

## pLEXY\_I-neo2

### LEXSinduce Expression Vector

Cat.-No.	Size
EGE-242	5 µg

#### Description:

Inducible LEXSY expression vector pLEXY\_I-neo2 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 (LT-110).

#### Storage conditions:

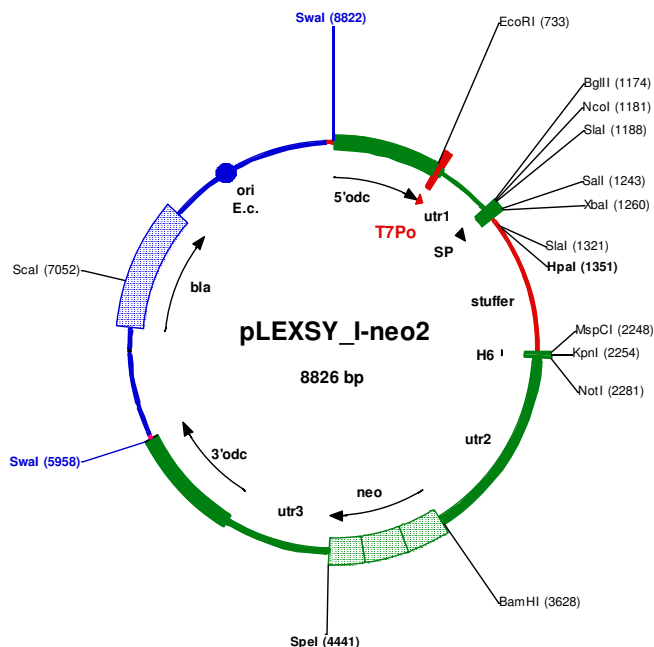
Store at -20°C.

Sequence information available for download from corporate WS [www.jenabioscience.com](http://www.jenabioscience.com)

**FOR RESEARCH USE ONLY.**

**NOT INTENDED FOR HUMAN OR ANIMAL**

**DIAGNOSTIC OR THERAPEUTIC USES.**



Find more information in the LEXSinduce User's Guide (download pdf from the LEXSY section of our website).