

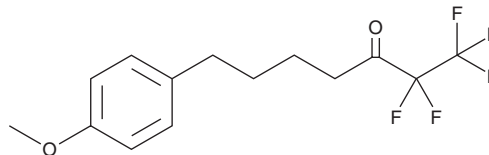
PRODUCT INFORMATION



GK187

Item No. 25130

CAS Registry No.: 1071001-50-7
Formal Name: 1,1,1,2,2-pentafluoro-7-(4-methoxyphenyl)-3-heptanone
MF: $C_{14}H_{15}F_5O_2$
FW: 310.3
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 279 nm
Supplied as: A solution in methyl acetate
Storage: $-20^{\circ}C$
Stability: ≥ 1 year



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

Laboratory Procedures

GK187 is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of GK187 in these solvents is approximately 10 mg/ml.

For maximum solubility in aqueous buffers, evaporate the methyl acetate and dissolve in ethanol. The ethanolic solution of GK187 should be diluted with the aqueous buffer of choice. GK187 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method.

Description

GK187 is an inhibitor of Group VIA (GVIA) calcium-independent phospholipase A_2 (iPLA $_2$), an enzyme involved in basal cell metabolism.¹ It inhibits GVIA iPLA $_2$ by 99.8% at 0.091 mole fraction in a mixed micelle activity assay and is selective for GVIA iPLA $_2$ over GIVA cPLA $_2$ and GV sPLA $_2$ where it shows less than 25% and 32.8% inhibition, respectively.

Reference

1. Magrioti, V., Nikolaou, A., Smyrniotou, A., *et al.* New potent and selective polyfluoroalkyl ketone inhibitors of GVIA calcium-independent phospholipase A2. *Bioorg. Med. Chem.* **21(18)**, 5823-5829 (2013).

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, 05/22/2019

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