

LEXSY BHI – Powder Media Kit, S pack

brain-heart infusion-based; recommended for transfection and strain maintenance

Ordering Information

[Order Online »](#)

JBS-ML-412S	1Kit
-------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability Store components as indicated on individual labels.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes For the growth of LEXSY hosts P10 (Cat.-No. LT-101) and T7-TR (Cat.-No. LT-110).

Couple Type Buffer

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