

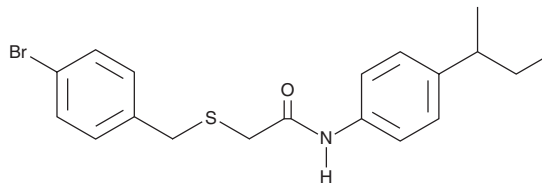
PRODUCT INFORMATION



AZ 12216052

Item No. 29460

CAS Registry No.: 1290628-31-7
Formal Name: 2-[[[(4-bromophenyl)methyl]thio]-N-[4-(1-methylpropyl)phenyl]-acetamide
MF: C₁₉H₂₂BrNOS
FW: 392.4
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 248 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

AZ 12216052 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZ 12216052 in the solvent of choice, which should be purged with an inert gas. AZ 12216052 is soluble in the organic solvent DMSO.

Description

AZ 12216052 is a positive allosteric modulator of metabotropic glutamate receptor 8 (mGluR8).¹ It potentiates glutamate-induced [³⁵S]GTPγS binding in GHEK cells expressing human mGluR8b (EC₅₀ = 1 μM). In a cell-based model of Parkinson's disease, AZ 12216052 inhibits MPP⁺-induced decreases in the viability of undifferentiated SH-SY5Y cells when used at concentrations ranging from 0.001 to 1 μM and in retinoic acid-differentiated SH-SY5Y cells at 0.1 and 1 μM.² AZ 12216052 (10 mg/kg) increases the time spent in the open areas of the elevated zero maze and reduces the acoustic startle response in mice, indicating anxiolytic-like activity in response to avoidable and unavoidable anxiogenic stimuli, respectively.¹

References

1. Duvoisin, R.M., Pfankuch, T., Wilson, J.M., *et al.* Acute pharmacological modulation of mGluR8 reduces measures of anxiety. *Behav. Brain Res.* **212**(2), 168-173 (2010).
2. Jantas, D., Greda, A., Golda, S., *et al.* Neuroprotective effects of metabotropic glutamate receptor group II and III activators against MPP(+)-induced cell death in human neuroblastoma SH-SY5Y cells: The impact of cell differentiation state. *Neuropharmacology* **83**, 36-53 (2014).

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, 12/06/2022

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