

pLEXY-sat2

LEXYcon2 Expression Vector

Cat.-No.	Size
EGE-234	5 µg

Description:

General expression vector pLEXY-sat2 with *sat* marker gene (encoding streptomycin acetyltransferase) allowing selection of recombinant LEXSY strains with the antibiotic LEXSY NTC (Cat.-No. AB-101 and AB-102).

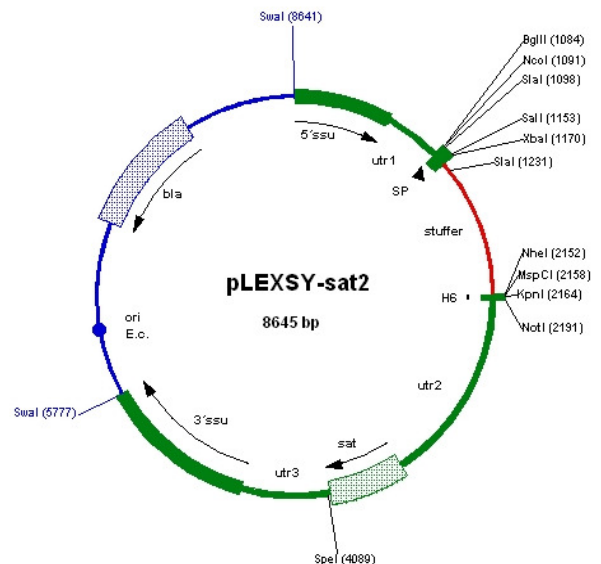
Storage conditions:

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).