

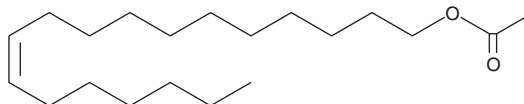
# PRODUCT INFORMATION



## 11(Z)-Vaccenyl Acetate

Item No. 10010101

**CAS Registry No.:** 6186-98-7  
**Formal Name:** (11Z)-11-octadecen-1-ol acetate  
**Synonym:** 11-cis Vaccenyl Acetate  
**MF:** C<sub>20</sub>H<sub>38</sub>O<sub>2</sub>  
**FW:** 310.5  
**Purity:** ≥98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

11(Z)-Vaccenyl acetate is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 11(Z)-vaccenyl acetate 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. The solubility of 11(Z)-vaccenyl acetate in PBS, pH 7.2, is approximately 2.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

The mating and social behaviors of insects are largely orchestrated by a suite of volatile hydrocarbon pheromones. 11(Z)-Vaccenyl Acetate is the male-specific mating pheromone of the fruit fly *D. melanogaster*. It acts selectively through the Or67d odorant receptor to control mating behavior in both male and female fruit flies.<sup>1</sup>

### Reference

1. Kurtovic, A., Widmer, A., and Dicksom, B.J. A single class of olfactory neurons mediates behavioural responses to a *Drosophila* sex pheromone. *Nature Letters* **446**, 542-546 (2007).

#### 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, 04/02/2024

#### 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