

pLEXY_I-neo3

LEXSinduce Expression Vector

Cat.-No.	Size
EGE-245	5 µg

Description:

Inducible LEXSY expression vector pLEXY_I-neo3 with the *neo* marker gene (encoding aminoglycoside phosphotransferase 3' (I) or APH(3')I) allowing selection of recombinant LEXSY strains with the antibiotic LEXSY Neo (Cat.-No. AB-105). Designed for inducible expression of target genes in LEXSY host T7-TR (Cat.-No.LT-110).

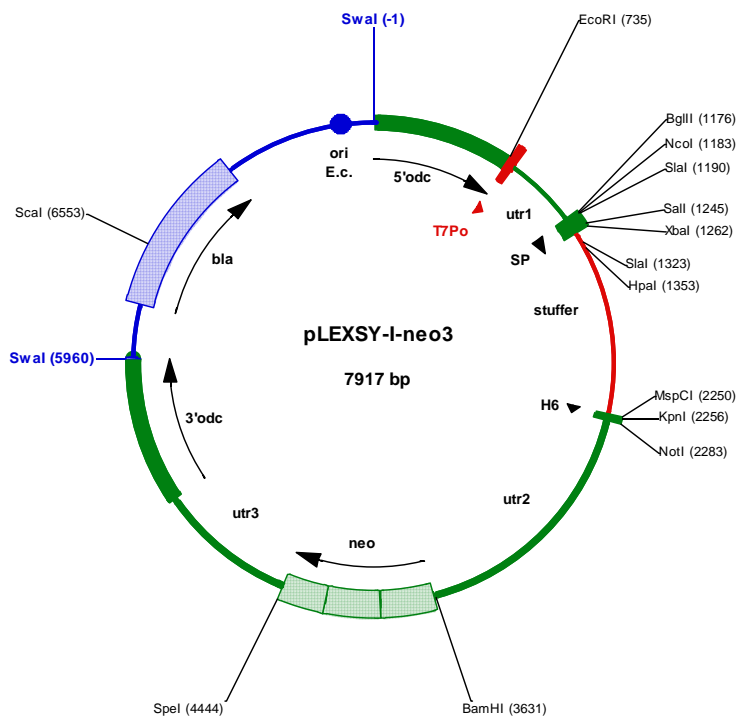
Please note:

The antibiotic LEXSY Neo is different from the anti-bacterial antibiotic Neomycin. The latter MUST NOT BE USED for LEXSY selections.

Storage conditions:

Store at -20°C.

Sequence information is available for download from http://www.jenabioscience.com/cms/en/1/catalog/1116_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.