

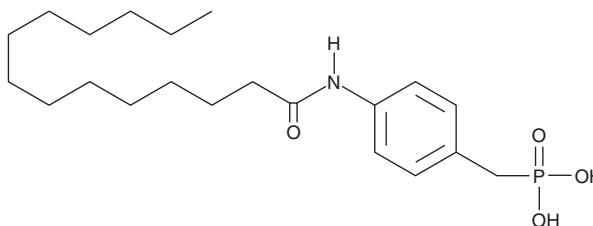
PRODUCT INFORMATION



S32826

Item No. 13664

CAS Registry No.: 1096770-84-1
Formal Name: P-[[4-[(1-oxotetradecyl)amino]phenyl]methyl]-phosphonic acid
MF: C₂₁H₃₆NO₄P
FW: 397.5
Purity: ≥98%
UV/Vis.: λ_{max}: 250 nm
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

S32826 is supplied as a crystalline solid. A stock solution may be made by dissolving the S32826 in the solvent of choice, which should be purged with an inert gas. S32826 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of S32826 in these solvents is approximately 2 and 5 mg/ml, respectively.

Description

Autotaxin is a lysophospholipase D, cleaving lysophosphatidylcholine to produce choline and lysophosphatidic acid (LPA). S32826 is a potent and selective inhibitor of autotaxin, inhibiting recombinant autotaxin b with an IC₅₀ value of 8.8 nM.¹ With similar effectiveness against the a and g autotaxin isoforms. It has no effect on a variety of receptors and enzymes, with the exceptions of src (IC₅₀ = 6 μM) and phosphatase 1B (IC₅₀ = 15 μM).¹ S32826 inhibits LPA release from adipocytes (IC₅₀ = 90 nM) and reduces plasma autotaxin activity by 57% when used at 5 μM.¹

Reference

1. Ferry, G., Moulharat, N., Pradère, J.-P., *et al.* S32826, a nanomolar inhibitor of autotaxin: Discovery, synthesis and applications as a pharmacological tool. *J. Pharmacol. Exp. Ther.* **327**, 809-819 (2008).

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, 11/14/2022

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