

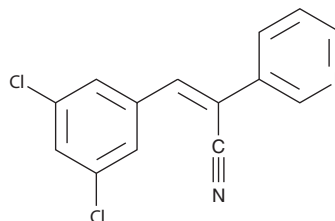
PRODUCT INFORMATION



RG-14620

Item No. 10010310

CAS Registry No.: 136831-49-7
Formal Name: α -[(3,5-dichlorophenyl)methylene]-3-pyridineacetonitrile
MF: C₁₄H₈Cl₂N₂
FW: 275.1
Purity: \geq 98%
UV/Vis.: λ_{max} : 206, 309 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

RG-14620 is supplied as a crystalline solid. A stock solution may be made by dissolving the RG-14620 in the solvent of choice, which should be purged with an inert gas. RG-14620 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of RG-14620 in ethanol is approximately 1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

Protein tyrosine kinase (PTK) inhibitors are potential antiproliferative agents for diseases caused by the hyperactivity of PTKs. Tyrophostins are a class of antiproliferative compounds that act as PTK blockers. PTK inhibitors that preferentially inhibit the epidermal growth factor (EGF) receptor kinase block EGF-dependent cell proliferation.¹ RG-14620 is an inhibitor of EGF receptor kinase with an IC₅₀ value of 3 μ M in HT-22 cells.²

References

1. Gazit, A., Yaish, P., Gilon, C., *et al.* Tyrophostins I: Synthesis and biological activity of protein tyrosine kinase inhibitors. *J. Med. Chem.* **32(10)**, 2344-2352 (1989).
2. Sagara, Y., Ishige, K., Tsai, C., *et al.* Tyrophostins protect neuronal cells from oxidative stress. *J. Biol. Chem.* **277(39)**, 36204-36215 (2002).

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, 05/09/2024

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