

PRODUCT INFORMATION



CGP 77675 (hydrate)

Item No. 21089

Formal Name: 1-[2-[4-[4-amino-5-(3-methoxyphenyl)-7H-pyrrolo[2,3-d]pyrimidin-7-yl]phenyl]ethyl]-4-piperidinol, hydrate

MF: C₂₆H₂₉N₅O₂ • XH₂O

FW: 443.5

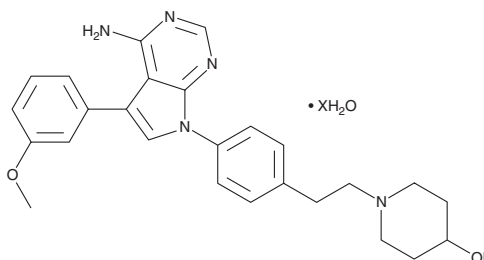
Purity: ≥98%

UV/Vis.: λ_{max}: 246, 291 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

CGP 77675 (hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the CGP 77675 (hydrate) in the solvent of choice, which should be purged with an inert gas. CGP 77675 (hydrate) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

CGP 77675 is an inhibitor of Src family kinases (SFKs) that blocks the phosphorylation of peptide substrates and autophosphorylation of purified c-Src (IC₅₀s = 5-20 and 40 nM, respectively).¹ It also inhibits the SFKs Lck and c-Yes with low nanomolar IC₅₀ values.¹⁻³ CGP 77675 is used to elucidate the role of SFKs in such processes as bone resorption, tight junction formation in epithelial cells, and allograft survival.^{1,2,4} It can also be used in combination with the GSK3 inhibitor CHIR99021 (Item No. 13122) to maintain mouse embryonic stem cells.⁵

References

1. Missbach, M., Jeschke, M., Feyen, J., *et al.* A novel inhibitor of the tyrosine kinase Src suppresses phosphorylation of its major cellular substrates and reduces bone resorption *in vitro* and in rodent models *in vivo*. *Bone* **24(5)**, 437-449 (1999).
2. Chen, Y.-H., Goodenough, D.A., and Jeansonne, B. Nonreceptor tyrosine kinase c-Yes interacts with occludin during tight junction formation in canine kidney epithelial cells. *Mol. Biol. Cell* **13(4)**, 1227-1237 (2002).
3. M. Šušća, M., N.-H. Luong-Nguyen, N.-H., J. Crespo, J., *et al.* Active recombinant human tyrosine kinase c-Yes: Expression in baculovirus system, purification, comparison to c-Src, and inhibition by a c-Src inhibitor. *Protein Expr. Purif.* **19(1)**, 99-106 (2000).
4. Zhang, Q., Fairchild, R.L., Reich, M.B., *et al.* Inhibition of Src kinases combined with CD40 ligand blockade prolongs murine cardiac allograft survival. *Transplantation* **80(8)**, 1112-1120 (2005).
5. Shimizu, T., Ueda, J., Ho, J.C., *et al.* Dual inhibition of Src and GSK3 maintains mouse embryonic stem cells, whose differentiation is mechanically regulated by Src signaling. *Stem Cells* **30(7)**, 1394-1404 (2012).

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, 12/19/2022

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