

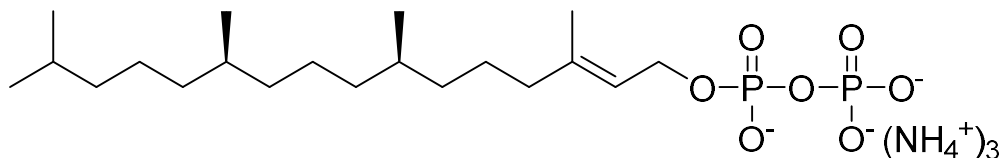
## Phytyl Diphosphate (PhytylPP)

Catalog number: I-0190

**Molecular Formula:** C<sub>20</sub>H<sub>51</sub>N<sub>3</sub>O<sub>7</sub>P<sub>2</sub>

**MW:** 507.59

**CAS:** 38476-27-6



**Alternate Name:** Phytyl pyrophosphate

**Solubility:** water (>1 mg/mL)

**Storage and Handling:** Store dry at -20 °C. Stock solutions should be stored frozen (-20 °C or below).

**Background:** Phytyl diphosphate is the precursor in the biosynthesis of tocopherol (Vitamin E) and phylloquinol in plants. It is formed by hydrolysis of the phytol side chain on chlorophyll to phytol followed by phosphorylation by the kinases VTE5 & VTE6.

**References:** 1) Ischebeck, T, et al. (2006) "A Salvage Pathway for Phytol Metabolism in *Arabidopsis*" J. Biol. Chem., 281(5), 2470-2477

2) vom Dorp, K., Hoelzl, G., Plohm, C., Eisenhut, M., Abraham, Marion; Weber, A. P. M., Hanson, A. D., Doermann, P. (2015) "Remobilization of phytol from chlorophyll degradation is essential for tocopherol synthesis and growth of *Arabidopsis*" Plant Cell, 27(10), 2846-2859.

3) Gutbrod, K., Romer, J., Doermann, P. (2019) "Phytyl Metabolism in Plants" Progress in Lipid Research, 74, 1-17

**Hazardous Properties and Cautions:** The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product 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.

