

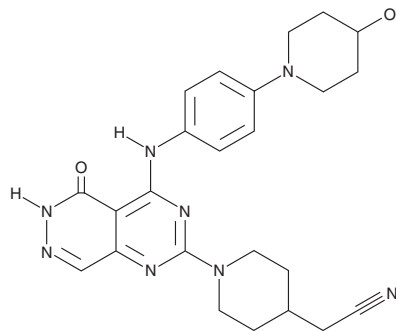
PRODUCT INFORMATION



Gusacitinib

Item No. 36554

CAS Registry No.: 1425381-60-7
Formal Name: 1-[5,6-dihydro-4-[[4-(4-hydroxy-1-piperidiny)phenyl]amino]-5-oxopyrimido[4,5-d]pyridazin-2-yl]-4-
Synonym: ASN-002
MF: C₂₄H₂₈N₈O₂
FW: 460.5
Purity: ≥98%
UV/Vis.: λ_{max}: 267 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

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

Description

Gusacitinib is a dual inhibitor of JAKs (IC₅₀s = 46, 4, and 11 nM for JAK1-3, respectively) and Syk family kinases (IC₅₀s = 5 and 8 nM for SyK and TYK2, respectively).¹

Reference

1. Bissonnette, R., Maari, C., Forman, S., *et al.* The oral Janus kinase/spleen tyrosine kinase inhibitor ASN002 demonstrates efficacy and improves associated systemic inflammation in patients with moderate-to-severe atopic dermatitis: Results from a randomized double-blind placebo-controlled study. *Br. J. Dermatol.* **181**(4), 733-742 (2019).

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