

## Glycerophosphoinositol (GroPI)

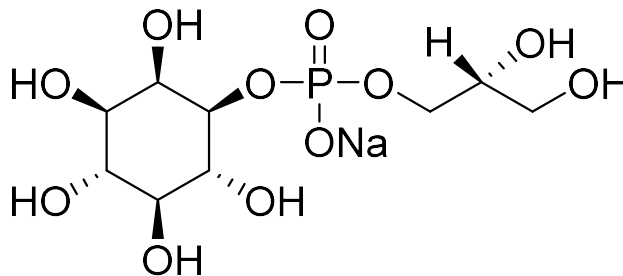
**Catalog number:** P-0000

**Molecular Formula:** C<sub>9</sub>H<sub>18</sub>NaO<sub>11</sub>P

**MW:** 356.2

**CAS#:** 129830-95-1

**Solubility:** >5mg/ml H<sub>2</sub>O



**Storage and Handling:** Glycerophosphoinositols polyphosphates (GroPIPs) are stable for at least one year when stored as a solid, protected from moisture, at -20 °C. Reconstitute with water or neutral pH, buffered salt solutions, i.e. PBS, TBS, etc. Storage in basic (pH > 9) or acidic (pH < 4) buffers may cause decomposition. After reconstitution, solutions of GroPIPs should be frozen and stored at -20 °C between uses. GroPIPs are stable for at least three months when handled in this way. Repeated freeze/thaw cycles do not affect GroPIPs. Do not store reconstituted GroPIPs at 4°C for more than 2-3 days.

**Background:** Phosphoinositides (PIPns) are minor components of cellular membranes but are integral signaling molecules for cellular communication. Phosphatidylinositol (PI or PtdIns) is biosynthesized from CDP-diacylglycerol and inositol by phosphatidylinositol synthase. PI is the building block for phosphoinositide polyphosphates through phosphorylation by PI-kinases and is a substrate for PI-PLC yielding diacylglycerol.

**References:** 1) R. Buccione, *et al.* "Glycerophosphoinositols inhibit the ability of tumour cells to invade the extracellular matrix." *Eur. J. Cancer*, **2005**, *41*, 470-476.

2) Zizza, P., C. Iurisci, *et al.* (2012). "Phospholipase A2 IValpha regulates phagocytosis independent of its enzymatic activity." *J Biol Chem* 287(20): 16849-16859.

3) Vessichelli, M., *et al.* (2017). "The natural phosphoinositide derivative glycerophosphoinositol inhibits the lipopolysaccharide-induced inflammatory and thrombotic responses." *Journal of Biological Chemistry* 292(31): 12828-12841.

**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.

Technical Data Sheet, Rev 2a, 1/7/2025 – **For research use only.** Not intended for diagnostic or therapeutic use.