

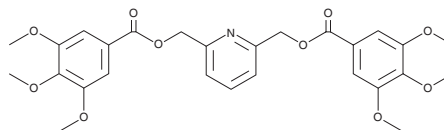
PRODUCT INFORMATION



Pirozadil

Item No. 37292

CAS Registry No.: 54110-25-7
Formal Name: 3,4,5-trimethoxy-benzoic acid, 1,1'-[2,6-pyridinediylbis(methylene)] ester
MF: C₂₇H₂₉NO₁₀
FW: 527.5
Purity: ≥98%
UV/Vis.: λ_{max}: 215 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

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

Description

Pirozadil is a lipid-lowering agent.¹ It reduces serum and aortic cholesterol levels, serum LDL levels, and atherosclerotic lesion areas on the aorta and coronary lumen in a rabbit model of atherosclerosis induced by a high-cholesterol diet when administered at a dose of 50 mg/kg per day.

Reference

1. Parwaresch, M.R., Haacke, H., and Mäder, C. Efficacy of hypolipidemic treatment in inhibition of experimental atherosclerosis: The effect of nicotinic acid and related compounds. *Atherosclerosis* **31(4)**, 395-401 (1978).

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, 10/05/2023

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