

pLEXSY_I-ble3

LEXSinduce Expression Vector

Cat.-No.	Size
EGE-244	5 µg

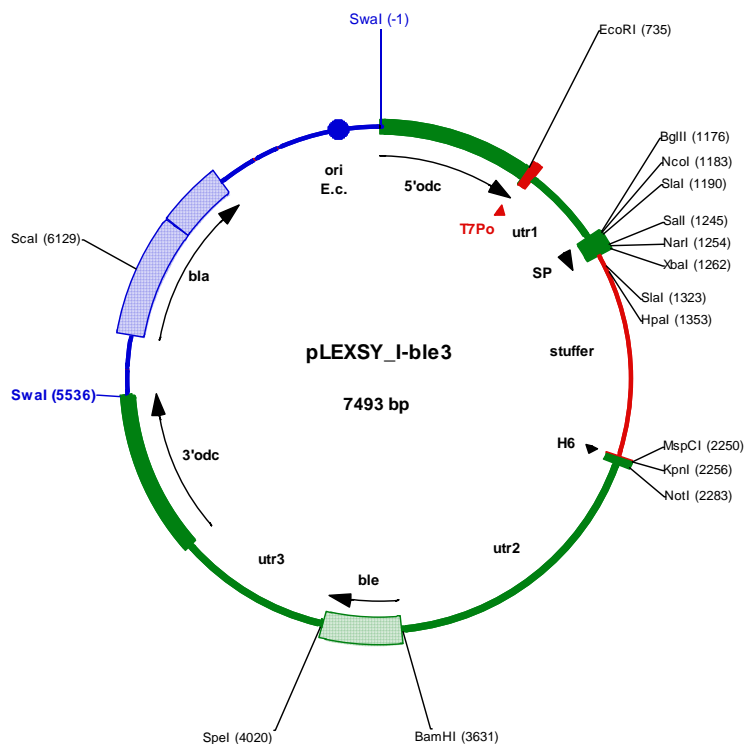
Description:

Inducible LEXSY expression vector pLEXSY_I-ble3 with the *ble* marker gene (*bleomycin resistance* gene from *S. hindustans*), allowing selection of recombinant LEXSY strains with the antibiotic LEXSY Bleo (Cat.-No. AB-103). 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



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.