

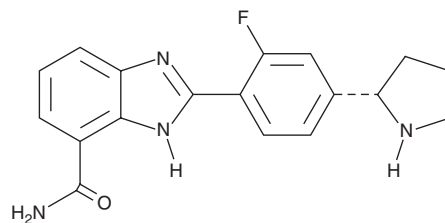
PRODUCT INFORMATION



A-966492

Item No. 26378

CAS Registry No.: 934162-61-5
Formal Name: 2-[2-fluoro-4-[(2S)-2-pyrrolidinyl]phenyl]-1H-benzimidazole-7-carboxamide
MF: C₁₈H₁₇FN₄O
FW: 324.4
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

A-966492 is supplied as a solid. A stock solution may be made by dissolving the A-966492 in the solvent of choice, which should be purged with an inert gas. A-966492 is soluble in DMSO.

Description

A-966492 is a dual inhibitor of poly(ADP-ribose) polymerase 1 (PARP1) and PARP2 ($K_s = 1$ and 1.5 nM, respectively).¹ It reduces tumor growth by 46 and 92% in an MX-1 mouse xenograft model when administered at doses of 100 and 200 mg/kg per day, respectively. A-966492 also reduces tumor growth in a B16/F10 murine melanoma model when administered in combination with temozolomide (Item No. 14163).

Reference

1. Penning, T.D., Zhu, G.D., Gong, J., *et al.* Optimization of phenyl-substituted benzimidazole carboxamide poly(ADP-ribose) polymerase inhibitors: Identification of (S)-2-(2-fluoro-4-(pyrrolidin-2-yl)phenyl)-1H-benzimidazole-4-carboxamide (A-966492), a highly potent and efficacious inhibitor. *J. Med. Chem.* **53**(8), 3142-3153 (2010).

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