

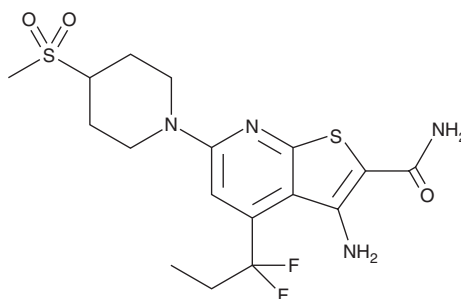
PRODUCT INFORMATION



BI-605906

Item No. 19184

CAS Registry No.: 960293-88-3
Formal Name: 3-amino-4-(1,1-difluoropropyl)-6-[4-(methylsulfonyl)-1-piperidiny]-thieno[2,3-b]pyridine-2-carboxamide
MF: C₁₇H₂₂F₂N₄O₃S₂
FW: 432.5
Purity: ≥98%
UV/Vis.: λ_{max}: 244, 306, 347, 360 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

BI-605906 is supplied as a crystalline solid. A stock solution may be made by dissolving the BI-605906 in the solvent of choice, which should be purged with an inert gas. BI-605906 is soluble in organic solvents such as DMSO and acetonitrile. The solubility of BI-605906 in these solvents is approximately 30 and 1 mg/ml, respectively.

Description

BI-605906 is an inhibitor of inhibitor of kappa B kinase subunit β (IKKβ; IC₅₀ = 380 nM).¹ It is selective for IKKβ over insulin-like growth factor 1 (IGF1; IC₅₀ = 7.6 μM) and over 100 kinases in a panel up to 10 μM. BI-605906 partially inhibits IL-1-stimulated activation of IKKε/TBK1 *in vitro*. BI-605906 also inhibits TNF-α-dependent IκB degradation and expression of the proinflammatory cytokines IL-6, IL-1β, and C-X-C motif ligand 1/2 (CXCL1/2) without affecting lipogenic gene expression or glucose production in murine primary hepatocytes.²

References

1. Clark, K., Peggie, M., Plater, L., *et al.* Novel cross-talk within the IKK family controls innate immunity. *Biochem. J.* **434**(1), 93-104 (2011).
2. Cameron, A.R., Morrison, V.L., Levin, D., *et al.* Anti-inflammatory effects of metformin irrespective of diabetes status. *Circ. Res.* **119**(5), 652-665 (2016).

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, 11/29/2022

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