

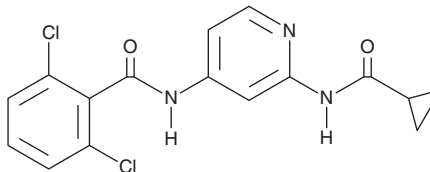
PRODUCT INFORMATION



GDC-046

Item No. 36601

CAS Registry No.: 1258292-64-6
Formal Name: 2,6-dichloro-N-[2-[(cyclopropylcarbonyl)amino]-4-pyridinyl]-benzamide
MF: C₁₆H₁₃Cl₂N₃O₂
FW: 350.2
Purity: ≥98%
UV/Vis.: λ_{max}: 242 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

GDC-046 is supplied as a solid. A stock solution may be made by dissolving the GDC-046 in the solvent of choice, which should be purged with an inert gas. GDC-046 is soluble in DMSO.

Description

GDC-046 is a tyrosine kinase 2 (TYK2) inhibitor ($K_i = 4.8$ nM).¹ It is selective for TYK2 over JAK1, -2, and -3 (K_i s = 83.8, 27.6, and 253 nM, respectively). GDC-046 selectively inhibits TYK2-mediated IL-12-induced phosphorylation of STAT4 over JAK1-mediated IL-6-induced phosphorylation of STAT3 and JAK2-mediated erythropoietin-induced phosphorylation of STAT5 in cell-based assays (EC_{50} s = 380, 2,000, and 1,700 nM, respectively).

Reference

1. Liang, J., van Abbema, A., Balazs, M., et al. Lead optimization of a 4-aminopyridine benzamide scaffold to identify potent, selective, and orally bioavailable TYK2 inhibitors. *J. Med. Chem.* **56**(11), 4521-4536 (2013).

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, 12/20/2022

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