

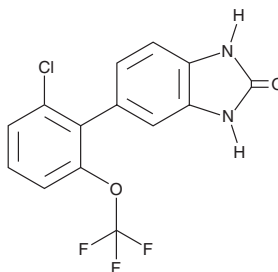
PRODUCT INFORMATION



JNJ-55511118

Item No. 29212

CAS Registry No.: 2036081-86-2
Formal Name: 5-[2-chloro-6-(trifluoromethoxy)phenyl]-1,3-dihydro-2H-benzimidazol-2-one
MF: C₁₄H₈ClF₃N₂O₂
FW: 328.7
Purity: ≥98%
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

JNJ-55511118 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-55511118 in the solvent of choice, which should be purged with an inert gas. JNJ-55511118 is soluble in the organic solvent DMSO.

Description

JNJ-55511118 is a negative modulator of AMPA receptors containing transmembrane AMPA receptor regulatory protein $\gamma 8$ (TARP- $\gamma 8$).¹ It inhibits glutamate-induced calcium flux in HEK293F cells co-transfected with TARP- $\gamma 8$ and either GluA1o, GluA1i, GluA2i, GluA3o, or GluA4o (IC_{50} s = 11.22, 12.3, 7.41, 38.02, and 15.85 nM, respectively). JNJ-55511118 is selective for GluA1o receptors containing TARP- $\gamma 8$ over GluA1o receptors containing TARP- $\gamma 2$, - $\gamma 3$, - $\gamma 4$, or - $\gamma 7$ (IC_{50} s = >10 μ M for all). It reduces peak glutamate-induced currents in acutely dissociated murine hippocampal neurons to 60.7% of control cells when used at a concentration of 1 μ M. JNJ-55511118 (1 μ M) reduces field excitatory postsynaptic potentials (fEPSPs) in the hippocampal CA1 region from wild-type, but not TARP- $\gamma 8$ knockout, mice. It also inhibits corneal kindling-induced seizures in mice (ED_{50} = 3.7 mg/kg).

Reference

1. Maher, M.P., Wu, N., Ravula, S., *et al.* Discovery and characterization of AMPA receptor modulators selective for TARP- $\gamma 8$. *J. Pharmacol. Exp. Ther.* **357**(2), 394-414 (2016).

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, 12/13/2022

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