

PRODUCT INFORMATION



Hentriacontane

Item No. 29776

CAS Registry No.: 630-04-6

Synonym: Untriacontane

MF: $C_{31}H_{64}$

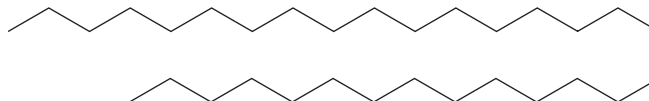
FW: 436.8

Purity: $\geq 95\%$

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

Hentriacontane is supplied as a crystalline solid. A stock solution may be made by dissolving the hentriacontane in the solvent of choice, which should be purged with an inert gas. Hentriacontane is soluble in the organic solvent heptane at a concentration of approximately 5 mg/ml.

Description

Hentriacontane is an alkane that has anti-inflammatory activity.¹ It reduces LPS-induced TNF- α , IL-6, and IL-1 β production in RAW 264.7 cells. Hentriacontane (2 and 5 mg/kg) reduces LPS-induced TNF- α , IL-6, and IL-1 β production in whole blood and carrageenan-induced paw edema in mice.

Reference

1. Khajuria, V., Gupta, S., Sharma, N., *et al.* Anti-inflammatory potential of hentriacontane in LPS stimulated RAW 264.7 cells and mice model. *Biomed. Pharmacother.* **92**, 175-186 (2017).

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, 12/08/2022

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