

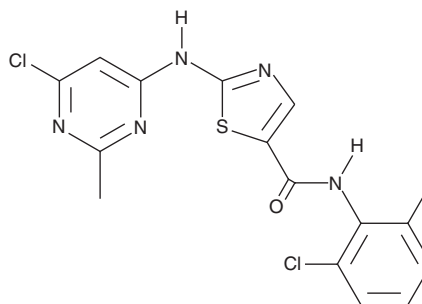
PRODUCT INFORMATION



CAY10697

Item No. 17806

CAS Registry No.: 302964-08-5
Formal Name: N-(2-chloro-6-methylphenyl)-2-
[[6-chloro-2-methyl-4-pyrimidinyl]
amino]-5-thiazolecarboxamide
MF: C₁₆H₁₃Cl₂N₅OS
FW: 394.3
Purity: ≥98%
UV/Vis.: λ_{max}: 319 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

CAY10697 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10697 in the solvent of choice, which should be purged with an inert gas. CAY10697 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10697 in these solvents is approximately 2 and 5 mg/ml, respectively.

CAY10697 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Dasatinib (Item No. 11498) is a potent inhibitor of receptor and non-receptor tyrosine kinases, including some drug resistant mutant forms.¹⁻⁴ It has potential therapeutic value in diseases that are characterized by elevated levels of these kinases, including some forms of cancer and fibrotic disease.^{1,2,5,6} CAY10697 is an intermediate in the synthesis of dasatinib.⁷

References

1. Lombardo, L.J., Lee, F.Y., Chen, P., *et al. J. Med. Chem.* **47**(27), 6658-6661 (2004).
2. Das, J., Chen, P., Norris, D., *et al. J. Med. Chem.* **49**(23), 6819-6823 (2006).
3. Davis, M.I., Hunt, J.P., Herrgard, S., *et al. Nat. Biotechnol.* **29**(11), 1046-1051 (2011).
4. Carter, T.A., Wodicka, L.M., Shah, N.P., *et al. Proc. Natl. Acad. Sci. USA* **102**(31), 11011-11016 (2005).
5. Distler, J.H.W. and Distler, O. *Rheumatology* **47**, 10-11 (2008).
6. McFarland, K.L. and Wetzstein, G.A. *Cancer Control* **16**(2), 132-140 (2009).
7. Deadman, B.J., Hopkin, M.D., Baxendale, I.R., *et al. Org. Biomol. Chem.* **11**, 1766-1800 (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, 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