

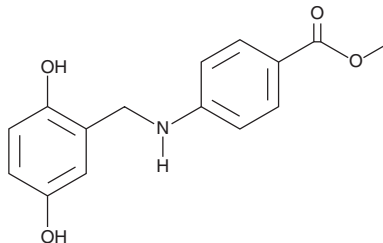
PRODUCT INFORMATION



AG957

Item No. 17259

CAS Registry No.: 140674-76-6
Formal Name: 4-[[[(2,5-dihydroxyphenyl)methyl]amino]-benzoic acid, methyl ester
Synonyms: NSC 654705, Tyrphostin AG957
MF: C₁₅H₁₅NO₄
FW: 273.3
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 308 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

AG957 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG957 in the solvent of choice, which should be purged with an inert gas. AG957 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

AG957 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

Tyrphostins are tyrosine phosphorylation inhibitors that act by inhibiting tyrosine kinases.¹ AG957 is a tyrphostin that targets transforming Bcr-Abl fusion proteins (p185^{Bcr-Abl}, p210^{Bcr-Abl}), as well as normal c-Abl (IC₅₀s = 4.3, 1, and 7.1 μM, respectively, for human proteins).² It also inhibits epidermal growth factor receptor (IC₅₀ = 0.25 μM).² As the constitutively-active 210 kDa Bcr-Abl fusion protein commonly occurs in chronic myelogenous leukemia (CML) cells, AG957 is commonly used to study Bcr-Abl signaling in the CML K562 cell line.^{3,4} AG957 is also used to study signaling through c-Abl.⁵

References

1. Levitzki, A. and Gazit, A. *Science* **267**(5205), 1782-1788 (1995).
2. Anafi, M., Gazit, A., Gilon, C., et al. *J. Biol. Chem.* **267**(8), 4518-4523 (1992).
3. Kaur, G., Gazit, A., Levitzki, A., et al. *Anticancer Drugs* **5**, 213-222 (1994).
4. Jamieson, L., Carpenter, L., Biden, T.J., et al. *J. Biol. Chem.* **274**(7), 3927-3930 (1999).
5. Fu, P., Usatyuk, P.V., Lele, A., et al. *Am. J. Physiol. Regul. Integr. Comp. Physiol.* **308**, L1025-L1038 (2015).

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, 10/24/2022

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