

**Poly (A) carrier RNA**

Polyadenylic acid, potassium salt

Cat. No.	Amount
RNT-017-S	100 mg
RNT-017-L	1 g

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** Short term exposure (up to 1 week cumulative) to ambient temperature possible. Store dark.**Shelf Life:** 12 months**Molecular Weight:** 100 – 500 kDa**Form:** lyophilised solid**Color:** white to off-white**Description:**

Poly (A) carrier RNA is used for quantitative precipitation/purification of RNA and DNA. It improves recovery of short fragments (< 200 bp) or low amounts of nucleic acids.

Chain length: 2100 to 10000 nucleotides

Recommended working concentration: 0.5 mg/ml

Quality control:

Poly (A) in relation to one mononucleotide unit:

2.3 - 2.5 $\mu\text{mol}/\text{mg}$ (A_{257}^* , $\epsilon = 10.7 \text{ L mmol}^{-1} \text{ cm}^{-1}$) A_{290}/A_{260}^* : 0.03 - 0.05 A_{280}/A_{260}^* : 0.25 - 0.35 A_{250}/A_{260}^* : 0.82 - 0.92

DNase/RNase contamination: not detectable (FRET assay)

* absorbance measurement in DNase/RNase free water