

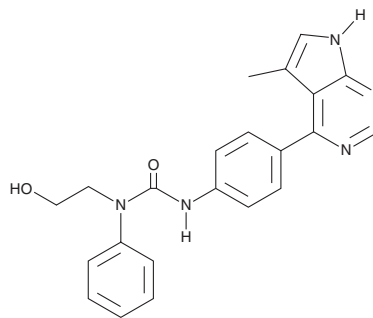
PRODUCT INFORMATION



SR 7826

Item No. 32976

CAS Registry No.: 1219728-20-7
Formal Name: N-(2-hydroxyethyl)-N'-[4-(5-methyl-7H-pyrrolo[2,3-d]pyrimidin-4-yl)phenyl]-N-phenyl-urea
MF: C₂₂H₂₁N₅O₂
FW: 387.4
Purity: ≥98%
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

SR 7826 is supplied as a solid. A stock solution may be made by dissolving the SR 7826 in the solvent of choice, which should be purged with an inert gas. SR 7826 is soluble (≥10 mg/ml) in DMSO.

Description

SR 7826 is an inhibitor of LIM kinase 1 (LIMK1; IC₅₀ = 43 nM).¹ It is selective for LIMK1 over Rho-associated kinase 1 (ROCK1) and ROCK2 (IC₅₀s = 6,565 and 5,536 nM, respectively) and a panel of 60 kinases at 1 μM but does inhibit serine/threonine kinase 16 (STK16) by greater than 80% at 1 μM. SR 7826 inhibits the phosphorylation of cofilin induced by hepatocyte growth factor (HGF) in A7r5 cells (IC₅₀ = 470 nM). It reduces the migration and invasion of PC3 prostate cancer cells when used at a concentration of 1 μM. SR 7826 (1 μM) inhibits electric field stimulation-, phenylephrine-, methoxamine-, norepinephrine-, or U-46619-induced contractions of prostate strips isolated from patients with prostate cancer.² *In vivo*, SR 7826 (10 mg/kg, p.o.) increases apical and basal thin spine density of CA1 pyramidal neurons in the hAPP-J20 mouse model of Alzheimer's disease.³

References

1. Yin, Y., Zheng, K., Eid, N., *et al.* Bis-aryl urea derivatives as potent and selective LIM kinase (Limk) inhibitors. *J. Med. Chem.* **58**(4), 1846-1861 (2015).
2. Yu, Q., Gratzke, C., Wang, Y., *et al.* Inhibition of human prostate smooth muscle contraction by the LIM kinase inhibitors, SR7826 and LIMKi3. *Br. J. Pharmacol.* **175**(11), 2077-2096 (2018).
3. Henderson, B.W., Greathouse, K., Ramdas, R., *et al.* Pharmacologic inhibition of LIMK1 provides dendritic spine resilience against β-amyloid. *Sci. Signal* **12**(587), eaaw9318 (2019).

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, 08/11/2025

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