

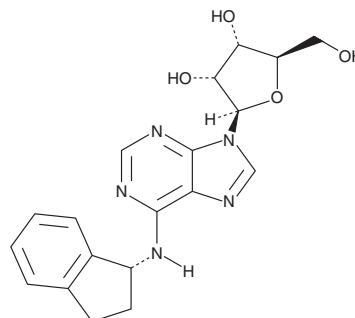
# PRODUCT INFORMATION



**PD 117519**

Item No. 28444

**CAS Registry No.:** 96392-15-3  
**Formal Name:** N-[(1R)-2,3-dihydro-1H-inden-1-yl]-adenosine  
**Synonym:** CI-947  
**MF:** C<sub>19</sub>H<sub>21</sub>N<sub>5</sub>O<sub>4</sub>  
**FW:** 383.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 211, 273 nm  
**Supplied as:** A 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

PD 117519 is supplied as a solid. A stock solution may be made by dissolving the PD 117519 in the solvent of choice, which should be purged with an inert gas. PD 117519 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PD 117519 in these solvents is approximately 0.1, 5, and 10 mg/ml, respectively.

PD 117519 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PD 117519 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PD 117519 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

PD 117519 is an adenosine receptor agonist.<sup>1</sup> It selectively binds to adenosine A<sub>2</sub> receptors (IC<sub>50</sub> = 30 μM) over A<sub>1</sub> receptors (IC<sub>50</sub> = 810 μM) in rat brain membranes. PD 117519 increases heart rate and decreases systolic blood pressure in normotensive dogs when administered at doses of 2 or 10 mg/kg.<sup>2</sup>

## References

1. Steffen, R.P. Method of treating heart failure and medicaments therefor. *Warner-Lambert Company*. **4,693,996** (1987).
2. Enerson, B.E., Lin, A., Lu, B., *et al.* Acute drug-induced vascular injury in beagle dogs: Pathology and correlating genomic expression. *Toxicol. Pathol.* **34**(1), 27-32 (2006).

### 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, 12/05/2022

## 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](http://WWW.CAYMANCHEM.COM)