

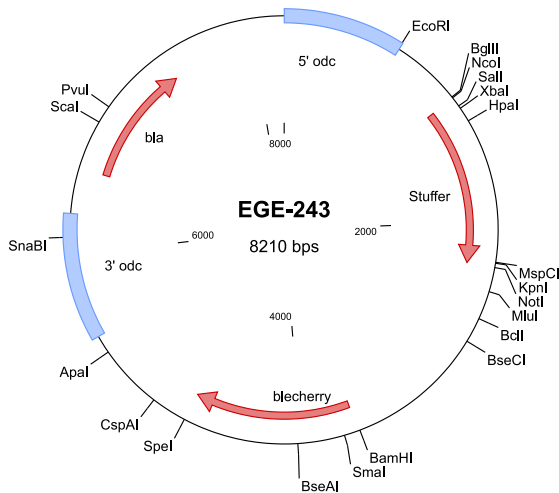


pLEXY_I-blecherry3

Integrative inducible expression vector

Antibiotic selection of transfectants with LEXSY Bleo and expression monitoring with Cherry fluorescence

Cat. No.	Amount
EGE-243	5 µg



Description:

Inducible LEXSY expression vector pLEXY_I-blecherry3 with the *blecherry* marker and reporter gene consisting of *bleomycin* resistance gene from *S. hindustans* and *cherry fluorescence* gene, enabling selection of recombinant LEXSY strains with the antibiotic LEXSY Bleo (Cat.-No. AB-103) and monitoring induction by cherry fluorescence. Designed for inducible expression of target genes in LEXSY host T7-TR (Cat.-No. LT-110).

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