

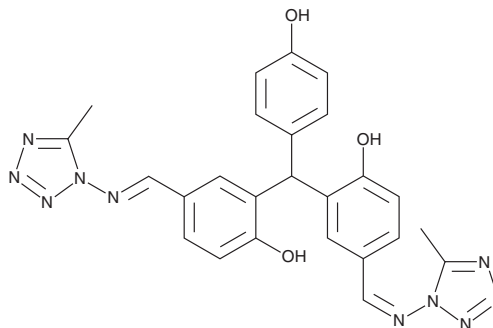
PRODUCT INFORMATION



VP-14637

Item No. 17805

CAS Registry No.: 235106-62-4
Formal Name: 2,2'-[(4-hydroxyphenyl)methylene]bis[4-[[[(5-methyl-1H-tetrazol-1-yl)imino]methyl]-phenol]
MF: C₂₅H₂₂N₁₀O₃
FW: 510.5
Purity: ≥98%
UV/Vis.: λ_{max}: 238, 326 nm
Supplied as: A crystalline 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

VP-14637 is supplied as a crystalline solid. A stock solution may be made by dissolving the VP-14637 in the solvent of choice, which should be purged with an inert gas. VP-14637 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of VP-14637 in these solvents is approximately 10 mg/ml.

VP-14637 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, VP-14637 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. VP-14637 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

VP-14637 is a small molecule inhibitor of RSV that demonstrates antiviral activity with an EC₅₀ value of 1.4 nM.^{1,2} It interferes directly with RSV fusion (EC₅₀ = 5.4 nM), binding to the viral F protein, which allows simultaneous interaction with both of the heptad repeat domains of the F protein trimer.^{1,2}

References

1. Douglas, J.L., Panis, M.L., Ho, E., *et al.* Inhibition of respiratory syncytial virus fusion by the small molecule VP-14637 via specific interactions with F protein. *J. Virol.* **77**(9), 5054-5064 (2003).
2. Douglas, J.L., Panis, M.L., Ho, E., *et al.* Small molecules VP-14637 and JNJ-2408068 inhibit respiratory syncytial virus fusion by similar mechanisms. *Antimicrob. Agents Chemother.* **49**(6), 2460-2466 (2005).

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, 11/17/2022

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