

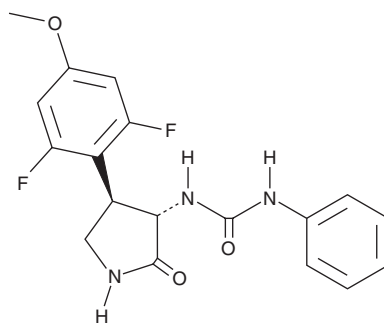
PRODUCT INFORMATION



BMS 986235

Item No. 45369

CAS Registry No.: 2253947-47-4
Formal Name: N-[(3S,4R)-4-(2,6-difluoro-4-methoxyphenyl)-2-oxo-3-pyrrolidinyl]-N'-phenyl-urea
Synonym: LAR-1219
MF: C₁₈H₁₇F₂N₃O₃
FW: 361.3
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

BMS 986235 is supplied as a solid. A stock solution may be made by dissolving the BMS 986235 in the solvent of choice, which should be purged with an inert gas. BMS 986235 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

BMS 986235 is a formyl peptide receptor 2 (FPR2) agonist.¹ It selectively induces calcium mobilization in HEK293 cells overexpressing human FPR2 over FPR1 (EC₅₀s = 0.41 and 2,800 nM, respectively). BMS 986235 inhibits chemotaxis of HL-60 human promyeloblast cells and induces phagocytosis of opsonized zymosan by isolated mouse peritoneal macrophages. *In vivo*, BMS 986235 inhibits neutrophil infiltration into the lungs in a mouse model of LPS-induced lung inflammation. BMS 986235 reduces left ventricle remodeling, scar length, and heart mass in a mouse model of myocardial infarction when administered at a dose of 0.3 mg/kg on days 4-28 post-occlusion. It also reduces right and left ventricle damage and inflammation in a mouse model of Duchenne muscular dystrophy.²

References

1. Asahina, Y., Wurtz, N.R., Arakawa, K., *et al.* Discovery of BMS-986235/LAR-1219: A potent formyl peptide receptor 2 (FPR2) selective agonist for the prevention of heart failure. *J. Med. Chem.* **63**(17), 9003-9019 (2020).
2. Novak, J.S., Lischin, A., Uapinyoying, P., *et al.* Failure to resolve inflammation contributes to juvenile onset cardiac damage in a mouse model of Duchenne muscular dystrophy. *Cell Death Dis.* **16**(1), 505 (2025).

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, 04/22/2026

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