

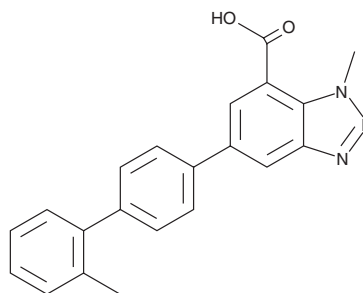
PRODUCT INFORMATION



AG-636

Item No. 36122

CAS Registry No.: 1623416-31-8
Formal Name: 1-methyl-5-(2'-methyl[1,1'-biphenyl]-4-yl)-1H-benzotriazole-7-carboxylic acid
MF: $C_{21}H_{17}N_3O_2$
FW: 343.4
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 259 nm
Supplied as: A 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

AG-636 is supplied as a solid. A stock solution may be made by dissolving the AG-636 in the solvent of choice, which should be purged with an inert gas. AG-636 is soluble in the organic solvent DMSO. AG-636 is slightly soluble in water.

Description

AG-636 is a dihydroorotate dehydrogenase (DHODH) inhibitor ($IC_{50} = 17 \text{ nM}$).¹ It increases levels of ureidosuccinic acid (Item No. 34126) and DHO, pyrimidine biosynthesis metabolites upstream of DHODH, in a panel of six cell lines when used at a concentration of $1 \mu\text{M}$. AG-636 (100 mg/kg) induces complete tumor stasis in an OCI-Ly19 diffuse large B cell lymphoma (DLBCL) mouse xenograft model and complete tumor regression in a Z-138 ibrutinib-resistant mantle cell lymphoma mouse xenograft model. It also increases DHO levels and reduces pyrimidine levels in Z-138 and JeKo-1 lymphoma mouse xenograft models.

Reference

1. McDonald, G., Chubukov, V., Coco, J., *et al.* Selective vulnerability to pyrimidine starvation in hematologic malignancies revealed by AG-636, a novel clinical-stage inhibitor of dihydroorotate dehydrogenase. *Mol. Cancer Ther.* **19**(12), 2502-2515 (2020).

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, 12/19/2022

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