

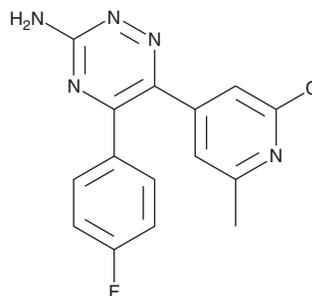
PRODUCT INFORMATION



AZD 4635

Item No. 26322

CAS Registry No.: 1321514-06-0
Formal Name: 6-(2-chloro-6-methyl-4-pyridinyl)-5-(4-fluorophenyl)-1,2,4-triazin-3-amine
Synonyms: HTL-1071, Imaradenant
MF: C₁₅H₁₁ClFN₅
FW: 315.7
Purity: ≥98%
UV/Vis.: λ_{max}: 273, 304 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

AZD 4635 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 4635 in the solvent of choice, which should be purged with an inert gas. AZD 4635 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AZD 4635 in these solvents is approximately 30 mg/ml.

AZD 4635 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AZD 4635 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AZD 4635 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AZD 4635 is an adenosine A₂ receptor antagonist (K_i = 1.7 nM for the human receptor).¹ It inhibits adenosine-induced cAMP production in CHO cells expressing the human A₂ receptor (IC₅₀ = 0.79 μM) and reverses suppression of IFN-γ secretion induced by 5'-N-ethylcarboxamidoadenosine (NECA; Item No. 21420) in CD8⁺ T cells. AZD 4635 reduces tumor growth when administered alone and in combination with checkpoint inhibitors in syngeneic mouse tumor models.

Reference

1. Borodovsky, A., Wang, Y., Ye, M., *et al.* Abstract 5580: Preclinical pharmacodynamics and antitumor activity of AZD4635, a novel adenosine 2A receptor inhibitor that reverses adenosine mediated T cell suppression. *Cancer Res.* **77(13 Supp.)**, 5580 (2017).

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, 10/10/2022

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