

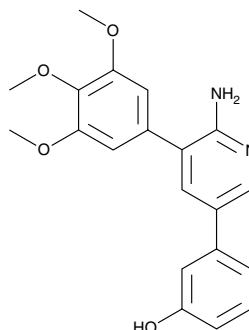
Product Information



K02288

Item No. 16678

CAS Registry No.: 1431985-92-0
Formal Name: 3-[6-amino-5-(3,4,5-trimethoxyphenyl)-3-pyridinyl]-phenol
MF: C₂₀H₂₀N₂O₄
FW: 352.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 255, 325 nm nm



Laboratory Procedures

For long term storage, we suggest that K02288 be stored as supplied at -20°C. It should be stable for at least two years.

K02288 is supplied as a crystalline solid. A stock solution may be made by dissolving the K02288 in the solvent of choice. K02288 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of K02288 in ethanol is approximately 0.1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

K02288 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, K02288 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. K02288 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Activation of bone morphogenetic protein (BMP) type I receptors, also known as activin receptor-like kinases (ALK1-7), leads to the assembly of SMAD complexes, which translocate to the nucleus to induce transcriptional activation important for normal development and tissue repair.¹ K02288 is a 2-aminopyridine-based inhibitor of ALK1 and ALK2 with IC₅₀ values of 1.8 and 1.1 nM, respectively.² It is less selective for ALK3, 4, 5, and 6 subtypes and the type II BMP receptor ActRIIA, demonstrating IC₅₀ values of 34.4, 302, 321, 6.4, and 220 nM, respectively.² K02288 can prevent BMP4-induced SMAD1/5/8 pathway activation *in vitro* (IC₅₀ = 100 nM) without affecting TGF-β signaling. Furthermore, at 8-10 μM, K02288 has been used to induce the dorsalization of zebrafish embryos. Through specific inhibition of BMP signaling, this compound can be used to research stem cell biology and disease models of musculoskeletal dysplasia and cancer.²

References

1. Brivanlou, A.H. and Darnell, J.E., Jr. Signal transduction and the control of gene expression. *Science* **295**, 813-818 (2002).
2. Sanvitale, C.E., Kerr, G., Chaikuad, A., *et al.* A new class of small molecule inhibitor of BMP signaling. *PLoS One* **8**(4), 62721 (2013).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/16678

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 10/28/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com