

Expression of Fusion Proteins: How to Get Started with the HaloTag® Technology

ABSTRACT

Appending fusion proteins such as affinity tags and fluorescent proteins is important for the study of protein structure and function. The HaloTag® Technology gives the functionalities of different fusion partners with a single fusion construct. Here we discuss basic considerations and steps to creating a HaloTag® fusion construct.

Promega Notes 100, 13–15.

Melissa McCornack, Trista Schagat, and Michael Slater

Promega Corporation

Publication Date: 2008



Download Article

(103 KB)
