

PRODUCT INFORMATION



TG101209

Item No. 14696

CAS Registry No.: 936091-14-4
Formal Name: N-(1,1-dimethylethyl)-3-[[5-methyl-2-[[4-(4-methyl-1-piperazinyl)phenyl]amino]-4-pyrimidinyl]amino]-benzenesulfonamide

MF: C₂₆H₃₅N₇O₂S

FW: 509.7

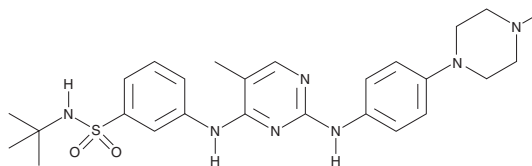
Purity: ≥98%

UV/Vis.: λ_{max}: 282 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

TG101209 is supplied as a crystalline solid. A stock solution may be made by dissolving the TG101209 in the solvent of choice. TG101209 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of TG101209 in these solvents is approximately 0.12, 12.5, and 16 mg/ml, respectively.

TG101209 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TG101209 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TG101209 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

TG101209 is a potent inhibitor of the tyrosine kinase janus kinase (JAK) 2, FMS-like tyrosine kinase 3, proto-oncogene RET, and JAK3 (IC₅₀ = 6, 25, 17, and 169 nM, respectively).¹ Through these effects, it induces cell cycle arrest and apoptosis in leukemic cell lines and CD45⁺ myeloma cells.^{1,2}

References

1. Pardanani, A., Lasho, T., Levine, R.L., *et al.* TG101209, a small molecule JAK2-selective kinase inhibitor potently inhibits myeloproliferative disorder-associated JAK2V617F and MPLW515L/K mutations. *Leukemia* **21**(8), 1658-1668 (2007).
2. Ramakrishnan, V., Kimlinger, T., Haug, J., *et al.* TG101209, a novel JAK2 inhibitor, has significant *in-vitro* activity in multiple myeloma and displays preferential cytotoxicity for CD45⁺ myeloma cells. *Am. J. Hematol.* **85**(9), 675-686 (2010).

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