

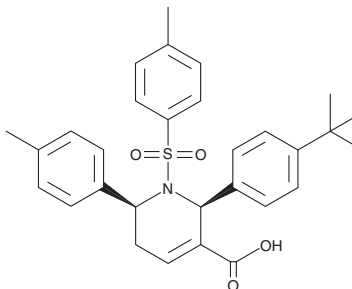
PRODUCT INFORMATION



P11

Item No. 17507

CAS Registry No.: 942285-55-4
Formal Name: (2R,6R)-rel-2-[4-(1,1-dimethylethyl)phenyl]-1,2,5,6-tetrahydro-6-(4-methylphenyl)-1-[(4-methylphenyl)sulfonyl]-3-pyridinecarboxylic acid
MF: C₃₀H₃₃NO₄S
FW: 503.7
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

P11 is supplied as a crystalline solid. A stock solution may be made by dissolving the P11 in the solvent of choice. P11 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of P11 in these solvents is approximately 30 mg/ml.

P11 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, P11 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. P11 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Lipoprotein-associated phospholipase A₂, also known as platelet-activating factor acetylhydrolase (PAF-AH) or phospholipase A₂ group 7, hydrolyzes glycerophospholipids to produce lyso-PAF/lyso-phosphatidylcholine and short and/or oxidized fatty acids, many of which have pro-inflammatory or pro-oxidative activities. PAF-AH1b2 and 1b3 form a complex with 1b1 to regulate brain development, spermatogenesis, and cancer pathogenesis. P11 is an inhibitor of PAF-AH1b2- and 1b3-mediated hydrolysis of PAF to lyso-PAF (IC₅₀s = 37 and 880 nM, respectively) without significant effect on other brain serine hydrolases.¹ At 10 μM, it has been shown to impair Neuro2a and PC3 tumor cell survival.¹

Reference

1. Chang, J.W., Zuhl, A.M., Speers, A.E., *et al.* Selective inhibitor of platelet-activating factor acetylhydrolases 1b2 and 1b3 that impairs cancer cell survival. *ACS Chem. Biol.* A-H (2015).

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, 11/17/2022

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