

Achieve the Protein Expression Level You Need with the Mammalian HaloTag 7 Flexi Vectors

ABSTRACT

A series of vectors containing deletions of human cytomegalovirus (CMV) immediate-early enhancer/promoter were created to provide different constitutive expression levels of HaloTag 7 fusion proteins, providing greater control over the level of protein expression. This allows expression of enough fusion protein to image or capture but not so much to disturb normal function of the protein partner fused to the HaloTag 7 protein.

Promega Notes 100, 16–18.

Michael Slater, Danette Hartzell, Jim Hartnett, Sarah Wheeler, Pete Stecha, and Natasha Karassina

Promega Corporation

Publication Date: 2008



Download Article

(142 KB)
