

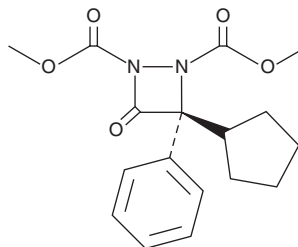
# PRODUCT INFORMATION



## ABL 127

Item No. 10833

**CAS Registry No.:** 1073529-41-5  
**Formal Name:** (3R)-3-cyclopentyl-4-oxo-3-phenyl-1,2-diazetidone-1,2-dicarboxylic acid, 1,2-dimethyl ester  
**MF:** C<sub>17</sub>H<sub>20</sub>N<sub>2</sub>O<sub>5</sub>  
**FW:** 332.4  
**Purity:** ≥98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

ABL 127 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO purged with an inert gas can be used. The solubility of ABL 127 in DMSO is approximately 5 mg/ml.

### Description

ABL 127 is a covalent aza-β-lactam inhibitor of protein phosphatase methylesterase-1 (PME-1; IC<sub>50</sub> = 4.2 nM).<sup>1</sup> It selectively inhibits PME-1 over 50 other serine hydrolases in MDA-MB-231 and HEK293T cell lysates at 10 μM. ABL 127 decreases proliferation of Ishikawa cells and migration of ECC-1 endometrial carcinoma cells when used at a concentration of 50 nM.<sup>2</sup>

### References

1. Bachovchin, D.A., Mohr, J.T., Speers, A.E., *et al.* Academic cross-fertilization by public screening yield a remarkable class of protein phosphatase methylesterase-1 inhibitors. *Proc. Natl. Acad. Sci. U.S.A.* **108**(17), 6811-6 (2011).
2. Pusey, M., Bail, S., Xu, Y., *et al.* Inhibition of protein methylesterase 1 decreased cancerous phenotypes in endometrial adenocarcinoma cell lines and xenograft tumor models. *Tumour Biol.* **37**(9), 11835-11842 (2016).

#### 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, 11/08/2022

#### 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)