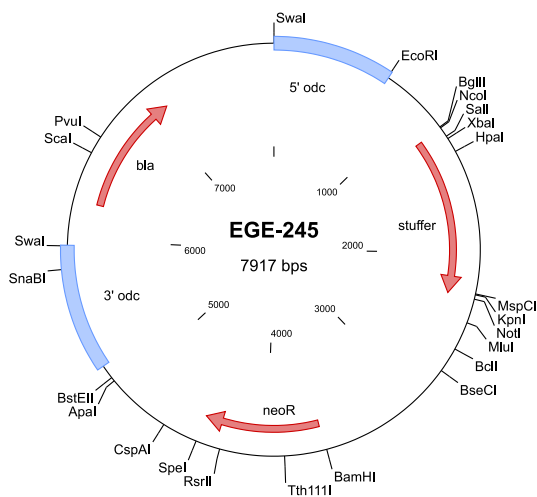




## pLEXY\_I-neo3

Integrative inducible expression vector  
Antibiotic selection of transfectants with LEXSY Neo

Cat. No.	Amount
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 selection.

**For general laboratory use.** Not intended for human or animal diagnostic or therapeutic uses.

**Shipping:** shipped on gel packs

**Storage Conditions:** Store at -20 °C

**Shelf Life:** 12 months