

PRODUCT INFORMATION



Y-27632 (hydrochloride) Ancillary Material

Item No. 44885

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

MF: C₁₄H₂₁N₃O • 2HCl

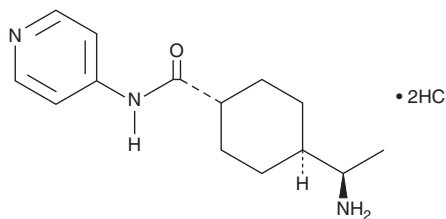
FW: 320.0

Purity: ≥98%

Supplied as: An amorphous solid

Storage: -20°C

Stability: *batch specific Please see the Certificate of Analysis for retest dating.*



Description

Y-27632 is a potent, ATP-competitive inhibitor of Rho-associated kinases (ROCKs).¹ It has been used to enhance cell survival and cloning efficiency, as well as to prevent dissociation-induced apoptosis, also known as anoikis, in pluripotent stem cells.²⁻⁴ Y-27632 (hydrochloride) Ancillary Material is intended for use in cell, gene, and tissue engineering applications.

This product has been tested in accordance with GMP principles and meets the expectations for clinical cell, gene, and tissue-engineered products in accordance with USP <1043> "Ancillary Materials for Cell, Gene, and Tissue-engineered Products".⁵

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. Wang, X., Lin, G., Martins-Taylor, K., *et al.* Inhibition of caspase-mediated anoikis is critical for basic fibroblast growth factor-sustained culture of human pluripotent stem cells. *J. Biol. Chem.* **284**(49), 34054-34064 (2009).
3. Gandham, V.D., Maddala, R.L., Rao, V., *et al.* Effects of Y27632 on keratinocyte procurement and wound healing. *Clin. Exp. Dermatol.* **38**(7), 782-786 (2013).
4. Fuentes, E.O., Leemhuis, J., Stark, G.B., *et al.* Rho kinase inhibitors Y27632 and H1152 augment neurite extension in the presence of cultured Schwann cells. *J. Brachial Plex. Peripher. Nerve Inj.* **3**(19) (2008).
5. **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 for use in manufacturing, processing, or packaging of cell-based or gene-engineered products. **Not intended for direct clinical administration.**

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, 03/06/2026

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