

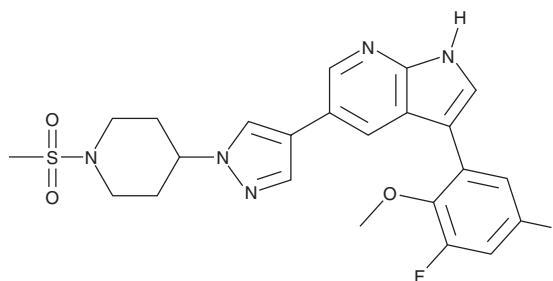
PRODUCT INFORMATION



ARN3261

Item No. 45729

CAS Registry No.: 1613711-28-6
Formal Name: 3-(3,5-difluoro-2-methoxyphenyl)-5-[1[1-(methylsulfonyl)-4-piperidinyl]-1H-pyrazol-4-yl]-1H-pyrrolo[2,3-b]pyridine
Synonyms: GRN-300
MF: C₂₃H₂₃F₂N₅O₃S
FW: 487.5
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

ARN3261 is supplied as a solid. A stock solution may be made by dissolving the ARN3261 in the solvent of choice, which should be purged with an inert gas. ARN3261 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

ARN3261 is an inhibitor of salt-inducible kinase 2 (SIK2).¹ It induces apoptosis in OVCAR-8, SKOV3, and OC 316 ovarian cancer cells and inhibits growth in ten ovarian cancer cell lines (IC₅₀s = 0.8-2.6 μM). ARN3261 enhances sensitivity to carboplatin (Item No. 13112) in both carboplatin-sensitive and -resistant ovarian cancer cells and increases carboplatin-induced DNA damage and apoptosis in OVCAR-8, SKOV3, and OC 316 cells. It reduces tumor weight in a SKOV3 mouse xenograft model when administered at a dose of 40 mg/kg with carboplatin.

Reference

1. Fan, D., Yang, H., Mao, W., *et al.* A novel salt inducible kinase 2 inhibitor, ARN-3261, sensitizes ovarian cancer cell lines and xenografts to carboplatin. *Cancers* **13**, 446 (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, 05/27/2026

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