

# Clone and Express Protein-Coding Regions Using the Flexi® Vector Systems

---

## ABSTRACT

*The Flexi® Vector Systems provide a rapid, efficient and high-fidelity method to transfer protein-coding regions between Flexi® Vectors with various expression or peptide-tag options while maintaining insert orientation and reading frame. A lethal gene in the Flexi® Vector cloning region allows positive selection of clones containing insert. Antibiotic resistance genes simplify selection of the desired clone during subsequent transfers to other Flexi® Vectors.*

*Cell Notes* **11**, 13–16.

Michael Slater, Jim Harnett, Natalie Betz, Jami English, Ethan Strauss, Becky Pferdehirt, and Elaine Schenborn  
Promega Corporation  
Publication Date: 2005



## Download Article

(312 KB)

---