

[U-13C, 15N] Protein Labeling Using *Escherichia coli* Strain...

[U-13C, 15N] Protein Labeling Using *Escherichia coli* Strain KRX

ABSTRACT

We have developed a protocol for efficient [U-13C, 15N]labeled protein production in the highly competent *E. coli* strain, KRX, using Silantes CN OD2 medium. This procedure can facilitate protein analysis by NMR spectroscopy and other methodologies that require 13C and 15N protein labeling.

Promega Notes 98, 6–7.

Kate Qin Zhao, Marco Tonelli, Jim Hartnett, and Michael Slater

¹Promega Corporation, ²NMRFAM, University of Wisconsin-Madison

Publication Date: 2008



Download Article

(448 KB)
