

pLEXY-ble2

LEXYcon2 Expression Vector

Cat.-No.	Size
EGE-231	5 µg

Description:

General expression vector pLEXY-ble2 with the *ble* marker gene (*bleomycin resistance gene from S. hindustans*), allowing selection of recombinant LEXSY strains with the antibiotic LEXSY Bleo (Cat.-No.AB-103).

Storage conditions:

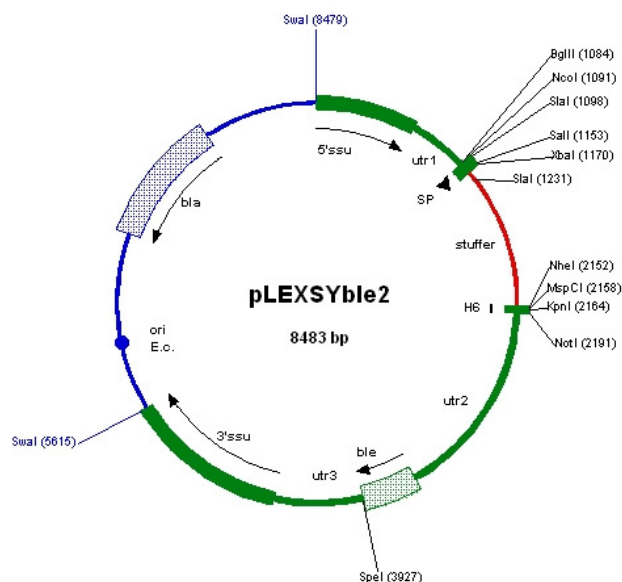
Store at -20°C.

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).