

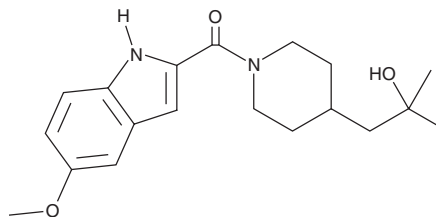
PRODUCT INFORMATION



ASP9521

Item No. 27308

CAS Registry No.: 1126084-37-4
Formal Name: [4-(2-hydroxy-2-methylpropyl)-1-piperidiny] (5-methoxy-1H-indol-2-yl)-methanone
MF: C₁₉H₂₆N₂O₃
FW: 330.4
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 294 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

ASP9521 is supplied as a crystalline solid. A stock solution may be made by dissolving the ASP9521 in the solvent of choice. ASP9521 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of ASP9521 in these solvents is approximately 10 mg/ml.

ASP9521 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ASP9521 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ASP9521 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ASP9521 is an inhibitor of 17β-hydroxysteroid dehydrogenase type 5/aldo-keto reductase 1C3 (17β-HSD5/AKR1C3; IC₅₀ = 120 nM).¹ It is selective for 17β-HSD5/AKR1C3 over AKR1C2 (IC₅₀ = 20 μM). ASP9521 inhibits conversion of androstenedione into testosterone by 17β-HSD5/AKR1C3 (IC₅₀ = 11 and 49 nM for human and cynomolgus monkey enzymes, respectively). It inhibits androstenedione-dependent production of prostate specific androgen (PSA) in and proliferation of LNCaP cells expressing 17β-HSD5/AKR1C3. ASP9521 (3 and 10 mg/kg) inhibits androstenedione-induced intratumor testosterone production in a CWR22R prostate cancer mouse xenograft model.

Reference

1. Kikuchi, A., Furutani, T., Azami, H., *et al.* In vitro and in vivo characterisation of ASP9521: A novel, selective, orally bioavailable inhibitor of 17β-hydroxysteroid dehydrogenase type 5 (17βHSD5; AKR1C3). *Invest New Drugs* **32**(5), 860-870 (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, 09/30/2022

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