

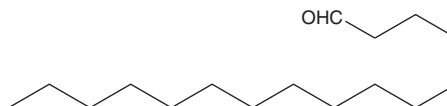
# PRODUCT INFORMATION



## Hexadecanal

Item No. 9001996

**CAS Registry No.:** 629-80-1  
**Formal Name:** hexadecanal  
**Synonyms:** 1-Hexadecanal, Palmitaldehyde  
**MF:** C<sub>16</sub>H<sub>32</sub>O  
**FW:** 240.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 207, 293 nm  
**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

Hexadecanal is supplied as a crystalline solid. A stock solution may be made by dissolving the hexadecanal in the solvent of choice, which should be purged with an inert gas. Hexadecanal is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of hexadecanal in ethanol and DMF is approximately 30 mg/ml and approximately 10 mg/ml in DMSO.

Hexadecanal is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, hexadecanal should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Hexadecanal has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Hexadecanal is the 16-carbon free fatty aldehyde analog of palmitic acid that, in conjunction with NAD<sup>+</sup>, acts as a substrate for hexadecanal:NAD<sup>+</sup> oxidoreductase (fatty aldehyde dehydrogenase).<sup>1</sup>

### Reference

1. Rizzo, W.B. and Craft, D.A. Sjögren-Larsson syndrome. Deficient activity of the fatty aldehyde dehydrogenase component of fatty alcohol:NAD<sup>+</sup> oxidoreductase in cultured fibroblasts. *J. Clin. Invest.* **88**(5), 1643-1648 (1991).

#### 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, 12/13/2022

#### 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