

## pLEXS<sub>Y</sub>\_I-ble2

### LEXS<sub>Y</sub>induce Expression Vector

Cat.-No.	Size
EGE-241	5 µg

#### Description:

Inducible LEXSY expression vector pLEXS<sub>Y</sub>\_I-ble2 with the *ble* marker gene (*bleomycin resistance gene* from *S. hindustans*), allowing selection of recombinant LEXSY strains with the antibiotic LEXSY Bleo (AB-103). Designed for inducible expression of target genes in LEXSY host T7-TR (LT-110).

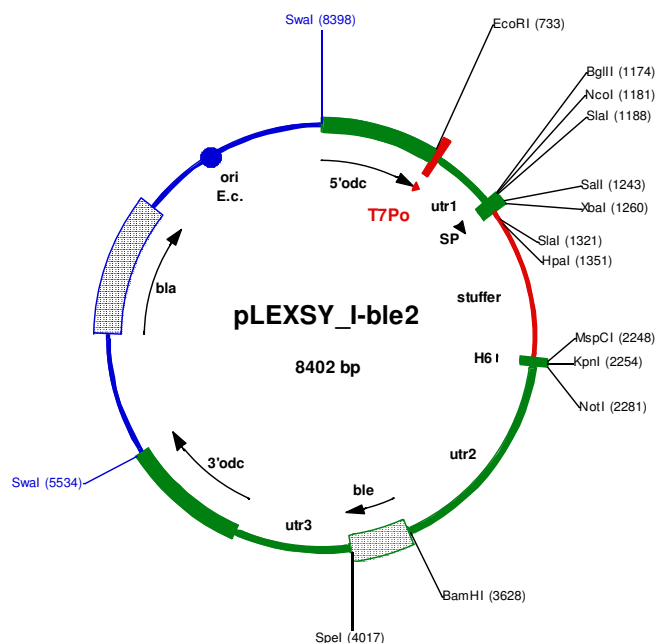
#### Storage conditions:

Store at -20°C.

Sequence information available for download from corporate WS [www.jenabioscience.com](http://www.jenabioscience.com)

**FOR RESEARCH USE ONLY.**

**NOT INTENDED FOR HUMAN OR ANIMAL  
 DIAGNOSTIC OR THERAPEUTIC USES.**



Find more information in the LEXS<sub>Y</sub>induce User's Guide (download pdf from the LEXSY section of our website).