

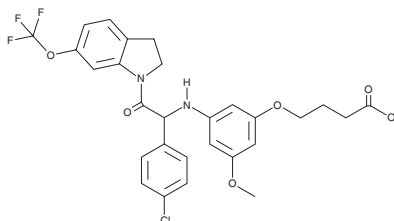
PRODUCT INFORMATION



JNJ-A07

Item No. 43714

CAS Registry No.: 2135640-93-4
Formal Name: (+)-4-[3-[[1-(4-chlorophenyl)-2-[2,3-dihydro-6-(trifluoromethoxy)-1H-indol-1-yl]-2-oxoethyl]amino]-5-methoxyphenoxy]-butanoic acid
MF: C₂₈H₂₆ClF₃N₂O₆
FW: 579.0
Purity: ≥98%
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

JNJ-A07 is supplied as a solid. A stock solution may be made by dissolving the JNJ-A07 in the solvent of choice, which should be purged with an inert gas. JNJ-A07 is sparingly soluble (1-10 mg/ml) in acetonitrile and DMSO.

Description

JNJ-A07 is an inhibitor of the protein-protein interaction between dengue virus non-structural protein 3 (NS3) and NS4B (EC₅₀ = 35 nM in a pull-down assay).¹ It inhibits replication of 21 dengue virus clinical isolates in Vero E6 cells (EC₅₀s = 0.07-0.3 nM). JNJ-A07 reduces dengue virus RNA levels in mosquitoes (*A. aegypti*) when fed blood meals containing it at a concentration of at 2 μM.² It decreases blood dengue virus RNA levels and increases survival in a mouse model of dengue virus infection when administered at doses of 1, 3, 10, or 30 mg/kg.¹

References

1. Kaptein, S.J.F., Goethals, O., Kiemel, D., *et al.* A pan-serotype dengue virus inhibitor targeting the NS3-NS4B interaction. *Nature* **597(7881)**, 504-509 (2021).
2. Rosales-Rosas, A.L., Goossens, S., Chiu, W., *et al.* The antiviral JNJ-A07 significantly reduces dengue virus transmission by *Aedes aegypti* mosquitoes when delivered via blood-feeding. *Sci. Adv.* **10(48)**, eadr8338 (2024).

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, 07/24/2025

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