

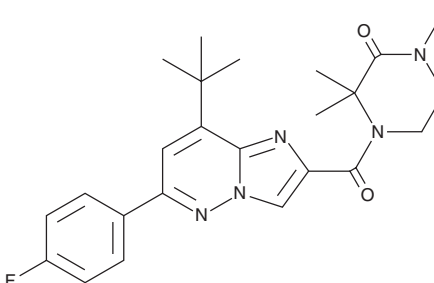
PRODUCT INFORMATION



I-191

Item No. 37553

CAS Registry No.: 1690172-25-8
Formal Name: 4-[[8-(1,1-dimethylethyl)-6-(4-fluorophenyl)imidazo[1,2-b]pyridazin-2-yl]carbonyl]-3,3-dimethyl-2-piperazinone
MF: C₂₃H₂₆FN₅O₂
FW: 423.5
Purity: ≥98%
UV/Vis.: λ_{max}: 245, 283 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

I-191 is supplied as a solid. A stock solution may be made by dissolving the I-191 in the solvent of choice, which should be purged with an inert gas. I-191 is soluble in DMSO and methanol.

Description

I-191 is an antagonist of proteinase-activated receptor 2 (PAR2; IC₅₀ = 72.4 nM in HT-29 cells expressing the human receptor).¹ It inhibits trypsin-induced calcium release in HT-29 cells (IC₅₀ = 200 nM for the human receptor). I-191 (0.1, 1, and 10 μM) reduces the migration of HT-29 cells in a scratch assay. It decreases the expression of mRNA encoding chemokine (C-X-C motif) ligand 8 (CXCL8) in, and secreted IL-8 levels from, HT-29 cells induced by the PAR2 agonist 2f-LIGRL-NH₂ when used at a concentration of 1 and 10 μM. *Ex vivo*, I-191 (10 μM) inhibits tissue factor-induced clotting using isolated human whole blood.²

References

1. Jiang, Y., Yau, M.K., Lim, J., *et al.* A potent antagonist of protease-activated receptor 2 that inhibits multiple signaling functions in human cancer cells. *J. Pharmacol. Exp. Ther.* **364**(2), 246-257 (2018).
2. Iyer, A., Humphries, T.L.R., Owens, E.P., *et al.* PAR2 activation on human kidney tubular epithelial cells induces tissue factor synthesis, that enhances blood clotting. *Front. Physiol.* **12**, 615428 (2021).

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, 10/10/2023

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