

## pLEXY\_I-blecherry3 LEXSinduce3 Expression Vector

Cat.-No.	Size
EGE-243	5 µg

### Description:

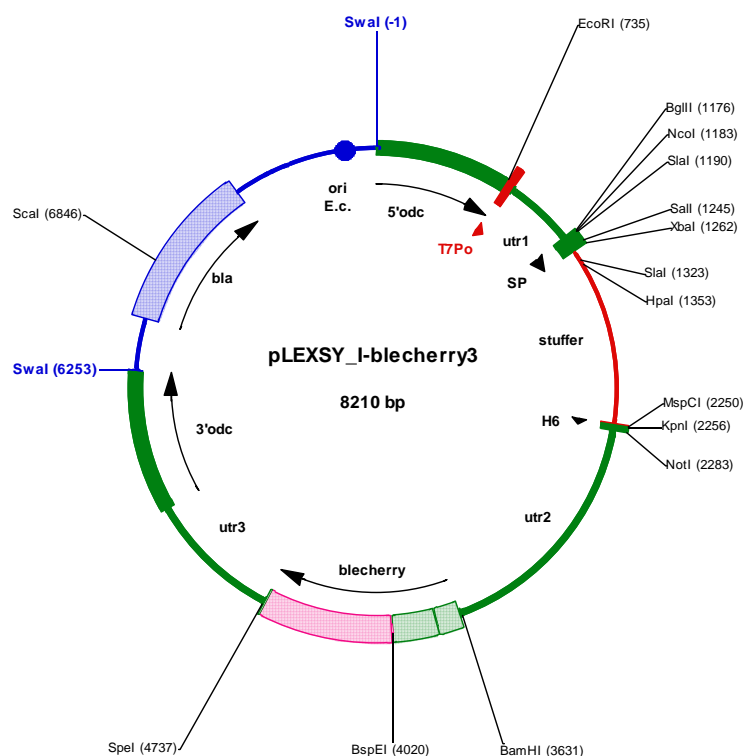
Inducible LEXSY expression vector pLEXY\_I-blecherry3 with the *blecherry* marker and reporter gene consisting of *bleomycin* resistance gene from *S. hindustans* and *cherry fluorescence* gene, enabling selection of recombinant LEXSY strains with the antibiotic LEXSY Bleo (Cat.-No. AB-103) and monitoring induction by cherry fluorescence. Designed for inducible expression of target genes in LEXSY host T7-TR (Cat.-No. LT-110).

### Storage conditions:

Store at -20°C.

Sequence information is available for download from [http://www.jenabioscience.com/cms/en/1/catalog/11\\_16\\_expression\\_vectors.html](http://www.jenabioscience.com/cms/en/1/catalog/11_16_expression_vectors.html)

**FOR RESEARCH USE ONLY.  
 NOT INTENDED FOR HUMAN OR ANIMAL  
 DIAGNOSTIC OR THERAPEUTIC USES.**



Please, find more information in the LEXSinduce3 User's Guide; download .pdf from the LEXSY section of our website [www.jenabioscience.com/LEXY](http://www.jenabioscience.com/LEXY).