

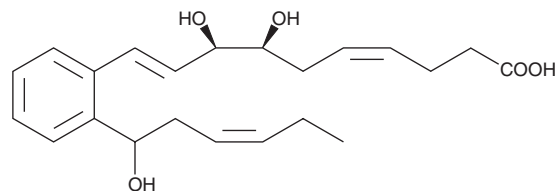
PRODUCT INFORMATION



17(R,S)-Benzo-Resolvin D1

Item No. 10623

Formal Name: (4Z,7S,8R,9E)-7,8-dihydroxy-10-(2-((Z)-1-hydroxyhex-3-en-1-yl)phenyl)deca-4,9-dienoic acid
Synonyms: Benzo-RvD1, 17(R,S)-Benzo-RvD1
MF: C₂₀H₃₀O₅
FW: 374.5
Purity: ≥95%
UV/Vis.: λ_{max}: 250 nm
Supplied as: A solution in ethanol
Storage: -80°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

17(R,S)-Benzo-resolvin D1 (benzo-RvD1) is a derivative of the specialized pro-resolving mediator (SPM) RvD1 (Item No. 10012554).¹ It reduces PDGF-BB-induced cytoskeletal changes in, and PDGF-induced migration of, isolated human vascular smooth muscle cells (VSMCs) when used at concentrations of 10 and 100 nM. Benzo-RvD1 (10 nM) inhibits p65 nuclear translocation in human umbilical vein endothelial cells (HUVECs). It also increases RAW 264.7 phagocytosis of *S. aureus* and zymosan particles.

Reference

1. Kim, A.S., Werlin, E.C., Kagaya, H., *et al.* 17R/S-Benzo-RvD1, a synthetic resolvin D1 analogue, attenuates neointimal hyperplasia in a rat model of acute vascular injury. *PLoS Biol.* **17**(2), e0264217 (2022).

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, 06/21/2023

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