

Cord Factor (endotoxin-free grade)

Immunomodulator

Immunomodulatory compound / adjuvant. Induces activation of macrophages independent of TLR2 or TLR4.

Citations: 19

[View Online »](#)

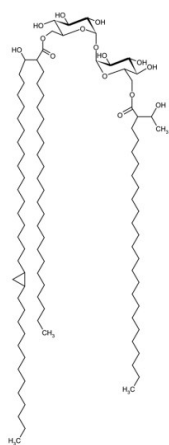
Ordering Information

[Order Online »](#)

ALX-581-210-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Trehalose 6,6'-dimycolate, TDM, 6,6'-Di-O-mycoloyl- α,α -trehalose
CAS	61512-20-7
Formulation	Lyophilized.
Purity	Single spot (TLC)
Solubility	Soluble in chloroform:methanol:water (90:10:1) (5mg/ml), hexane or isopropanol.
Source	Isolated from <i>Mycobacterium bovis</i> , Bacillus Calmette Guérin.

Last modified: November 25, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)