclohexanecarboxamide, dihydrochloride

MF: $C_{14}H_{21}N_3O \bullet 2HCl$

FW: 320.0

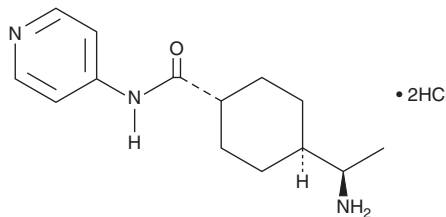
Purity: $\geq 98\%$

Supplied as: An amorphous solid

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

Stability: **batch specific** Please see certificate of analysis for retest dating

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



Description

Y-27632 is a potent, ATP-competitive inhibitor of ROCKs.¹ In cell and gene engineering applications, it has been shown to facilitate cell proliferation, migration, differentiation, and viability.

This product has been manufactured and tested in accordance with GMP principles set forth by the ICH and FDA and meets the expectations for commercial cell, gene, and tissue-engineered products in accordance with USP <1043> "Ancillary Materials for Cell, Gene, and Tissue-engineered Products."²

References

1. Davies, S.P., Reddy, H., Caivano, M., *et al.* Specificity and mechanism of action of some commonly used protein kinase inhibitors. *Biochem. J.* **351**(1), 95-105 (2000).
2. **United States Pharmacopeia.** <1043> Ancillary Materials for Cell, Gene, and Tissue-Engineered Products. In: *USP-NF*. Rockville, MD: United States Pharmacopeial Convention

WARNING

Ancillary Material intended for further processing in accordance with USP <1043>

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, 06/04/2026

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