

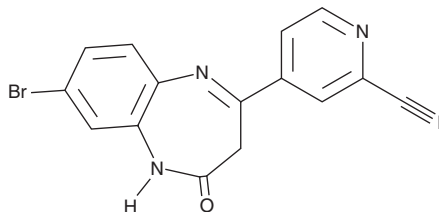
PRODUCT INFORMATION



MNI137

Item No. 45486

CAS Registry No.: 946619-21-2
Formal Name: 4-(8-bromo-2,3-dihydro-2-oxo-1H-1,5-benzodiazepin-4-yl)-2-pyridinecarbonitrile
MF: $C_{15}H_9BrN_4O$
FW: 341.2
Purity: $\geq 98\%$
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

MNI137 is supplied as a solid. A stock solution may be made by dissolving the MNI137 in the solvent of choice, which should be purged with an inert gas. MNI137 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

MNI137 is a negative allosteric modulator of metabotropic glutamate receptor 2 (mGluR2).¹ It selectively inhibits glutamate-induced calcium mobilization in HEK293 cells expressing human mGluR2 ($IC_{50} = 11.4$ nM) over HEK293 cells expressing rat mGluR1, -5, -7, or -8 and those expressing human mGluR4 at 10 μM .

Reference

1. Hemstapat, K., Da Costa, H., Nong, Y., et al. A novel family of potent negative allosteric modulators of group II metabotropic glutamate receptors. *J. Pharmacol. Exp. Ther.* **322**(1), 254-264 (2007).

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, 05/04/2026

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