

PRODUCT INFORMATION



Pyr3

Item No. 16888

CAS Registry No.: 1160514-60-2
Formal Name: 1-[4-[(2,3,3-trichloro-1-oxo-2-propen-1-yl)amino]phenyl]-5-(trifluoromethyl)-1H-pyrazole-4-carboxylic acid, ethyl ester

MF: C₁₆H₁₁Cl₃F₃N₃O₃

FW: 456.6

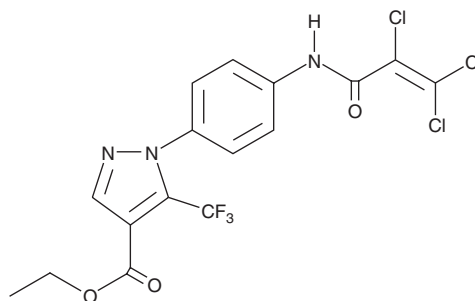
Purity: ≥98%

UV/Vis.: λ_{max}: 210, 270 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

Pyr3 is supplied as a crystalline solid. A stock solution may be made by dissolving the Pyr3 in the solvent of choice, which should be purged with an inert gas. Pyr3 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Pyr3 in these solvents is approximately 10, 20, and 50 mg/ml, respectively.

Description

Pyr3 is a pyrazole compound that selectively antagonizes the transient receptor potential canonical channel 3 (TRPC3).¹ It inhibits TRPC3-mediated Ca²⁺ influx with an IC₅₀ value of 0.7 μM without effect on other TRPC members and suppresses activation of nuclear factor of activated T cells with an IC₅₀ value of 0.05 μM.¹ Pyr3 has been shown to suppress cardiac hypertrophy in mice at 0.1 mg/kg/day.¹

Reference

1. Kiyonaka, S., Kato, K., Nishida, M., *et al.* Selective and direct inhibition of TRPC3 channels underlies biological activities of a pyrazole compound. *Proc. Natl. Acad. Sci. USA* **106**(13), 5400-5405 (2009).

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, 11/29/2022

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