

# LEXSYcon2.1 Expression Kit (contains the integrative vector pLEXSY-hyg2.1)

Ordering Information

[Order Online »](#)

JBS-EGE-1310hyg	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 constitutive cytosolic or secretory protein expression

**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)