

# PRODUCT INFORMATION

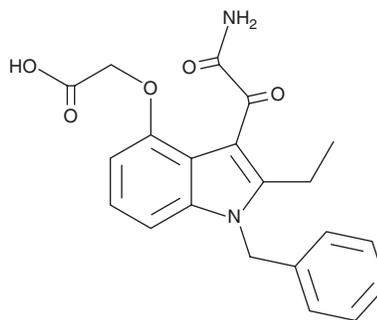


**LY315920**

Item No. 18267

**CAS Registry No.:** 172732-68-2  
**Formal Name:** 2-[[3-(2-amino-2-oxoacetyl)-2-ethyl-1-(phenylmethyl)-1H-indol-4-yl]oxy]-acetic acid

**Synonym:** Varespladib  
**MF:** C<sub>21</sub>H<sub>20</sub>N<sub>2</sub>O<sub>5</sub>  
**FW:** 380.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 217, 256, 331 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

LY315920 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY315920 in the solvent of choice, which should be purged with an inert gas. LY315920 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LY315920 in these solvents is approximately 3 and 2 mg/ml, respectively.

LY315920 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LY315920 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. LY315920 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

LY315920 is an inhibitor of human group IIA (hGIIA) secretory phospholipase A<sub>2</sub> (sPLA<sub>2</sub>; IC<sub>50</sub> = 9 nM).<sup>1</sup> It also potently inhibits human GII E and GX isoforms, as well as the corresponding mouse isoforms, but poorly inhibits other PLA<sub>2</sub> enzymes.<sup>1-4</sup> LY315920 inhibits sPLA<sub>2</sub>-induced release of thromboxane A<sub>2</sub> from isolated guinea pig lung bronchoalveolar lavage cells (IC<sub>50</sub> = 0.79 μM).<sup>2</sup> Intravenous or oral administration of LY315920 to transgenic mice expressing human sPLA<sub>2</sub> inhibits serum sPLA<sub>2</sub> activity in a dose-related manner over four hours.<sup>2</sup>

## References

1. Draheim, S.E., Bach, N.J., Dillard, R.D., *et al.* Indole inhibitors of human nonpancreatic secretory phospholipase A<sub>2</sub>. 3. Indole-3-glyoxamides. *J. Med. Chem.* **39(26)**, 5159-5175 (1996).
2. Snyder, D.W., Bach, N.J., Dillard, R.D., *et al.* Pharmacology of LY315920/S-5920, [[3-(aminooxoacetyl)-2-ethyl-1-(phenylmethyl)-1H-indol-4-yl]oxy] acetate, a potent and selective secretory phospholipase A<sub>2</sub> inhibitor: A new class of anti-inflammatory drugs, SPI. *J. Pharmacol. Exp. Ther.* **288(3)**, 1117-1124 (1999).
3. Oslund, R.C., Cermak, N., and Gelb, M.H. Highly specific and broadly potent inhibitors of mammalian secreted phospholipases A<sub>2</sub>. *J. Med. Chem.* **51(15)**, 4708-4714 (2008).
4. Smart, B.P., Pan, Y.H., Weeks, A.K., *et al.* Inhibition of the complete set of mammalian secreted phospholipases A<sub>2</sub> by indole analogues: A structure-guided study. *Bioorg. Med. Chem.* **12(7)**, 1737-1749 (2004).

### 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, 10/06/2022

## 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