

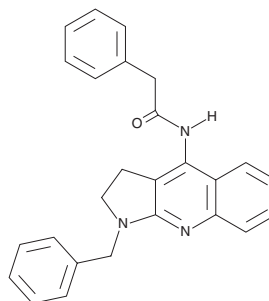
PRODUCT INFORMATION



PGP-4008

Item No. 20901

CAS Registry No.: 365565-02-2
Formal Name: N-[2,3-dihydro-1-(phenylmethyl)-1H-pyrrolo[2,3-b]quinolin-4-yl]-benzeneacetamide
MF: C₂₆H₂₃N₃O
FW: 393.5
Purity: ≥95%
UV/Vis.: λ_{max}: 210, 256, 347 nm
Supplied as: A crystalline 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

PGP-4008 is supplied as a crystalline solid. A stock solution may be made by dissolving the PGP-4008 in the solvent of choice, which should be purged with an inert gas. PGP-4008 is soluble in organic solvents such as ethanol and DMSO. The solubility of PGP-4008 in these solvents is approximately 5 and 50 mM, respectively.

Description

PGP-4008 is a selective inhibitor of P-glycoprotein (P-gp) that does not affect the activity of multidrug resistance-related protein 1 (MRP1).^{1,2} It is effective *in vitro* and, when delivered intraperitoneally, *in vivo*, inhibiting tumor growth when given with doxorubicin (Item No. 15007).^{1,2} PGP-4008 also enhances the ability of docetaxel (Item No. 11637) to inhibit growth and drive apoptosis in ovarian cancer cells.³

References

1. Lee, B.D., French, K.J., Zhuang, Y., *et al.* Development of a syngeneic *in vivo* tumor model and its use in evaluating a novel P-glycoprotein modulator, PGP-4008. *Oncol. Res.* **14(1)**, 49-60 (2003).
2. Lee, B.D., Li, Z., French, K.J., *et al.* Synthesis and evaluation of dihydropyrroloquinolines that selectively antagonize P-glycoprotein. *J. Med. Chem.* **47(6)**, 1413-1422 (2004).
3. Miettinen, S., Gr nman, S. and Ylikomi, T. Inhibition of P-glycoprotein-mediated docetaxel efflux sensitizes ovarian cancer cells to concomitant docetaxel and SN-38 exposure. *Anticancer Drugs* **20(4)**, 267-276 (2009).

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, 11/14/2022

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