# **Dodecylphosphate**

### **LPA** receptor ligand

Lysophosphatidic acid (LPA) receptor ligand. A specific agonist at  ${\rm LPA}_2$  (EC  $_{50}$  =700 nM) and antagonist at LPA

Citations: 1

View Online »

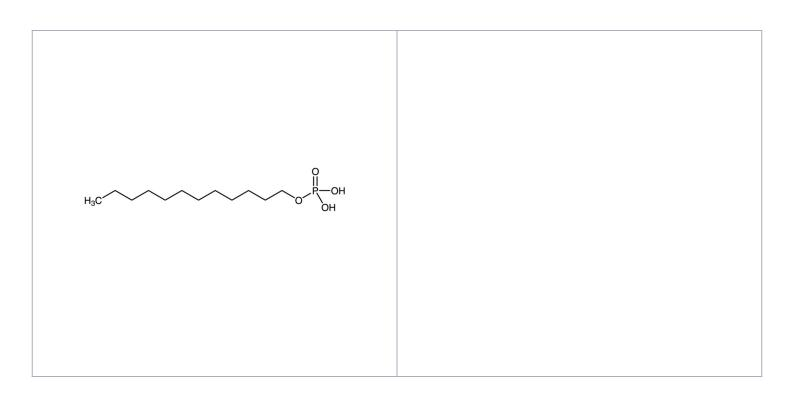
## **Ordering Information**

Order Online »

BML-LP105-0005	5mg
BML-LP105-0025	25mg

Manuals, SDS & CofA

View Online »



#### **Handling & Storage**

Use/Stability As indicated on product label or CoA when stored as recommended. Store, as supplied,

at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.

Long Term Storage -20°C

**Shipping** Ambient Temperature

### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name Dihydrogen monolauryl phosphate

**Appearance** White to off-white solid.

**CAS** 2627-35-2

Couple Target LPA receptor

Couple Type Ligand

Formula  $C_{12}H_{27}O_4P$ 

MW 266.3

Purity ≥98% (TLC)

Soluble in DMSO (15 mg/ml warm) or ethanol (2 mg/ml

warm).

**Source** Synthetic.

Last modified: May 29, 2024

