

RNasin® Ribonuclease Inhibitor Part I: Characterization of the Protein

ABSTRACT

Ribonucleases (RNases) are ubiquitous enzymes that can detrimentally affect RNA isolation in the laboratory. Many applications such as RT-PCR and in vitro transcription and translation may be affected as well. RNase can renature after autoclaving and is difficult to inactivate. Many RNase inhibitors are available commercially. Promega's RNasin® Ribonuclease Inhibitor, one of the first introduced, is exceptionally effective at controlling RNase activity during molecular biology procedures.

Promega Notes 77, 8–11.

John Shultz, Robin Hurst and Natalie Betz Promega Corporation Publication Date: 2001



Download Article

(164 KB)