



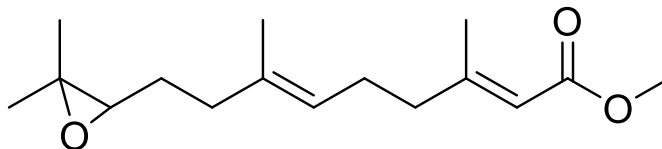
Echelon Biosciences Inc.  
675 Arapeen Drive, Suite 302  
Salt Lake City, UT 84108  
Telephone 866-588-0455  
Fax 801-588-0497  
echelon@echelon-inc.com  
www.echelon-inc.com

## Technical Data Sheet

**For research use only**  
Not intended or approved for  
diagnostic or therapeutic use.

**Product Name:** Juvenile Hormone III (*trans-trans*-10,11-Epoxyfarnesenic Acid methyl ester)

**Product Number:** S-0155



**Molecular Formula:** C<sub>16</sub>H<sub>26</sub>O<sub>3</sub>

**MW:** 266.38

**CAS:** 24198-95-6

**Solubility:** Chloroform, ethyl acetate, ethanol, most organic solvents, >1 mg/mL

**Storage:** -20 °C or below

**Field of Interest:** Juvenile Hormone III (JH III) is a sesquiterpanoid found in insects which regulates important physiological processes such as reproduction and metamorphosis. JH III is biosynthesized via the mevalonate pathway and is metabolized by JH esterase and JH epoxide hydrolase terminating the activity of the hormone. JH III has also been found in certain plants such as *Cyperus iria* and *C. Aromaticus* potentially acting as a protectant against insect herbivory.

**References:** 1) C. Helvig, *et al.* "CYP15A1, the cytochrome P450 that catalyzes epoxidation of methyl farnesoate to juvenile hormone III in cockroach corpora allata" *Proc. Nat. Sci. Acad.* 2004, 101(12), 4024–4029, doi: 10.1073/pnas.0306980101. 2) J.A. Tillman, S.J. Seybold, R.A. Jurenka, R.A., *et al.* "Insect pheromones - an overview of biosynthesis and endocrine regulation." *Insect Biochemistry and Molecular Biology*, 1999, 29(6), 481-514. 3) Y.C. Toong, D.A. Schooley, & F.C. Baker. "Isolation of Insect Juvenile Hormone III from a plant" *Nature*, 1988, 333, 170-171.

**Hazardous Properties and Cautions:** The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. Juvenile Hormone III is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

**Warranty and Disclaimer:** Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.