

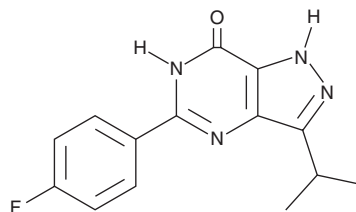
PRODUCT INFORMATION



NPD-2975

Item No. 39466

CAS Registry No.: 3032452-65-3
Formal Name: 5-(4-fluorophenyl)-1,6-dihydro-3-(1-methylethyl)-7H-pyrazolo[4,3-d]pyrimidin-7-one
MF: C₁₄H₁₃FN₄O
FW: 272.3
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 292 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

NPD-2975 is supplied as a solid. A stock solution may be made by dissolving the NPD-2975 in the solvent of choice, which should be purged with an inert gas. NPD-2975 is soluble in methanol and DMSO.

Description

NPD-2975 is an antiparasitic agent.¹ It reduces the proliferation of *T. brucei* (IC₅₀ = 63.1 nM) but not *T. cruzi*, *L. infantum*, *P. falciparum*, or human MRC-5 fibroblasts (IC₅₀s = >6,000 nM for all). *In vivo*, NPD-2975 (50 mg/kg twice per day) reduces peripheral blood parasite levels below detection limits and increases survival in a mouse model of *T. brucei* infection.

Reference

1. Zheng, Y., van den Kerkhof, M., van der Meer, T., *et al.* Discovery of 5-phenylpyrazolopyrimidinone analogs as potent antitrypanosomal agents with in vivo efficacy. *J. Med. Chem.* **66(15)**, 10252-10264 (2023).

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, 10/20/2025

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