

T4 DNA Ligase, S pack

E. coli lambda lysogen NM 989

Ordering Information

[Order Online »](#)

JBS-EN-149S	400U
-------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Concentration 2.5 Weiss units/μl (500 CE units/μl)

Formulation liquid (Supplied in 10 mM Tris-HCl pH 7.4, 50 mM KCl, 0.1 mM EDTA, 1 mM DTT, 200 μg/ml BSA and 50 % [v/v] glycerol)

Quantity 400 Weiss units (80000 CE units). One Weiss unit is defined as the amount of enzyme required to catalyze the exchange of 1 nmol of 32P from pyrophosphate to ATP, into Norit-adsorbable material in 20 minutes at 37°C.

Source E. coli lambda lysogen NM 989

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