

PRODUCT INFORMATION



DQP1105

Item No. 31121

CAS Registry No.: 380560-89-4
Formal Name: 5-(4-bromophenyl)-3-(1,2-dihydro-6-methyl-2-oxo-4-phenyl-3-quinolinyl)-4,5-dihydro- γ -oxo-1H-pyrazole-1-butanoic acid

MF: C₂₉H₂₄BrN₃O₄

FW: 558.4

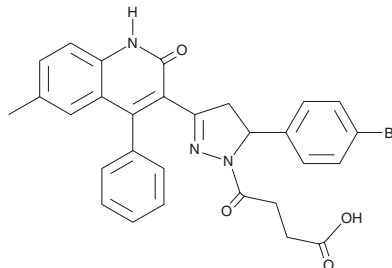
Purity: \geq 98%

UV/Vis.: λ_{max} : 232 nm

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

DQP1105 is supplied as a solid. A stock solution may be made by dissolving the DQP1105 in the solvent of choice, which should be purged with an inert gas. DQP1105 is soluble in the organic solvent DMSO at a concentration of approximately 50 mg/ml.

Description

DQP1105 is an NMDA receptor antagonist.¹ It selectively inhibits glutamate-induced currents in *Xenopus* oocytes expressing rat NR2C and NR2D subunit-containing NMDA receptors (IC₅₀s = 8.5 and 2.7 μ M, respectively) over NR2A, NR2B, NRA1, and NRK2 subunit-containing receptors (IC₅₀s = 206, 121, 198, and 153 μ M, respectively).

Reference

1. Acker, T.M., Yuan, H., Hansen, K.B., *et al.* Mechanism for noncompetitive inhibition by novel GluN2C/D N-methyl-D-aspartate receptor subunit-selective modulators. *Mol. Pharmacol.* **80(5)**, 782-795 (2011).

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, 11/30/2022

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