

Moniliformin . sodium salt

Mycotoxin

Mycotoxin.

Citations: 10

[View Online »](#)

Ordering Information

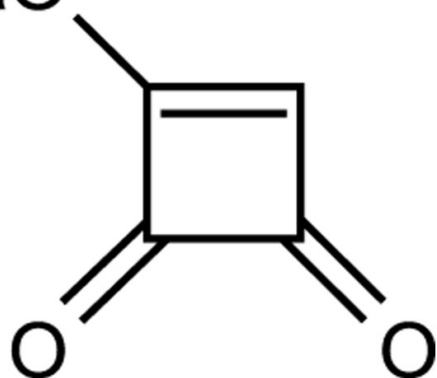
[Order Online »](#)

ALX-630-111-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

NaO



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-Hydroxycyclobut-1-ene-3,4-dione . sodium salt

Appearance Light yellow powder.

CAS 71376-34-6

Formula $C_4H_3O_3 \cdot Na$

Identity Identity determined by NMR.

MW 97.1 . 23.0

Purity ≥98% (TLC)

RTECS GU1815000

Solubility Soluble in water, methanol or DMSO.

Source Isolated from *Fusarium moniliforme*.

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-
be@enzolifesciences.com](mailto:info-be@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)