

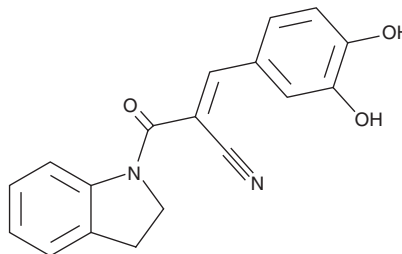
PRODUCT INFORMATION



AG-528

Item No. 29666

CAS Registry No.: 133550-49-9
Formal Name: α E-[(3,4-dihydroxyphenyl)methylene]-2,3-dihydro- β -oxo-1H-indole-1-propanenitrile
Synonym: Tyrphostin AG-528
MF: C₁₈H₁₄N₂O₃
FW: 306.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 248, 363 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

AG-528 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-528 in the solvent of choice, which should be purged with an inert gas. AG-528 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AG-528 in these solvents is approximately 1, 20, and 25 mg/ml, respectively.

Description

AG-528 is a tyrphostin that inhibits EGFR and HER2 (IC₅₀s = 12 and 4.9 μ M, respectively, for inhibition of autophosphorylation).¹ It also inhibits pantothenate kinase 3 (Pank3; IC₅₀ = 3 μ M).² AG-528 inhibits EGF-dependent proliferation of HER14 cells (IC₅₀ = 25 μ M).¹

References

1. Gazit, A., Osherov, N., Posner, I., *et al.* Tyrphostins. 2. Heterocyclic and α -substituted benzylidenemalononitrile tyrphostins as potent inhibitors of EGF receptor and ErbB2/neu tyrosine kinases. *J. Med. Chem.* **34**(6), 1896-1907 (1991).
2. Leonardi, R., Zhang, Y.-M., Yun, M.-K., *et al.* Modulation of pantothenate kinase 3 activity by small molecules that interact with the substrate/allosteric regulatory domain. *Chem. Biol.* **17**(8), 892-902 (2010).

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, 10/17/2022

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