

pLEXSY-hyg2

LEXSYcon2 Expression Vector

| Cat.-No. | Size |
|----------|------|
| EGE-232 | 5 µg |

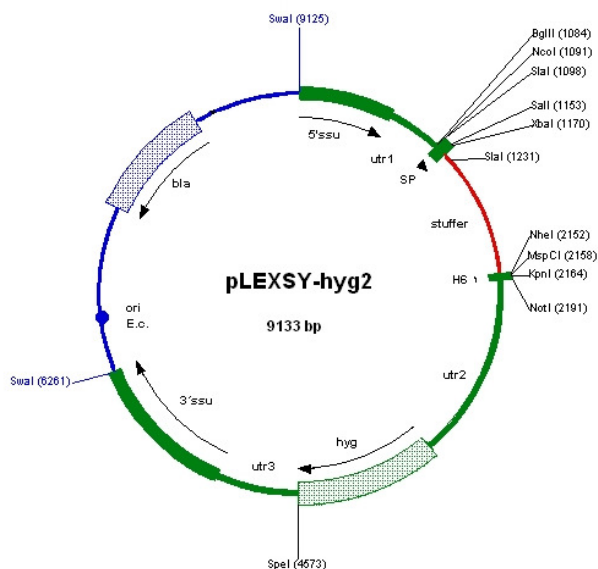
Description:

General expression vector pLEXSY-hyg2 with the *hyg* marker gene (encoding Hygromycin phosphotransferase, HPT) allowing selection of recombinant LEXSY strains with the antibiotic LEXSY Hygro (Cat.-No. AB-104).

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).