

AgarACE™ Agarose-Digesting Enzyme Part Three: DNA/RNA Recovery and Applications Testing...

AgarACE™ Agarose-Digesting Enzyme Part Three: DNA/RNA Recovery and Applications Testing for Use in Translation Transcription/Translation Prime-a-Gene® Labeling and Fluorescent DNA Sequencing

ABSTRACT

AgarACE™ Enzyme is a unique agarase that enables simple, quantitative recovery of nucleic acids from agarose gels. Previously, we presented data describing the use of AgarACE™ Enzyme in cloning, sequencing and PCR-related applications. In this article, we present data on recovery of supercoiled plasmid DNA, high molecular weight linear DNA and RNA, and show that DNA and RNA isolated using AgarACE™ Enzyme perform well in four additional popular applications.

Promega Notes 61, 21.

Mark Knuth, Kimberly Knoche and Paul Otto

Promega Corporation

Publication Date: 1997



Download Article

(316 KB)
