

Nourseothricin – Powder (100 g)

powder (non-sterile)

Ordering Information

[Order Online »](#)

JBS-AB-102-100G	100g
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NTC

Appearance beige powder

CAS 96736-11-7

Couple Type Inhibitor

Formula $C_{19}H_{34}N_8O_8 \cdot H_2SO_4$ (Streptothricin F)

Formulation non-sterile powder

MW 600.6 (Streptothricin F)

Solubility 1000 g/l in water

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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