

# PRODUCT INFORMATION



## DPP-9 Inhibitor 42

Item No. 39881

**Formal Name:** tert-butyl ((7S)-3r-((2-(5R,6-difluoroisoindolin-2-yl)-2-oxoethyl)amino)adamantan-1s-yl)carbamate

**Synonym:** Dipeptidyl Peptidase 9 Inhibitor 42

**MF:** C<sub>25</sub>H<sub>33</sub>F<sub>2</sub>N<sub>3</sub>O<sub>3</sub>

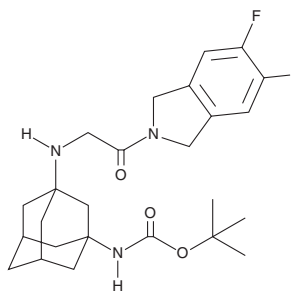
**FW:** 461.7

**Purity:** ≥95%

**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

DPP-9 inhibitor 42 is supplied as a solid. A stock solution may be made by dissolving the DPP-9 inhibitor 42 in the solvent of choice, which should be purged with an inert gas. DPP-9 inhibitor 42 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

### Description

DPP-9 inhibitor 42 is an inhibitor of dipeptidyl peptidase 9 (DDP-9; IC<sub>50</sub> = 0.0034 μM).<sup>1</sup> It is selective for DDP-9 over DPP-8, DPP-4, DPP-2, fibroblast activation protein (FAP), and prolyl endopeptidase (IC<sub>50</sub>s = 0.6, >10, >5, 10, and >10 μM, respectively). DPP-9 inhibitor 42 (10 μM) induces lactate dehydrogenase (LDH) release from THP-1 acute monocytic and HL-60 promyeloblast leukemia cells.

### Reference

1. Benramdane, S., De Loose, J., Filippi, N., *et al.* Highly selective inhibitors of dipeptidyl peptidase 9 (DPP9) derived from the clinically used DPP4-inhibitor vildagliptin. *J. Med. Chem.* **66(18)**, 12717-12738 (2023).

#### 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, 05/16/2024

#### 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