

RNasin® Ribonuclease Inhibitor Part II: A Tale of Two Proteins

ABSTRACT

In this article, we present data that compares RNasin® Ribonuclease Inhibitor to SUPERaseIn™, as RNase inhibitor recently introduced by Ambion. The two RNase inhibitors are compared for their ability to protect RNA from degradation under a variety of different reaction conditions. In addition, we tested both inhibitors for their ability to protect RNA during in vitro translation and transcription.

Promega Notes 77, 12–15.

John Shultz, Robin Hurst, Ken Lewis, Natalie Betz and Ray Williams

Promega Corporation

Publication Date: 2001



Download Article

(220 KB)
