

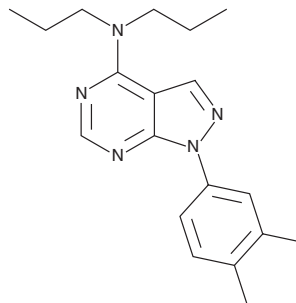
PRODUCT INFORMATION



S3QEL-2

Item No. 18556

CAS Registry No.: 890888-12-7
Formal Name: 1-(3,4-dimethylphenyl)-N,N-dipropyl-1H-pyrazolo[3,4-d]pyrimidin-4-amine
MF: C₁₉H₂₅N₅
FW: 323.4
Purity: ≥98%
UV/Vis.: λ_{max}: 243, 298 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

S3QEL-2 is supplied as a crystalline solid. A stock solution may be made by dissolving the S3QEL-2 in the solvent of choice, which should be purged with an inert gas. S3QEL-2 is soluble in the organic solvent dimethyl formamide (DMF) at a concentration of approximately 10 mg/ml. S3QEL-2 is also slightly soluble in DMSO.

S3QEL-2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, S3QEL-2 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. S3QEL-2 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

S3QEL-2 is a cell-permeable, selective suppressor of electron leak from mitochondrial respiratory complex III, inhibiting site III_{Qo} superoxide production (IC₅₀ = 1.7 μM) without altering normal bioenergetics function.¹ It reduces production of H₂O₂ during oxidation of glutamate plus malate to 25% of the control level. S3QEL-2 also protects against ROS-induced cell stress in pancreatic β-cells and decreases HIF-1α induction in response to hypoxia.

Reference

1. Orr, A.L., Vargas, L., Turk, C.N., *et al.* Suppressors of superoxide production from mitochondrial complex III. *Nat. Chem. Biol.* **11**, 834-840 (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, 11/21/2022

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