

PRODUCT INFORMATION



β -L-Gulopyranosyl-caldarchaetidyl-glycerol

Item No. 27203

CAS Registry No.: 416880-23-4

Formal Name: [(2S,7R,11R,15S,19S,22S,26S,30R,34R,38S,43R,47R,51S,55S,58S,62S,66R,70R)-38-[[[(2,3-dihydroxypropoxy)hydroxyphosphinyl]oxy]methyl]-7,11,15,19,22,26,30,34,43,47,51,55,58,62,66,70-hexadecamethyl-1,4,37,40-tetraoxacyclodoheptacont-2-yl)methyl β -L-gulopyranoside

Synonyms: GuCGp, Gulopyranosyl-(β 1-1)-caldarchaetidylglycerol

MF: $C_{95}H_{189}O_{16}P$

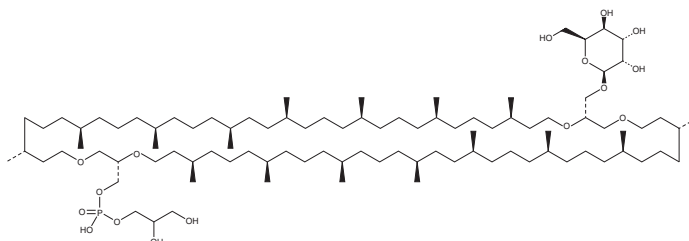
FW: 1,618.5

Purity: $\geq 95\%$

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

β -L-Gulopyranosyl-caldarchaetidyl-glycerol is supplied as a solid. A stock solution may be made by dissolving the β -L-gulopyranosyl-caldarchaetidyl-glycerol in the solvent of choice, which should be purged with an inert gas. β -L-Gulopyranosyl-caldarchaetidyl-glycerol is soluble in a 2:1 solution of chloroform:methanol.

Description

β -L-Gulopyranosyl-caldarchaetidyl-glycerol is the main polar lipid of the archaeon *T. acidophilum*.¹ Cellular levels of β -L-gulopyranosyl-caldarchaetidyl-glycerol in *T. acidophilum* HO-62 positively and negatively correlate with increases in culture pH and temperature, respectively. β -L-gulopyranosyl-caldarchaetidyl-glycerol has been used as a standard for the quantification of polar lipids extracted from *T. acidophilum* HO-62.

Reference

1. Shimada, H., Nemoto, N., Shida, Y., *et al.* Effects of pH and temperature on the composition of polar lipids in *Thermoplasma acidophilum* HO-62. *J. Bacteriol.* **190**(15), 5404-5411 (2008).

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, 11/30/2022

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