

LEXSY *in vitro* Translation Kit (for plasmid based cell- free protein synthesis)

The LEXSY cell extract for *in vitro* translation contains functional ribosomes and all essential components of the eukaryotic translation and folding machinery. For target mRNA generation the heterologous T7 RNA polymerase was added to the extracts. To ensure a low background the translation of endogenous host mRNAs is efficiently blocked by an antisense oligonucleotide making use of the unique gene organization of *Leishmania*.

DNA templates can be generated by cloning into a plasmid or by PCR amplification. The plasmid-based version was designed for high yield *in vitro* translation whereas the PCR-based version allows high-throughput applications.

Ordering Information

[Order Online »](#)

JBS-EGE-2002-15	30Reactions
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

For plasmid based cell-free protein synthesis

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