

# LEXSY BHI – Powder Media Kit, XL pack

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

## Ordering Information

[Order Online »](#)

JBS-ML-412XL	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](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-be@enzolifesciences.com](mailto:info-be@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)