

## GeneClip™ U1 Hairpin Cloning Systems for Expression of Short Hairpin RNAs in vivo

## **ABSTRACT**

The GeneClip<sup>TM</sup> U1 Hairpin Cloning Systems enable cloning and expression of short hairpin RNAs (shRNAs) for RNA interference (RNAi) experiments. The GeneClip<sup>TM</sup> Systems contain linearized vectors that facilitate cloning of complementary oligonucleotides downstream of a U1 promoter. Upon transfection with pGeneClip<sup>TM</sup> Vectors containing these cloned sequences, shRNA is produced in the cell, leading to reduction of target gene expression through RNAi. In this article we illustrate use of the GeneClip<sup>TM</sup> Systems to suppress Renilla luciferase expression in a range of cell types and to suppress p53 expression in both a transient and stable manner.

Promega Notes 89, 21-24.

Thomas Yeager, Cheryl Bailey, Jolanta Vidugiriene, Brian Almond and Doug Storts Promega Corporation

Publication Date: 2005



**Download Article** 

(247 KB)