

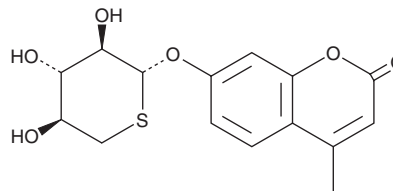
PRODUCT INFORMATION



Odiparcil

Item No. 44255

CAS Registry No.: 137215-12-4
Formal Name: 4-methyl-7-[(5-thio-β-D-xylopyranosyl)oxy]-2H-1-benzopyran-2-one
Synonym: SB-424323
MF: C₁₅H₁₆O₆S
FW: 324.3
Purity: ≥95%
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

Odiparcil is supplied as a solid. A stock solution may be made by dissolving the odiparcil in the solvent of choice, which should be purged with an inert gas. Odiparcil is slightly soluble (0.1-1 mg/ml) in DMSO and a 1:1 solution acetonitrile:water.

Description

Odiparcil is a β-D-thioxyloside derivative and driver of circulating glycosaminoglycan (GAG) synthesis.^{1,2} It increases the secretion of total sulfated GAGs in bovine aortic endothelial (BAE) cells when used at a concentration of 10 μM.¹ Odiparcil (10 μM) reduces intracellular chondroitin sulfate levels in primary skin fibroblasts isolated from patients with mucopolysaccharidosis type VI (MPS VI), an inborn error of metabolism characterized by a deficiency in the degradation of specific GAGs. Dietary administration of odiparcil (4.5 g/kg of chow) increases urinary excretion of sulfated GAGs and decreases hepatic and renal accumulation of sulfated GAGs in a mouse model of MPS VI induced by a homozygous gene mutation in the *Arsb* gene, which encodes the enzyme arylsulfatase B. Odiparcil reduces thrombus weight in a rat model of thrombosis (ED₅₀ = 3.2 mg/kg).²

References

1. Entchev, E., Jantzen, I., Masson, P., *et al.* Odiparcil, a potential glycosaminoglycans clearance therapy in mucopolysaccharidosis VI-Evidence from *in vitro* and *in vivo* models. *PLoS One* **15**(5), e0233032 (2020).
2. Toomey, J.R., Abboud, M.A., Valocik, R.E., *et al.* A comparison of the β-D-xyloside, odiparcil, to warfarin in a rat model of venous thrombosis. *J. Thromb. Haemost.* **4**(9), 1989-1996 (2006).

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, 09/10/2025

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