

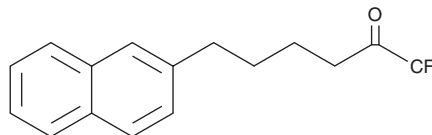
PRODUCT INFORMATION



FKGK 18

Item No. 13943

CAS Registry No.: 1071001-09-6
Formal Name: 1,1,1-trifluoro-6-(2-naphthalenyl)-2-hexanone
MF: C₁₆H₁₅F₃O
FW: 280.3
Purity: ≥95%
UV/Vis.: λ_{max}: 225, 271 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

FKGK 18 is supplied as a crystalline solid. A stock solution may be made by dissolving the FKGK 18 in the solvent of choice. FKGK 18 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of FKGK 18 in these solvents is approximately 20 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of FKGK 18 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of FKGK 18 in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

FKGK 18 is an inhibitor of group VIA (GVIA) calcium-independent phospholipase A₂ (iPLA₂).¹ It inhibits GVIA iPLA₂ by 99.9% at 0.091 mole fraction in a mixed micelle activity assay and is selective for GVIA iPLA₂ over GIVA cPLA₂ and GV sPLA₂ where it shows 80.8 and 36.8% inhibition, respectively. FKGK 18 inhibits iPLA₂β activity in cytosolic extracts from INS-1 cells overexpressing iPLA₂β (IC₅₀ = ~50 nM) as well as iPLA₂γ activity in mouse heart membrane fractions (IC₅₀s = ~1-3 μM).² It inhibits glucose-induced increases in prostaglandin E₂ (PGE₂; Item No. 14010) production and insulin secretion in human pancreatic islets when used at a concentration of 10 μM and inhibits thapsigargin-induced apoptosis in INS-1 cells overexpressing iPLA₂β in a concentration-dependent manner. FKGK 18 (20 mg/kg, 3 times per week) reduces blood glucose levels in an intraperitoneal glucose tolerance test, decreases the incidence of diabetes, and increases serum insulin levels in non-obese diabetic (NOD) mice.³

References

1. Kokotos, G., Hsu, Y.H., Burke, J.E., *et al.* Potent and selective fluoroketone inhibitors of group VIA calcium-independent phospholipase A₂. *J. Med. Chem.* **53**(9), 3602-3610 (2010).
2. Ali, T., Kokotos, G., Magrioti, V., *et al.* Characterization of FKGK18 as inhibitor of group VIA Ca²⁺-independent phospholipase A₂ (iPLA₂β): Candidate drug for preventing β-cell apoptosis and diabetes. *PLoS One* **8**(8):e71748, (2013).
3. Bone, R.N., Gai, Y., Magrioti, V., *et al.* Inhibition of Ca²⁺-independent phospholipase A₂β (iPLA₂β) ameliorates islet infiltration and incidence of diabetes in NOD mice. *Diabetes* **64**(2), 541-554 (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, 05/20/2019

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