

**RNase Inhibitor - cloned**

mice cell line

Cat. No.	Amount
PCR-504S	2.000 units
PCR-504L	10.000 units

**Unit Definition:** One unit of the protein inhibits the activity of 5 ng RNase A by 50 %. Inhibitor activity is assayed in: 40 mM Tris-HCl (pH 8.0, 25 °C), 10 mM MgSO<sub>4</sub>, 1 mM CaCl<sub>2</sub>.

**For *in vitro* use only!**

**Shipping:** shipped on blue ice

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles

**Shelf Life:** 12 months

**Form:** liquid (Supplied in 20 mM HEPES-KOH pH 7.5, 100 mM KCl, 0.1 mM EDTA and 50 % [v/v] glycerol)

**Concentration:** 40 units/μl

**Applications:**

RNase inhibitor should be applied in a working concentration of 40 units (1 μl) per 50 μl reaction volume. The application of RNase inhibitor is recommended for:

- cDNA synthesis / RT-PCR
- *in vitro* transcription
- *in vitro* translation
- isolation of mammalian cell fractions that contain mRNA-protein complex
- Separation and identification of specific ribonuclease activities
- Studies of tumor suppression

**Description:**

RNase Inhibitor completely inhibits the activity of RNases A, B and C by non-covalent binding. It binds the RNases in a 1:1 ratio with an association constant of approx. 10<sup>14</sup> M. It does not inhibit the RNases I, T1, T2, H, U1, U2 and CL3.