

pLEXSY_I-egfp-neo3 LEXSinduce Expression Vector

Cat.-No.	Size
EGE-248	5 µg

Description:

Inducible LEXSY control vector pLEXSY_I-egfp 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) and monitoring induction by EGFP fluorescence. Designed as transfection control for inducible expression in LEXSY host T7-TR (Cat.-No. LT-110).

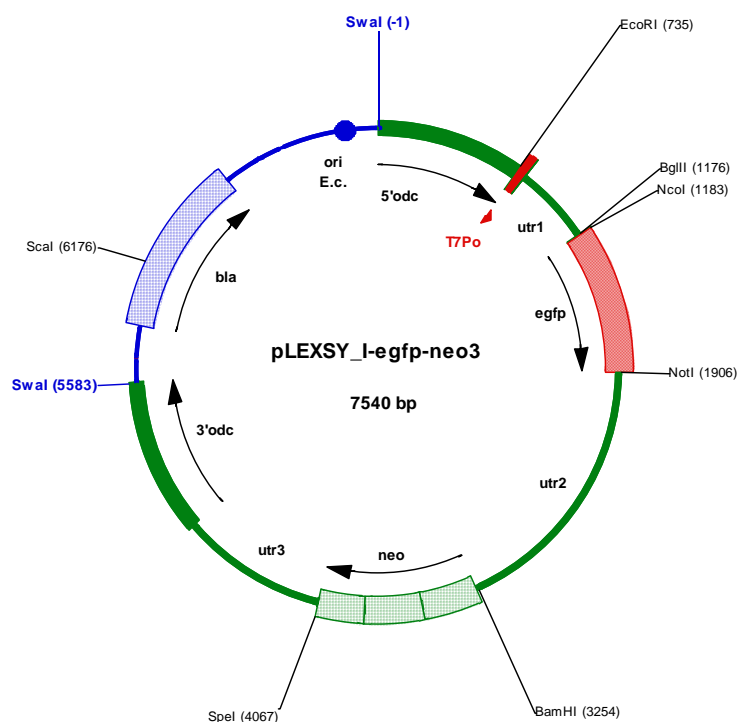
Please note:

The antibiotic LEXSY Neo is different from the antibacterial 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.