

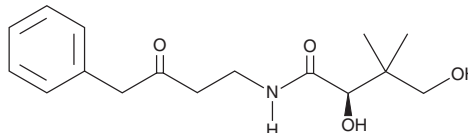
PRODUCT INFORMATION



RR6

Item No. 9001411

CAS Registry No.: 1351758-37-6
Formal Name: 2R-hydroxy-3-(hydroxymethyl)-3-methyl-N-(3-oxo-4-phenylbutyl)-butanamide
MF: C₁₆H₂₃NO₄
FW: 293.4
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

RR6 is supplied as a solid. A stock solution may be made by dissolving the RR6 in the solvent of choice, which should be purged with an inert gas. RR6 is soluble in dichloromethane and methanol.

Description

RR6 is an inhibitor of the pantetheinase vanin-1 ($IC_{50} = 0.54 \mu\text{M}$ for the recombinant human enzyme) and is a derivative of the CoA precursor D-pantothenic acid (Item No. 17288).¹ It is selective for vanin-1 over biotinidase, cathepsin B, cathepsin L, and papain (IC_{50} s = >200 μM for all). RR6 is active against *P. falciparum* ($IC_{50} = 14.5 \mu\text{M}$).² It reduces plasma levels of cholesterol and increases plasma levels of free fatty acids in fasted rats when administered in the drinking water at 3 mg/ml.¹ RR6 decreases neointima formation and macrophage infiltration in aortic allografts in a rat model of transplant vasculopathy.³

References

1. Jansen, P.A., van Diepen, J.A., Ritzen, B., *et al.* Discovery of small molecule vanin inhibitors: New tools to study metabolism and disease. *ACS Chem. Biol.* **8**(3), 530-534 (2013).
2. Pett, H.E., Jansen, P.A.M., Hermkens, P.H.H., *et al.* Novel pantothenate derivatives for anti-malarial chemotherapy. *Malar. J.* **14**, 169 (2015).
3. Wedel, J., Jansen, P.A.M., Botman, P.N.M., *et al.* Pharmacological inhibition of vanin activity attenuates transplant vasculopathy in rat aortic allografts. *Transplantation* **100**(8), 1656-1666 (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, 01/31/2024

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