

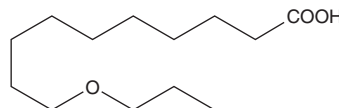
# PRODUCT INFORMATION



## O-11

Item No. 90510

**CAS Registry No.:** 119290-12-9  
**Formal Name:** 10-propoxy-decanoic acid  
**Synonyms:** FA 13:0;O, 11-Oxatetradecanoic Acid  
**MF:** C<sub>13</sub>H<sub>26</sub>O<sub>3</sub>  
**FW:** 230.3  
**Purity:** ≥98%  
**Supplied as:** A solution in methyl acetate  
**Storage:** -20°C  
**Stability:** ≥1 years



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

### Laboratory Procedures

O-11 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 (DMF) purged with an inert gas can be used. The solubility of O-11 in ethanol and DMF is approximately 30 mg/ml and approximately 20 mg/ml in DMSO.

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. If an organic solvent-free solution of O-11 is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of O-11 in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

O-11 is an analog of the fully saturated, 14-carbon fatty acid myristic acid, in which the methylene group at position 11 is replaced with oxygen. It is highly effective and selective at killing *Trypanosoma brucei*, the protozoan parasite responsible for African sleeping sickness, exhibiting an LD<sub>50</sub> of less than 1 μM in a cell culture assay.<sup>1,2</sup> The toxic effects of O-11 appear to be caused by its ability to inhibit the incorporation of a single myristate into the GPI anchor of the variant surface glycoprotein (VSG), a protein critical for evading the host immune response.<sup>1</sup> O-11 exhibits essentially no anti-fungal activity when assayed using *C. neoformans*, but does have a minor inhibitory effect on HIV-1 replication in T-lymphocytes.<sup>3</sup>

### References

1. Doering, T.L., Raper, J., Buxbaum, L.U., *et al.* An analog of myristic acid with selective toxicity for African trypanosomes. *Science* **252**, 1851-1854 (1991).
2. Doering, T.L., Lu, T., Werbovets, K.A. *et al.* Toxicity of myristic acid analogs toward African trypanosomes. *Proc. Natl. Acad. Sci. USA* **91**, 9735-9739 (1994).
3. Langner, C.A., Lodge, J.K., Travis, S.J., *et al.* 4-Oxatetradecanoic acid is fungicidal for *Cryptococcus neoformans* and inhibits replication of human immunodeficiency virus I. *J. Biol. Chem.* **267**, 17159-17169 (1992).

#### 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, 02/13/2024

#### 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