

15N Labeling of Proteins Overexpressed in the Escherichia coli Strain KRX

ABSTRACT

We describe a protocol for efficient 15N-labeled protein production in the E. coli strain, KRX. This procedure can help streamline the process of target cloning, protein expression, screening and preliminary protein folding/aggregation assessment by NMR spectroscopy and other methods that require 15N protein labeling.

Promega Notes 97, 28–29.

Kate Qin Zhao, Jim Hartnett and Michael R. Slater

Promega Corporation

Publication Date: 2009



Download Article

(416 KB)
