

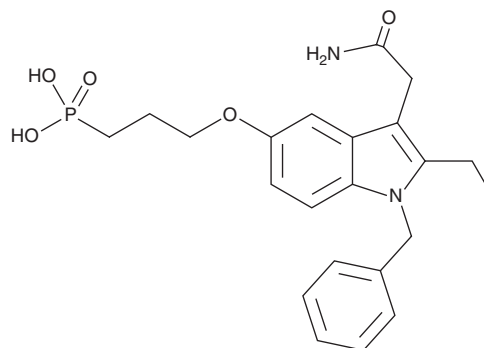
PRODUCT INFORMATION



LY311727

Item No. 17973

CAS Registry No.: 164083-84-5
Formal Name: P-[3-[[3-(2-amino-2-oxoethyl)-2-ethyl-1-(phenylmethyl)-1H-indol-5-yl]oxy]propyl]-phosphonic acid
MF: C₂₂H₂₇N₂O₅P
FW: 430.4
Purity: ≥95%
UV/Vis.: λ_{max}: 225, 283 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

LY311727 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY311727 in the solvent of choice. LY311727 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of LY311727 in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of LY311727 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of LY311727 in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Secreted phospholipase A₂ (sPLA₂) isoforms are low molecular weight, millimolar calcium-dependent enzymes that hydrolyze the fatty acid from the sn-2 position of membrane phospholipids.¹ Group IIA sPLA₂, also known as sPLA₂IIA and non-pancreatic sPLA₂, has roles in the regulation of eicosanoid synthesis and may modulate inflammatory signaling.^{1,2} LY311727 is an inhibitor of Group IIA sPLA₂ (IC₅₀ = 0.47 μM) that interacts with the active site of the enzyme in a non-covalent manner.^{2,3} It shows greater than 1,500-fold selectivity over pancreatic sPLA₂ (Group IB sPLA₂).³ LY311727 is commonly used to distinguish the actions of Group IIA sPLA₂ from those of other sPLA₂ isoforms in biological systems.⁴⁻⁷

References

1. Burke, J.E. and Dennis, E.A. *J. Lipid Res.* **50**, S237-S242 (2009).
2. Dong, C.-Z., Romieu, A., Heymans, F., et al. *Biochem. J.* **365**, 505-511 (2002).
3. Schevitz, R.W., Bach, N.J., Carlson, D.G., et al. *Nat. Struct. Biol.* **2(6)**, 458-465 (1995).
4. Hurt-Camejo, E., Andersen, S., Standal, R., et al. *Arterioscler. Thromb. Vasc. Biol.* **17**, 300-309 (1997).
5. Thommesen, L., Sjursen, W., Gåsvik, K., et al. *J. Immunol.* **161**, 3421-3430 (1998).
6. Shinohara, H., Balboa, M.A., Johnson, C.A., et al. *J. Biol. Chem.* **274(18)**, 12263-12268 (1999).
7. Balsinde, J., Shinohara, H., Lefkowitz, L.J., et al. *J. Biol. Chem.* **274**, 25967-25970 (1999).

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, 11/21/2022

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