

PRODUCT INFORMATION



DUN41875

Item No. 37151

CAS Registry No.: 1295541-87-5
Formal Name: 5-methyl-N-(1-methyl-1H-pyrazol-3-yl)-1-[[3-(trifluoromethyl)phenyl]methyl]-1H-imidazole-4-carboxamide

Synonyms: SCD1 Inhibitor, Stearyl-CoA Desaturase 1 Inhibitor, Stearoyl-Coenzyme A Desaturase 1 Inhibitor

MF: C₁₇H₁₆F₃N₅O

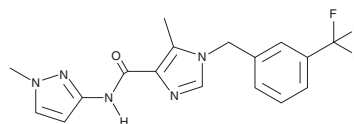
FW: 363.3

Purity: ≥98%

Supplied as: A 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

DUN41875 is supplied as a solid. A stock solution may be made by dissolving the DUN41875 in the solvent of choice, which should be purged with an inert gas. DUN41875 is soluble in methanol.

Description

DUN41875 is an inhibitor of stearyl-CoA desaturase 1 (SCD1; IC₅₀ = 11.8 nM for the rat enzyme).¹ It increases tetramerization of soluble α-synuclein and reduces mutant α-synuclein incorporation into tetramers in BE(2)-M17 neuroblastoma cells expressing a mutant α-synuclein containing three glutamate-to-lysine substitutions at positions 35, 46, and 61 (3K) when used at concentrations of 0.01, 0.1, and 1 μM.² DUN41875 (10 or 30 mg/kg) reduces the plasma ratio of oleic acid to stearic acid in rats fed a low essential fatty acid diet.¹ It reduces the time to turn and the number of falls in the pole test, increases time to fall from the rotarod, and improves endurance in the hanging wire test in the 3K-3798 transgenic mouse model of Parkinson's disease when administered in the drinking water at a concentration of 1.9% (v/v).² DUN41875 reduces cortex somatic aggregates of phosphorylated α-synuclein and cortex levels of lysosome-associated membrane protein 1 (LAMP-1), also known as CD107a, in the brains from the same mice.

References

1. Atkinson, K.A., Beretta, E.E., Brown, J.A., *et al.* N-Benzylimidazole carboxamides as potent, orally active stearylCoA desaturase-1 inhibitors. *Bioorg. Med. Chem. Lett.* **21(6)**, 1621-1625 (2011).
2. Nuber, S., Nam, A.Y., Rajsombath, M.M., *et al.* A stearyl-coenzyme A desaturase inhibitor prevents multiple Parkinson disease phenotypes in α-synuclein mice. *Ann. Neurol.* **89(1)**, 74-90 (2021).

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, 11/06/2023

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