

PRODUCT INFORMATION



RY785

Item No. 19813

CAS Registry No.: 1393748-80-5
Formal Name: 3-methoxy- α -methyl-N-[2-(4-thiazolyl)-1H-benzimidazol-6-yl]-benzenepropanamide

MF: C₂₁H₂₀N₄O₂S

FW: 392.5

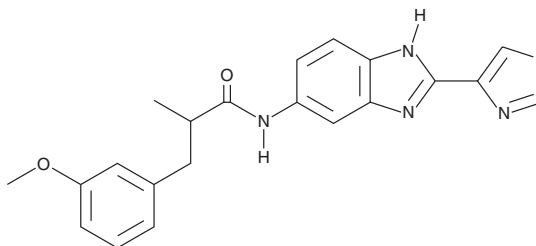
Purity: \geq 98%

UV/Vis.: λ_{\max} : 210, 245, 317 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

RY785 is supplied as a crystalline solid. A stock solution may be made by dissolving the RY785 in the solvent of choice. RY785 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of RY785 in ethanol and DMSO is approximately 15 mg/ml and the solubility in DMF is approximately 20 mg/ml.

Description

RY785 is a selective inhibitor of K_v2.1 voltage-gated potassium channels (IC₅₀ = 0.05 μ M) that does not affect Ca_v2 calcium channels.¹ It displays over 100-fold selectivity for K_v2.1 over Ca_v1.2 (IC₅₀ = 17 μ M).¹

Reference

- Herrington, J., Solly, K., Ratliff, K., *et al.* Identification of novel and selective K_v2 channel inhibitors. *Mol. Pharmacol.* **80**(6), 959-964 (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, 12/19/2022

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