

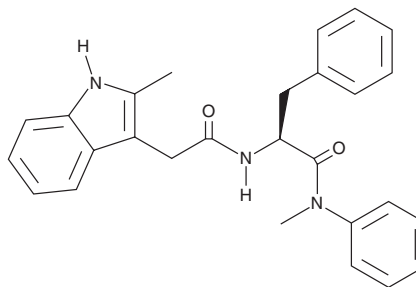
# PRODUCT INFORMATION



**PF-3450074**

Item No. 36107

**CAS Registry No.:** 1352879-65-2  
**Formal Name:** 2-methyl-N-[(1S)-2-(methylphenylamino)-2-oxo-1-(phenylmethyl)ethyl]-1H-indole-3-acetamide  
**Synonym:** PF74  
**MF:** C<sub>27</sub>H<sub>27</sub>N<sub>3</sub>O<sub>2</sub>  
**FW:** 425.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 220 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

PF-3450074 (PF74) is supplied as a solid. A stock solution may be made by dissolving the PF74 in the solvent of choice, which should be purged with an inert gas. PF74 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PF74 in these solvents is approximately 2 and 1 mg/ml, respectively.

## Description

PF74 is an inhibitor of HIV-1 capsid protein.<sup>1</sup> It reduces viral infectivity in HeLa cells infected with wild-type HIV-1 but not P38A, E45A, Q63A/Q67A, or Q219A capsid protein mutant HIV-1, simian immunodeficiency virus, or murine leukemia virus when used at a concentration of 10 μM. PF74 (0.62-10 μM) reduces capsid protein levels in HIV-1 viral cores.

## Reference

1. Shi, J., Zhou, J., Shah, V.B., *et al.* Small-molecule inhibition of human immunodeficiency virus type 1 infection by virus capsid destabilization. *J. Virol.* **85**(1), 542-549 (2011).

### 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, 12/19/2022

## 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)