

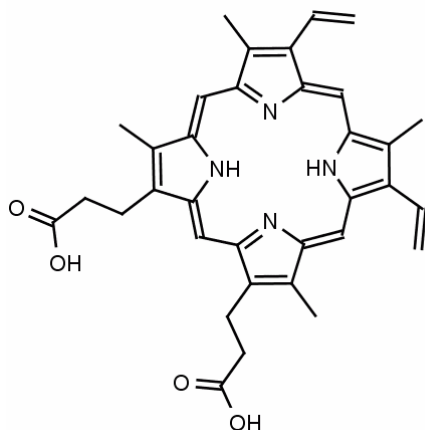


Echelon Biosciences Inc.  
675 Arapeen Drive, Suite 302  
Salt Lake City, UT 84108  
Telephone 866-588-0455  
Fax 801-588-0497  
echelon@echelon-inc.com  
www.echelon-inc.com

## Technical Data Sheet

**For research use only**

**Product Name:** Protoporphyrin IX



| Catalog Number | CAS Registry No. | MW (g/mol) | Chemical Formula  | Solubility                                   |
|----------------|------------------|------------|---|--|
| F-H050         | 533-12-8         | 562.66     | C <sub>34</sub> H <sub>34</sub> N <sub>4</sub> O <sub>4</sub> | Chloroform, glacial acetic acid or conc. HCl |

**Storage:** Protoporphyrin IX is stable for at least one year when stored as a solid, protected from moisture, at -20 °C. Protect from light.

**Field of Use:** Protoporphyrin IX is an intermediate in the Heme synthesis pathway and is formed from Protoporphyrinogen III in a reaction catalyzed by the enzyme Protoporphyrin III oxidase. Iron in the ferrous state is added to Protoporphyrin IX to form heme in a reaction catalyzed by the enzyme ferrochelatase. Protoporphyrin IX has been found to activate guanylate cyclase<sup>1</sup> and has been found to induce apoptosis in HeLa cells<sup>2</sup>.

**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.

### References:

1. Ignarro, L. J.; Wood, K. S.; Wolin, M. S. Activation of purified soluble guanylate cyclase by protoporphyrin IX. *Proc Natl Acad Sci U S A* 1982, 79, 2870-3.
2. Bednarz, N.; Zawacka-Pankau, J.; Kowalska, A. Protoporphyrin IX induces apoptosis in HeLa cells prior to photodynamic treatment. *Pharmacol Rep* 2007, 59, 474-9.

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.