

Certificate of Analysis

Ribo m⁷G Cap Analog:

Part No.	Size (units)
P171A	10 A ₂₅₄ units
P171B	25 A ₂₅₄ units

Description: 5' 7-methyl guanosine nucleotide (m⁷G(5')ppp(5')G) or cap structure is incorporated into RNA synthesized in vitro to mimic the capped structure of mRNA. This product is intended for use with Riboprobe® and RiboMAX™ Systems.

Formula: C₂₁H₂₇N₁₀O₁₈P₃Na₂.

Formula Weight: 846.

Storage Conditions: See the product information label for storage recommendations. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes.

Volume: Part# P171A contains at least 13.7µl Ribo m⁷G Cap Analog in Nuclease-Free Water. Part# P171B contains at least 34.3µl Ribo m⁷G Cap Analog in Nuclease-Free Water.

Part# 9PIP171

Revised 4/18



AF9PIP171 0418P171

Quality Control Assays

Activity Assays

Concentration: 40 ± 2mM, as determined by absorbance at 254nm.

Functional Assay: The Ribo m⁷G Cap Analog is tested for capped transcript synthesis using T7 RNA Polymerase and a 10:1 ratio of Ribo m⁷G Cap Analog to GTP. Capped and uncapped transcripts are separated by gel electrophoresis using a denaturing polyacrylamide gel. The minimum passing specification is > 50% capped transcripts.

References

1. Nakagawa, I. *et al.* (1980) A "capping" agent: P¹-S-Phenol P²-7-methylguanosine-5' pyrophosphorothioate. *Synthesis* 556.
2. Krieg, P.A. and Melton, D.A. (1987) In vitro RNA synthesis with SP6 RNA polymerase. *Meth. Enzymol.* **155**, 397-415.
3. Paterson, B.M. and Rosenberg, M. (1979) Efficient translation of prokaryotic mRNAs in a eukaryotic cell-free system requires addition of a cap structure. *Nature* **279**, 692-6.
4. Drummond, D.R., Armstrong, J. and Colman, A. (1985) The effect of capping and polyadenylation on the stability, movement and translation of synthetic messenger RNAs in *Xenopus oocytes*. *Nucl. Acids Res.* **13**, 7375-94.



Promega

Promega Corporation

2800 Woods Hollow Road	
Madison, WI 53711-5399 USA	
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	www.promega.com

PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

© 1999, 2001, 2005, 2006, 2010, 2013, 2016, 2018 Promega Corporation. All Rights Reserved.

pGEM, Riboprobe, RNasin and TNT are registered trademarks of Promega Corporation. RiboMAX is a trademark of Promega Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIP171
Printed in USA. Revised 4/18

Signed by:

R. Wheeler, Quality Assurance

1. Standard Applications: Synthesis in vitro of Capped RNA Transcripts

Materials to Be Supplied by the User

All materials can be found in Section 3.

(Solution compositions are provided in Section 2.)

- DNA template, linearized
- Nuclease-Free Water
- Recombinant RNasin® Ribonuclease Inhibitor
- rNTP capping mix
- Transcription Optimized 5X Buffer
- Phage RNA Polymerase

1. In a microcentrifuge tube, add the following reagents at room temperature in the order listed:

Transcription Optimized 5X Buffer	10µl
DTT, 100mM	5µl
Recombinant RNasin® Ribonuclease Inhibitor	50 units
rNTP capping mix (see Section 2)	5µl
Ribo m ⁷ G Cap Analog, 5mM	5µl
DNA template, linearized (in water or TE buffer)	5µg
Phage RNA Polymerase	40 units
Nuclease-Free Water to final volume of	50µl

2. Incubate for 1 hour at 37°C. To increase the yield of RNA, add an additional 40 units of Phage RNA Polymerase and incubate for 1 hour.

2. Composition of Buffers and Solutions

rNTP capping mix

5mM	rATP
5mM	rCTP
5mM	rUTP
0.5mM	rGTP

in Nuclease-Free Water

Transcription Optimized 5X Buffer

200mM	Tris-HCl (pH 7.9 at 25°C)
50mM	NaCl
30mM	MgCl ₂
10mM	spermidine

5mM Ribo m⁷G Cap Analog

5mM Ribo m⁷G Cap Analog

in Nuclease-Free Water

3. Related Products

A. Related Systems

Product	Cat.#
Riboprobe® System—SP6	P1420
Riboprobe® System—T3	P1430
Riboprobe® System—T7	P1440
Riboprobe® System Buffers	P1121
RiboMAX™ Large Scale RNA Production System—SP6	P1280
RiboMAX™ Large Scale RNA Production System—T7	P1300
TnT® SP6 Coupled Reticulocyte Lysate System	L4600
TnT® T3 Coupled Reticulocyte Lysate System	L4950
TnT® T7 Coupled Reticulocyte Lysate System	L4610
TnT® T7/SP6 Coupled Reticulocyte Lysate System	L5020
TnT® T7/T3 Coupled Reticulocyte Lysate System	L5010
TnT® SP6 Coupled Reticulocyte Lysate System, Trial Size	L4601
TnT® T7 Coupled Reticulocyte Lysate System, Trial Size	L4611
TnT® T7 Quick Coupled Transcription/Translation System	L1170
TnT® T7 Quick Coupled Transcription/Translation System, Trial Size	L1171
TnT® SP6 Quick Coupled Transcription/Translation System	L2080
TnT® SP6 Quick Coupled Transcription/Translation System, Trial Size	L2081

B. Related Products

Product	Size	Cat.#
SP6 Promoter Primer	2µg	Q5011
pGEM® Express Positive Control Template	10µg (2 × 5µg)	P2561
SP6 RNA Polymerase	5,000u	P1081
	1,000u	P1085
SP6 RNA Polymerase (HC)	2,500u	P4084
T7 RNA Polymerase	5,000u	P2077
	1,000u	P2075
T7 RNA Polymerase (HC)	10,000u	P4074
T3 RNA Polymerase	1,000u	P2083
T3 RNA Polymerase (HC)	2,500u	P4024
rATP, 10mM	0.5ml	P1132
rCTP, 10mM	0.5ml	P1142
rGTP, 10mM	0.5ml	P1152
rUTP, 10mM	0.5ml	P1162
Transcription Optimized 5X Buffer	200µl	P1181
rATP, rCTP, rGTP and rUTP, each at 10mM	0.5ml each	P1221
Nuclease-Free Water	50ml (2 × 25ml)	P1193

Product	Conc.	Size	Cat.#
Recombinant RNasin®			
Ribonuclease Inhibitor	20–40u/µl	2,500u	N2511
	20–40u/µl	10,000u	N2515