

## pLEXSY\_I-egfp-ble3

### LEXSinduce3 Control Vector

Cat.-No.	ble Size
EGE-247	5 µg

#### Description:

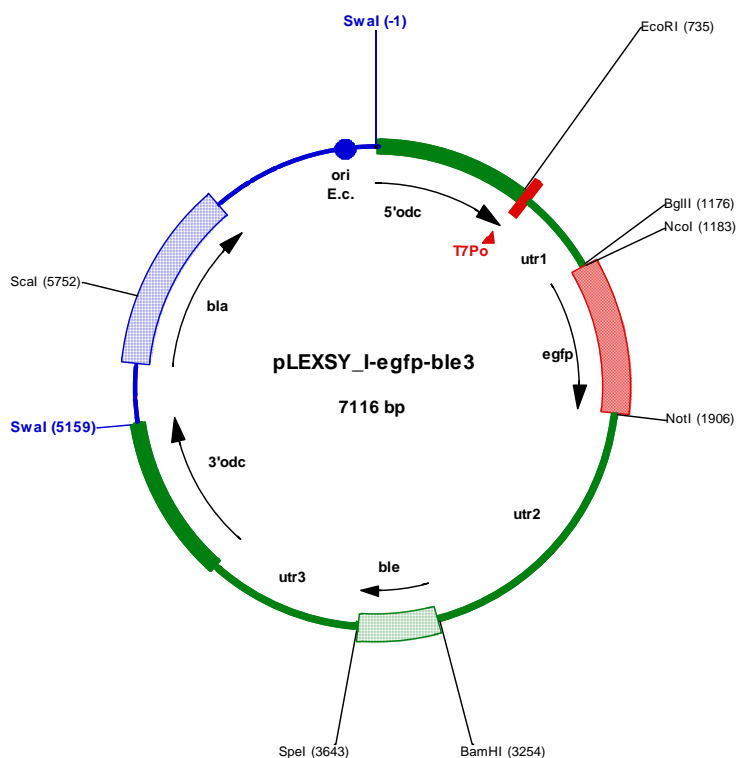
Inducible LEXSY control vector pLEXSY\_I-egfp-ble3 with *ble* marker gene encoding the *bleomycin* resistance gene from *S. hindustans*, enabling selection of recombinant LEXSY strains with the antibiotic LEXSY Bleo (Cat.-No. AB-103) and monitoring induction by EGFP fluorescence. Designed as transfection control for inducible expression 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/LEXSY](http://www.jenabioscience.com/LEXSY).