

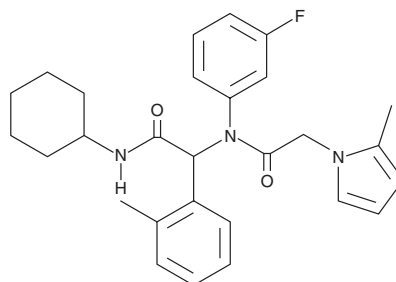
PRODUCT INFORMATION



AGI-5198

Item No. 14624

CAS Registry No.: 1355326-35-0
Formal Name: N-[2-(cyclohexylamino)-1-(2-methylphenyl)-2-oxoethyl]-N-(3-fluorophenyl)-2-methyl-1H-imidazole-1-acetamide
MF: C₂₇H₃₁FN₄O₂
FW: 462.6
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

AGI-5198 is supplied as a crystalline solid. A stock solution may be made by dissolving the AGI-5198 in the solvent of choice, which should be purged with an inert gas. AGI-5198 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AGI-5198 in these solvents is approximately 5, 10, and 20 mg/ml, respectively.

AGI-5198 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AGI-5198 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AGI-5198 has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Isocitrate dehydrogenases (IDHs) are nicotinamide adenine dinucleotide (NAD⁺) and NAD phosphate (NADP⁺)-dependent enzymes in the tricarboxylic acid cycle that catalyze oxidative decarboxylation of isocitrate producing α-ketoglutarate (2-OG) and carbon dioxide. IDH1 and IDH2 are mutated in >70% of lower grade gliomas.¹ The most common mutations map to arginine residues in the catalytic pockets of IDH1 (R132) or IDH2 (R140 and R172) and impart new gain of function catalytic activity leading to the NADPH-dependent conversion of 2-OG to 2-hydroxyglutarate (2-HG).²⁻⁴ AGI-5198 is a potent, selective inhibitor of IDH1 R132H and R132C mutants *in vitro* with IC₅₀ values of 0.07 and 0.16 μM, respectively, but not wild-type IDH1, wild-type IDH2, or IDH2 mutants (IC₅₀s > 100 μM).⁵ AGI-5198 has been shown to have anti-tumor efficacy in the TS603 glioma cell line and to lower tumor 2-HG production in HT1080 and U87MG cells with IC₅₀ values of 0.48 and 0.07 μM, respectively.⁶ In R132H-IDH1 glioma xenografts, AGI-5198 (450 mg/kg/day) caused 50-60% growth inhibition over a treatment period of three weeks with no affect in the growth of IDH1 wild-type glioma xenografts.⁵ Under conditions of near complete 2-HG inhibition, AGI-5198 can induce demethylation of histone H3K9me3 and expression of genes associated with gliogenic differentiation.⁵

References

1. Turcan, S., Rohle, D., Goenka, A., *et al. Nature* **483(7390)**, 479-483 (2012).
2. Reitman, Z.J. and Yan, H. *J. Natl. Cancer Inst.* **102(13)**, 932-941 (2010).
3. Dang, L., White, D.W., Gross, S., *et al. Nature* **462(7274)**, 739-744 (2009).
4. Koivunen, P., Lee, S., Duncan, C.G., *et al. Nature* **483(7390)**, 484-488 (2012).
5. Rohle, D., Popovici-Muller, J., Palaskas, N., *et al. Science* **340(6132)**, 626-630 (2013).
6. Popovici-Muller, J., Saunders, J.O., Salituro, F.G., *et al. ACS Med. Chem. Lett.* **3(10)**, 850-855 (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, 01/16/2024

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