

Monosodium urate (crystals)

NLRP3 inflammasome activator

NLRP3/NALP3 inflammasome activator. Stimulates the caspase-1 activating NLRP3/NALP3 inflammasome, resulting in the production of active IL-1 β and IL-18.

Citations: 13

[View Online »](#)

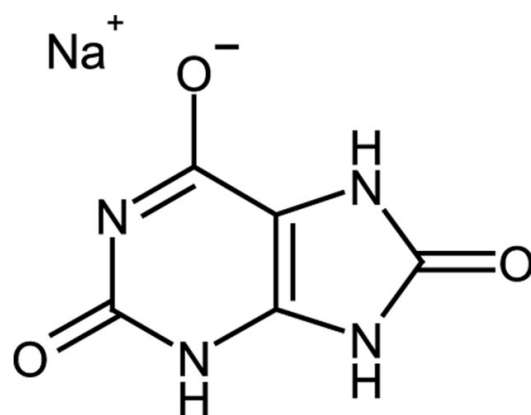
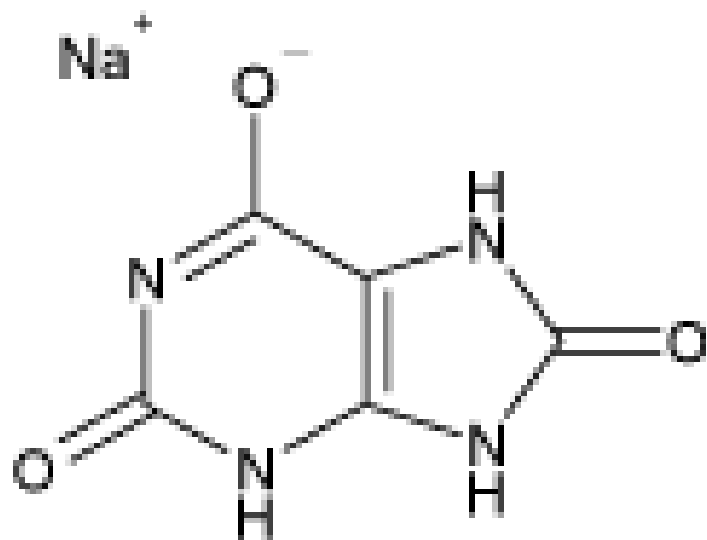
Ordering Information

[Order Online »](#)

ALX-400-047-M002	2mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Vortex before use.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	MSU (crystals), Uric Acid (crystals)
Appearance	White to off white solid.
CAS	1198-77-2
Couple Target	Inflammasome
Couple Type	Activator
Formula	$C_5H_3N_4O_3 \cdot Na$
Identity	Determined by IR.
MW	190.1
RTECS	YU7062000
Solubility	Soluble in PBS or 1M NaOH (25 mg/ml). Solution may appear as a slightly hazy suspension.

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)