

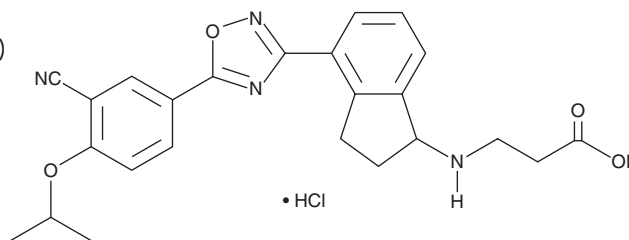
PRODUCT INFORMATION



RP 001 (hydrochloride)

Item No. 16636

CAS Registry No.: 1781880-34-9
Formal Name: N-[4-[5-[3-cyano-4-(1-methylethoxy)phenyl]-1,2,4-oxadiazol-3-yl]-2,3-dihydro-1H-inden-1-yl]-β-alanine, monohydrochloride
MF: C₂₄H₂₄N₄O₄ • HCl
FW: 468.9
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

RP 001 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the RP 001 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. RP 001 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

RP 001 is a potent sphingosine-1-phosphate receptor 1 (S1P₁) agonist (EC₅₀ = 9 pM).¹ It is competitive with W146 (Item No. 10009109), a selective S1P₁ antagonist, and induces internalization and polyubiquitination of S1P₁ in a dose-dependent manner. RP001 induces dose-dependent lymphopenia in mice (EC₅₀ = 0.03 mg/kg). It reduces S1P₁-EGFP expression *ex vivo* on lymphocytes in the isolated lymph node of *Edg1^{eGFP/eGFP}* mice but has no effect *in vivo*. RP 001 also completely inhibits *M. stadtmanae* (MSS) rechallenge-induced enlargement of B cell-rich tertiary lymphoid tissues (TLTs) and reduces neutrophil accumulation in the airways of a mouse model of experimental hypersensitivity pneumonitis.²

References

1. Cahalan, S.M., Gonzalez-Cabrera, P.J., Sarkisyan, G., *et al.* Actions of a picomolar short-acting S1P₁ agonist in S1P₁-eGFP knock-in mice. *Nat. Chem. Biol.* **7**(5), 254-256 (2011).
2. Huppé, C.A., Lecours, P.B., Lechasseur, A., *et al.* A sphingosine-1-phosphate receptor 1 agonist inhibits tertiary lymphoid tissue reactivation and hypersensitivity in the lung. *Mucosal. Immunol.* **11**(1), 112-119 (2018).

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