

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: $C_{16}H_{11}Cl_3F_3N_3O_3$

FW: 456.6

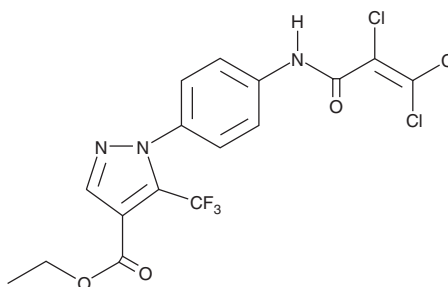
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 210, 270 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

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

Pyr3 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Pyr3 is a pyrazole compound that selectively antagonizes the transient receptor potential canonical channel 3 (TRPC3).¹ It inhibits TRPC3-mediated Ca^{2+} influx with an IC_{50} value of 0.7 μM without effect on other TRPC members and suppresses activation of nuclear factor of activated T cells with an IC_{50} 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, 04/25/2016

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