

pLEXYneo2

LEXYcon2 Expression Vector

Cat.-No.	Size
EGE-233	5 µg

Description:

General expression vector pLEXYneo2 with *neo* marker gene encoding aminoglycoside phosphotransferase 3' (I) syn. APH(3')I allowing selection of recombinant LEXSY strains with the antibiotic LEXSY Neo (Cat.-No. AB-105).

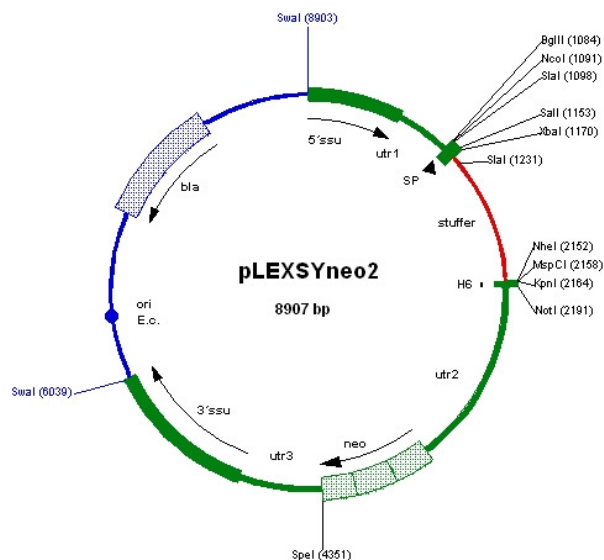
Storage conditions:

Store at -20°C.

Sequence information available for download from corporate WS www.jenabioscience.com

FOR RESEARCH USE ONLY.

**NOT INTENDED FOR HUMAN OR ANIMAL
 DIAGNOSTIC OR THERAPEUTIC USES.**



Find more information in the LEXSYcon2 Expression Kit User's Guide (download pdf from the LEXSY section of our website).