

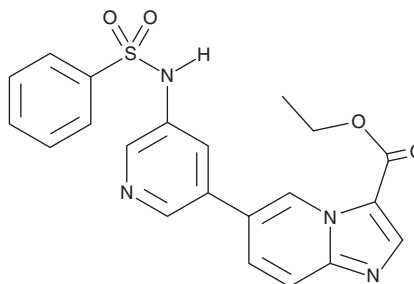
PRODUCT INFORMATION



HS-173

Item No. 19156

CAS Registry No.: 1276110-06-5
Formal Name: 6-[5-[(phenylsulfonyl)amino]-3-pyridinyl]-imidazo[1,2-a]pyridine-3-carboxylic acid, ethyl ester
MF: C₂₁H₁₈N₄O₄S
FW: 422.5
Purity: ≥98%
UV/Vis.: λ_{max}: 246, 297 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

HS-173 is supplied as a crystalline solid. A stock solution may be made by dissolving the HS-173 in the solvent of choice, which should be purged with an inert gas. HS-173 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HS-173 in ethanol is approximately 1 mg/ml and approximately 25 mg/ml in DMSO and DMF.

HS-173 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HS-173 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. HS-173 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

HS-173 is an imidazopyridine analog that acts as a PI3Kα inhibitor with an IC₅₀ value of 0.8 nM.¹ It demonstrates antiproliferative activity in T47D, SK-BR-3, and MCF-7 cells with IC₅₀ values of 0.6, 1.5, and 7.8 μM, respectively.¹ HS-173 is reported to induce apoptosis by arresting the cell cycle at the G₂/M phase and by activating caspases.^{2,3} It has also been shown to block VEGF-induced angiogenesis both *in vitro* and *in vivo*.²

References

1. Kim, O., Jeong, Y., Lee, H., *et al.* Design and synthesis of imidazopyridine analogues as inhibitors of phosphoinositide 3-kinase signaling and angiogenesis. *J. Med. Chem.* **54**(7), 2455-2466 (2011).
2. Lee, H., Jung, K.H., Jeong, Y., *et al.* HS-173, a novel phosphatidylinositol 3-kinase (PI3K) inhibitor, has anti-tumor activity through promoting apoptosis and inhibiting angiogenesis. *Cancer Lett.* **328**(1), 152-159 (2013).
3. Son, M.K., Ryu, Y.-L., Jung, K.H., *et al.* HS-173, a novel PI3K inhibitor, attenuates the activation of hepatic stellate cells in liver fibrosis. *Sci. Rep.* **3**:3470, (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, 10/05/2022

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