

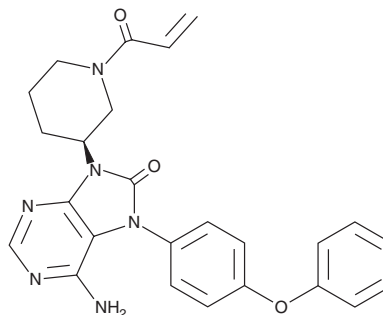
PRODUCT INFORMATION



ONO-4059 analog

Item No. 18924

CAS Registry No.: 1351635-67-0
Formal Name: 6-amino-7,9-dihydro-9-[(3S)-1-(1-oxo-2-propen-1-yl)-3-piperidinyl]-7-(4-phenoxyphenyl)-8H-purin-8-one
Synonym: GS-4059 analog
MF: C₂₅H₂₄N₆O₃
FW: 456.5
Purity: ≥98%
UV/Vis.: λ_{max}: 273 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

ONO-4059 analog is supplied as a crystalline solid. A stock solution may be made by dissolving the ONO-4059 analog in the solvent of choice, which should be purged with an inert gas. ONO-4059 analog is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of ONO-4059 analog in these solvents is approximately 20 and 25 mg/ml, respectively.

ONO-4059 analog is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ONO-4059 analog should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ONO-4059 analog has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ONO-4059 analog is a derivative of the Bruton's tyrosine kinase (BTK) inhibitor ONO-4059 (Item No. 32957).

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, 11/23/2022

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