



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Yeast tRNA

(Cat. # 786-058, 786-059)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

Transfer RNAs (tRNAs) are used as carriers or co-precipitants for helping in recovery of nucleic acids when alcohol precipitation method is used. Carrier or co-precipitants are inert substances that enhance recovery of nucleic acid when alcohol is used for precipitation. They are insoluble in ethanol or isopropanol solutions and form precipitate that helps to trap nucleic acids and seen as visible pellet after centrifugation. tRNA play significant role in quantitative recovery of small amounts of nucleic acids in dilute solutions. For small amounts of nucleic acid precipitation, ethanol is preferred over isopropanol.

It is also used as a blocking agent in hybridization reactions where RNA probes are used like northern blot. It may also be used in studies that involves use of natural RNA in an *in vivo* and *in vitro* protein synthesizing system

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-058	Yeast tRNA [10 mg/ml]	0.5ml
786-059	Yeast tRNA [10 mg/ml]	1ml

STORAGE CONDITIONS

Shipped on blue ice, store at -20°C upon arrival.

HANDLING

tRNA is sensitive to degradation by exogenous ribonucleases. Therefore one should work in RNAase free zone and always wear gloves when using this product. RNase-free reagents, tubes and filter tips should be used when handling RNA.

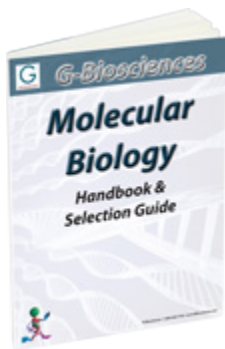
INSTRUCTIONS FOR USE

1. Select the appropriate monovalent cation salt and adjust the concentration of salt in nucleic acid containing solution as per protocol
For example: final concentration to 0.3M Sodium acetate, pH 5.2¹.
2. Add Yeast tRNA to the solution to a final concentration of 10-20 µg/ml and mix.
3. Add 2 volumes of ethanol, mix well and chill at -20°C for 15 minutes.
4. Centrifuge at $\geq 10,000 \times g$ for 15 minutes to precipitate the nucleic acid.
5. Discard the supernatant and dissolve the precipitate in DI water or appropriate buffer.

REFERENCES

1. Maniatis, T. et al (1982). Molecular Cloning. A Laboratory Manual. New York: Cold Spring Harbor Laboratory

RELATED PRODUCTS



Download our Molecular Biology Handbook. <http://info2.gbiosciences.com/complete-molecular-biology-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 7/29/2016 CMH



www.GBiosciences.com