

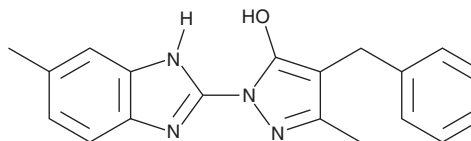
PRODUCT INFORMATION



HUHS015

Item No. 32915

CAS Registry No.: 1453097-13-6
Formal Name: 3-methyl-1-(6-methyl-1H-benzimidazol-2-yl)-4-(phenylmethyl)-1H-pyrazol-5-ol
MF: C₁₉H₁₈N₄O
FW: 318.4
Purity: ≥98%
UV/Vis.: λ_{max}: 301 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

HUHS015 is supplied as a crystalline solid. A stock solution may be made by dissolving the HUHS015 in the solvent of choice, which should be purged with an inert gas. HUHS015 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HUHS015 in DMSO is approximately 1 mg/ml. HUHS015 is slightly soluble in ethanol and DMF.

HUHS015 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HUHS015 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. HUHS015 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HUHS015 is an inhibitor of prostate cancer antigen-1 (PCA-1; IC₅₀ = 0.67 μM), also known as α-ketoglutarate-dependent dioxygenase alkB homolog 3 (ALKBH3).¹ It inhibits proliferation of DU145 human prostate cancer cells by 54 and 81% in 2D dish and 3D agar assays, respectively when used at a concentration of 10 μM. HUHS015 (32 mg/kg) reduces tumor growth in a DU145 mouse xenograft model.

Reference

1. Nakao, S., Mabuchi, M., Shimizu, T., *et al.* Design and synthesis of prostate cancer antigen-1 (PCA-1/ALKBH3) inhibitors as anti-prostate cancer drugs. *Bioorg. Med. Chem. Lett.* **24(4)**, 1071-1074 (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, 10/05/2022

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