



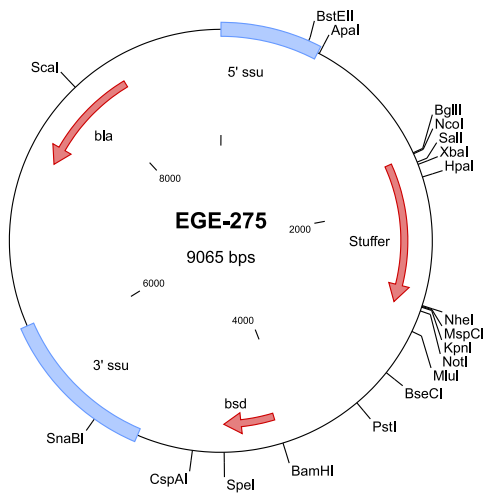
pLEXSY-bsd2.1

Integrative constitutive expression vector
Antibiotic selection of transfectants with LEXSY BSD

Cat. No.	Amount
EGE-275	5 µg

Description:

General expression vector pLEXSY-bsd2.1 with *bsd* marker gene (blasticidin-resistance gene of *Aspergillus terreus*) allowing selection of recombinant LEXSY strains with the antibiotic LEXSY BSD (Cat.-No. AB-107).



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