

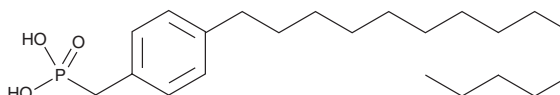
# PRODUCT INFORMATION



## BMP-22

Item No. 27264

CAS Registry No.: 1306684-90-1  
Formal Name: *p*-[[4-(pentadecylphenyl)methyl]-phosphonic acid  
MF: C<sub>22</sub>H<sub>39</sub>O<sub>3</sub>P  
FW: 382.5  
Purity: ≥95%  
UV/Vis.: λ<sub>max</sub>: 220, 265 nm  
Supplied as: A crystalline solid  
Storage: -20°C  
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

BMP-22 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMP-22 in the solvent of choice. BMP-22 is soluble in organic solvents such as ethanol and dimethyl formamide, which should be purged with an inert gas. The solubility of BMP-22 in these solvents is approximately 10 and 20 mg/ml, respectively.

### Description

BMP-22 is an inhibitor of autotaxin (IC<sub>50</sub> = 170 nM).<sup>1</sup> It is selective for autotaxin over the phosphodiesterases NPP6 and NPP7 at 10 μM. BMP-22 (0.1-1,000 nM) inhibits autotaxin-mediated production of lysophosphatidic acid (LPA) from lysophosphatidylcholine (LPC; Item No. 24331) *in vitro* in a concentration-dependent manner. It inhibits LPC-dependent MM1 cell invasion of a human umbilical vein endothelial cell (HUVEC) monolayer. BMP-22 (0.5 mg/kg per day) decreases the number of lung metastatic foci in a B16/F10 syngeneic mouse melanoma model of lung metastasis.

### Reference

1. Gupte, R., Patil, R., Liu, J., *et al.* Benzyl and naphthalene-methyl phosphonic acid inhibitors of autotaxin with anti-invasive and anti-metastatic actions. *ChemMedChem* **6**(5), 922-935 (2011).

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/30/2019

### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM