

pLEXSY_I-egfp-blecherry3

LEXSinduce3 Control Vector

Cat.-No.	Size
EGE-246	5 µg

Description:

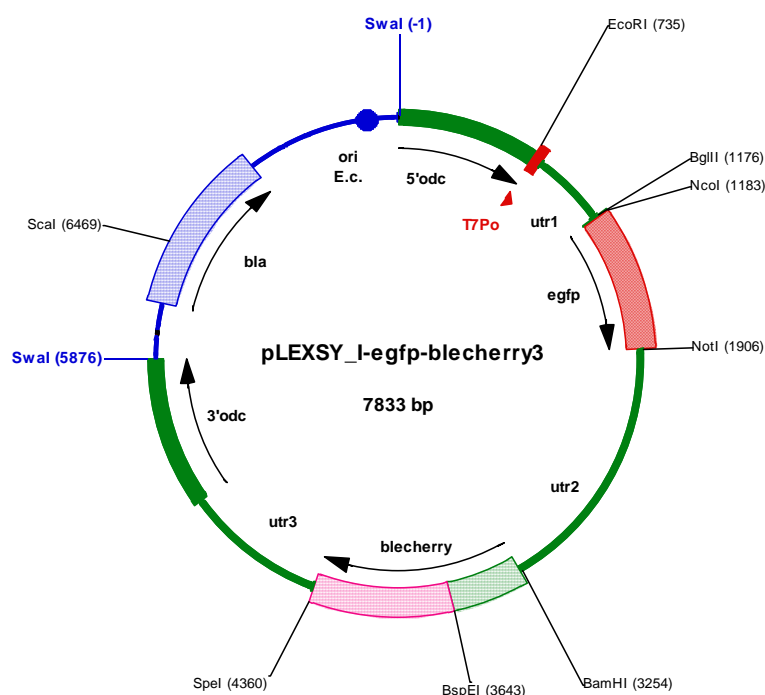
Inducible LEXSY control vector pLEXSY_I-egfp-blecherry3 with the *blecherry* marker and reporter gene consisting of *bleomycin* resistance gene from *S. hindustans* and *cherry fluorescence* gene, enabling selection of recombinant LEXSY strains with the antibiotic LEXSY Bleo (Cat.-No. AB-103) and monitoring induction by Cherry and EGFP fluorescences. 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

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.