

Fumonisin B₂

Ceramide synthase inhibitor

Mycotoxin, structurally similar to fumonisin B₁ (Prod. No. BML-SL220). Found as contaminant mostly in corn silage; was detected recently for the first time in the industrially important *Aspergillus niger*. Induces cancer. Inducer of apoptosis. Inhibits sphinganine N-acyltransferase (ceramide synthase) at IC₅₀ = 0.1 μM. Inhibits de novo sphingolipid biosynthesis.

Citations: 8

[View Online »](#)

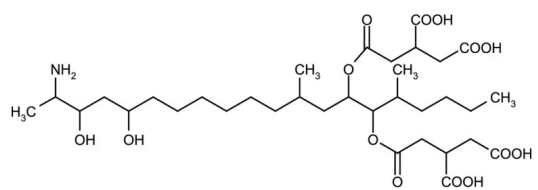
Ordering Information

[Order Online »](#)

BML-SL219-0001	1mg
BML-SL219-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Off-white solid.
CAS	116355-84-1
Couple Target	Ceramide synthase, Sphinganine N-acyltransferase
Couple Type	Inhibitor
Formula	$C_{34}H_{59}NO_{14}$
MI	14: 4289 (Fumonisin B ₁)
MW	705.8
Purity	≥98% (TLC)
RTECS	TZ8335000
Solubility	Soluble in water (25mg/ml).

Last modified: May 29, 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)