

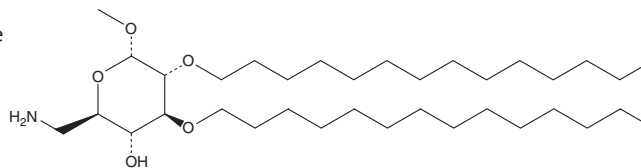
PRODUCT INFORMATION



IAXO-102

Item No. 34567

CAS Registry No.: 1115270-63-7
Formal Name: methyl 6-amino-6-deoxy-2,3-di-O-tetradecyl- α -D-glucopyranoside
MF: $C_{35}H_{71}NO_5$
FW: 585.9
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

IAXO-102 is supplied as a solid. A stock solution may be made by dissolving the IAXO-102 in the solvent of choice, which should be purged with an inert gas. IAXO-102 is soluble in methanol.

Description

IAXO-102 is a glycolipid and an antagonist of toll-like receptor 4 (TLR4).¹ It inhibits lipid A-induced TLR4 activation in a cell-based assay ($IC_{50} = 5.013 \mu\text{M}$). IAXO-102 (10 mg/kg) reduces LPS-induced lethality in mice. It decreases angiotensin II-induced aortic expansion and rupture, as well as the incidence of abdominal aortic aneurysms, in *ApoE*^{-/-} mice when administered at a dose of 3 mg/kg.²

References

1. Piazza, M., Rossini, C., Fiorentina, S.D., et al. Glycolipids and benzylammonium lipids as novel antiseptis agents: Synthesis and biological characterization. *J. Med. Chem.* **52**(4), 1209-1213 (2009).
2. Huggins, C., Pearce, S.F., Peri, F., et al. A novel small molecule TLR4 antagonist (IAXO-102) negatively regulates non-hematopoietic toll like receptor 4 signalling and inhibits aortic aneurysms development. *Atherosclerosis* **242**(2), 5630570 (2015).

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, 12/06/2022

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