

PRODUCT INFORMATION



ASK1 Inhibitor 10

Item No. 28512

CAS Registry No.: 1005775-56-3
Formal Name: 4-(1,1-dimethylethyl)-N-[6-(1H-imidazol-1-yl)imidazo[1,2-a]pyridin-2-yl]-benzamide, dihydrochloride

MF: C₂₁H₂₁N₅O • 2HCl
FW: 432.4

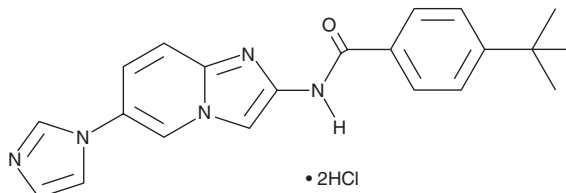
Purity: ≥95%

UV/Vis.: λ_{max}: 253, 335 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

ASK1 inhibitor 10 is supplied as a solid. A stock solution may be made by dissolving the ASK1 inhibitor 10 in the solvent of choice, which should be purged with an inert gas. ASK1 inhibitor 10 is soluble in the organic solvent DMSO at a concentration of approximately 40 mg/ml.

Description

ASK1 inhibitor 10 is an orally bioavailable inhibitor of apoptosis signal-regulating kinase 1 (ASK1; IC₅₀ = 14 nM).¹ It is selective for ASK1 over ASK2 (IC₅₀ = 510 nM) as well as MEKK1, TAK-1, IKKβ, ERK1, JNK1, p38α, GSK3β, PKCθ, and B-RAF (IC₅₀s = >10,000 nM for all). It inhibits streptozotocin-induced increases in JNK and p38 phosphorylation in INS-1 pancreatic β cells in a concentration-dependent manner.

Reference

1. Terao, Y., Suzuki, H., Yoshikawa, M., *et al.* Design and biological evaluation of imidazo[1,2-a]pyridines as novel and potent ASK1 inhibitors. *Bioorg. Med. Chem. Lett.* **22(24)**, 7326-7329 (2012).

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, 10/03/2022

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