

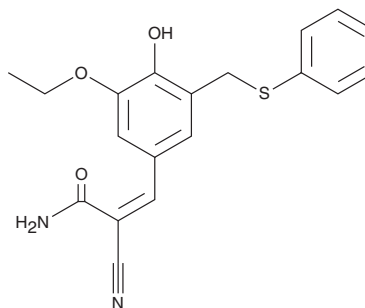
PRODUCT INFORMATION



ST638

Item No. 13337

CAS Registry No.: 107761-24-0
Formal Name: 2-cyano-3-[3-ethoxy-4-hydroxy-5-[(5-phenylthio)methyl]phenyl]-2-propenamide
MF: C₁₉H₁₈N₂O₃S
FW: 354.4
Purity: ≥98%
UV/Vis.: λ_{max}: 252, 363 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

ST638 is supplied as a crystalline solid. A stock solution may be made by dissolving the ST638 in the solvent of choice, which should be purged with an inert gas. ST638 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of ST638 in these solvents is approximately 30 and 50 mg/ml, respectively.

ST638 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ST638 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ST638 has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ST638 is a tyrosine kinase inhibitor (IC₅₀ = 370 nM). It suppresses tyrosine phosphorylation induced by tumor necrosis factor (TNF-α) and phorbol myristate acetate (PMA) in neutrophils and by angiotensin II in cardiac fibroblasts.¹⁻³

References

1. Fuortes, M., Jin, W.-W., and Nathan, C. Adhesion-dependent protein tyrosine phosphorylation in neutrophils treated with tumor necrosis factor. *J. Cell Biol.* **120**(3), 777-784 (1993).
2. Murasawa, S., Mori, Y., Nozawa, Y., *et al.* Role of calcium-sensitive tyrosine kinase Pyk2/CAKb/RAFTK in angiotensin II-induced Ras/ERK signaling. *Hypertension* **32**, 668-675 (1998).
3. Takemasa, H., Imagawa, N., Kawakami-Honda, N., *et al.* Effect of phosphotyrosine proteins on phorbol myristate acetate-induced NADPH oxidase activation in guinea pig peritoneal polymorphonuclear leukocytes. *Biol. Pharm. Bull.* **26**(7), 1009-1012 (2003).

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, 10/19/2022

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