

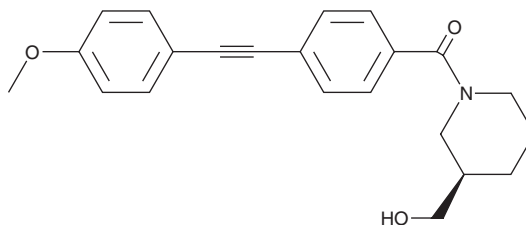
PRODUCT INFORMATION



ML-289

Item No. 15195

CAS Registry No.: 1382481-79-9
Formal Name: [3R-(hydroxymethyl)-1-piperidinyl]
[4-[2-(4-methoxyphenyl)ethynyl]
phenyl]-methanone
Synonym: VU0463597
MF: C₂₂H₂₃NO₃
FW: 349.4
Purity: ≥98%
UV/Vis.: λ_{max}: 303 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

ML-289 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-289 in the solvent of choice, which should be purged with an inert gas. ML-289 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

Glutamate, the major excitatory neurotransmitter in the CNS, activates eight known subtypes of metabotropic glutamate receptors (mGluRs). Highly selective modulators designed to act at allosteric sites within the receptor are being developed to preferentially target individual mGluR subtypes in order to determine their distinct pharmacological roles. ML-289 is a negative allosteric modulator of mGlu₃ (IC₅₀ = 0.66 μM) that displays 15-fold selectivity over mGlu₂ and inactivity against mGlu₅.¹

References

1. Sheffler, D.J., Wenthur, C.J., Bruner, J.A., *et al.* Development of a novel, CNS-penetrant, metabotropic glutamate receptor 3 (mGlu₃) NAM probe (ML289) derived from a closely related mGlu₅ PAM. *Bioorg. Med. Chem. Lett.* **22**(12), 3921-3925 (2012).

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, 12/14/2022

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