

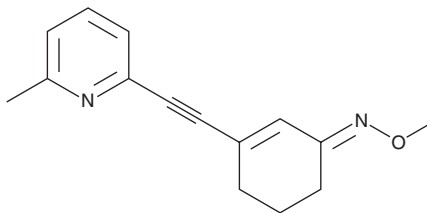
PRODUCT INFORMATION



ABP 688

Item No. 30929

CAS Registry No.: 924298-51-1
Formal Name: 3-[2-(6-methyl-2-pyridinyl)ethynyl]-2-cyclohexen-1-one, O-methyloxime
MF: C₁₅H₁₆N₂O
FW: 240.3
Purity: ≥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

ABP 688 is supplied as a solid. A stock solution may be made by dissolving the ABP 688 in the solvent of choice, which should be purged with an inert gas. ABP 688 is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

ABP 688 is a non-competitive antagonist of metabotropic glutamate receptor 5 (mGluR5; $K_i = 3.5$ nM in a radioligand binding assay using L(tk-) cell membranes expressing human receptors).¹ It is selective for mGluR5 receptors over a panel of CNS G protein-coupled receptors, ion channels, and transporters at 10 μ M. ABP 688 inhibits quisqualate-induced phosphoinositol accumulation in and glutamate-induced calcium release from L(tk-) cells expressing human mGluR5 (IC_{50} s = 2.4 and 2.3 nM, respectively). Radiolabeled forms of ABP 688 have been used as PET tracers for *in vivo* imaging of mGluR5 receptors in rodents.^{1,2}

References

- Hintermann, S., Vranesic, I., Allgeier, H., *et al.* ABP688, a novel selective and high affinity ligand for the labeling of mGlu5 receptors: Identification, *in vitro* pharmacology, pharmacokinetic and biodistribution studies. *Bioorg. Med. Chem.* **15**(2), 903-914 (2007).
- Elmenhorst, D., Minuzzi, L., Aliaga, A., *et al.* *In vivo* and *in vitro* validation of reference tissue models for the mGluR5 ligand [¹¹C]ABP688. *J. Cereb. Blood Flow Metab.* **30**(8), 1538-1549 (2010).

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