

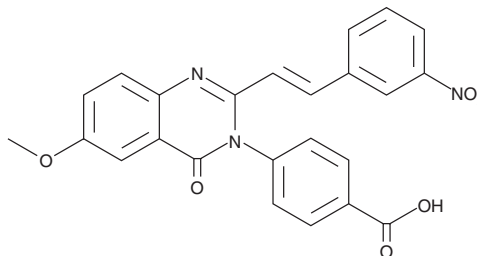
PRODUCT INFORMATION



QNZ 46

Item No. 14857

CAS Registry No.: 1237744-13-6
Formal Name: 4-[6-methoxy-2-[(1E)-2-(3-nitrophenyl)ethenyl]-4-oxo-3(4H)-quinazoliny]-benzoic acid
MF: C₂₄H₁₇N₃O₆
FW: 443.4
Purity: ≥98%
UV/Vis.: λ_{max}: 228, 346 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

QNZ 46 is supplied as a crystalline solid. A stock solution may be made by dissolving the QNZ 46 in the solvent of choice, which should be purged with an inert gas. QNZ 46 is soluble in 0.01 M DMSO.

Description

QNZ 46 is an antagonist of NR2C- and NR2D-subunit containing NMDA receptors (IC₅₀s = 6 and 3 μM, respectively, for receptors expressed in *Xenopus* oocytes).¹ It is selective for NR2C- and NR2D-subunit containing NMDA receptors over receptors containing NR2A and NR2B subunits as well as over the AMPA receptor GluR1 (IC₅₀s = 229, >300, and >300 μM, respectively). QNZ 46 acts in a noncompetitive and voltage-independent fashion that requires glutamate but not glycine binding for its activity.²

References

1. Mosley, C.A., Acker, T.M., Hansen, K.B., *et al.* Quinazolin-4-one derivatives: A novel class of noncompetitive NR2C/D subunit-selective N-methyl-D-aspartate receptor antagonists. *J. Med. Chem.* **53(15)**, 5476-5490 (2010).
2. Hansen, K.B. and Traynelis, S.F. Structural and mechanistic determinants of a novel site for noncompetitive inhibition of GluN2D-containing NMDA receptors. *J. Neurosci.* **31(10)**, 3650-3661 (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, 11/14/2022

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