

# Flexi® Rabbit Reticulocyte Lysate System

INSTRUCTIONS FOR USE OF PRODUCT L4540.

**Quick**  
PROTOCOL

## In Vitro Translation Protocol

### Before You Begin

Remove the reagents from  $-70^{\circ}\text{C}$  storage. Rapidly thaw the lysate by hand-warming, and immediately store on ice. Thaw other components at  $37^{\circ}\text{C}$ , and store on ice.

### Preparation of Template RNA

Denature the template mRNA at  $65^{\circ}\text{C}$  for 3 minutes, and immediately cool on ice. The template should be free of ethanol, calcium, RNases and salts.

### Translation Procedure

1. Assemble the reaction components, appropriate for the label being used, in a 0.5ml polypropylene microcentrifuge tube. After all components are added, gently mix the lysate by pipetting and stirring the reaction with the pipette tip. If necessary, centrifuge briefly to return the sample to the bottom of the tube.

Component	Reaction Using [ $^{35}\text{S}$ ]methionine	Reaction Using Transcend™ tRNA
Flexi® Rabbit Reticulocyte Lysate	33 $\mu\text{l}$	33 $\mu\text{l}$
Amino Acid Mixture, Minus Leucine, 1mM	—	0.5 $\mu\text{l}$
Amino Acid Mixture, Minus Methionine, 1mM	1 $\mu\text{l}$	0.5 $\mu\text{l}$
[ $^{35}\text{S}$ ]methionine (1,200Ci/mmol at 10mCi/ml)	2 $\mu\text{l}$	—
Magnesium Acetate, 25mM	0–4 $\mu\text{l}$	0–4 $\mu\text{l}$
Potassium Chloride, 2.5M	1.4 $\mu\text{l}$	1.4 $\mu\text{l}$
DTT, 100mM	0–1 $\mu\text{l}$	0–1 $\mu\text{l}$
RNasin® Ribonuclease Inhibitor (40u/ $\mu\text{l}$ )	1 $\mu\text{l}$	1 $\mu\text{l}$
RNA substrate	1–12 $\mu\text{l}$	1–12 $\mu\text{l}$
Transcend™ tRNA	—	1–2 $\mu\text{l}$
Nuclease-Free Water to a final volume of	50 $\mu\text{l}$	50 $\mu\text{l}$

2. Incubate the translation reaction at  $30^{\circ}\text{C}$  for 90 minutes.
3. Analyze the results of translation. See Technical Bulletin #TB127 for procedures for incorporation assays, gel analysis of translation products and an assay for luciferase production in the control reactions.

### Control Translation Reactions

1. We recommend including a control reaction containing no added mRNA. This allows measurement of any background incorporation of labeled amino acids.
2. The protocol for a positive control reaction using the Luciferase Control RNA is on the opposite side of this card.

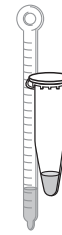
See additional protocol information in Technical Bulletin #TB127, available online at: [www.promega.com](http://www.promega.com)

### ORDERING/TECHNICAL INFORMATION:

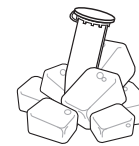
[www.promega.com](http://www.promega.com) • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601



Keep all components on ice.



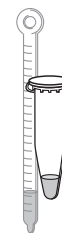
Incubate mRNA at  $65^{\circ}\text{C}$  for 3 minutes. Keep other reagents on ice.



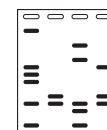
Immediately cool mRNA on ice.



Assemble reaction components. Gently mix. Return unused components to  $-70^{\circ}\text{C}$ .



Incubate at  $30^{\circ}\text{C}$  for 90 minutes.



Analyze results.

3061MA09\_0A



**Promega**

Printed in USA Revised 3/09.  
Part #9FB041

# Flexi® Rabbit Reticulocyte Lysate System

INSTRUCTIONS FOR USE OF PRODUCT L4540.

**Quick**  
PROTOCOL

## Luciferase Positive Control Translation Reaction

### Before You Begin

Remove the reagents from  $-70^{\circ}\text{C}$  storage. Rapidly thaw the lysate by hand-warming, and immediately store on ice. Thaw other components at  $37^{\circ}\text{C}$ , and store on ice.



Keep all components on ice.

### Non-Radioactive Luciferase Control Reaction

1. Assemble the following reaction:

Component	Volume
Flexi® Rabbit Reticulocyte Lysate	35 $\mu\text{l}$
Amino Acid Mixture, Minus Leucine, 1mM	0.5 $\mu\text{l}$
Amino Acid Mixture, Minus Methionine, 1mM	0.5 $\mu\text{l}$
Potassium Chloride, 2.5M	1.4 $\mu\text{l}$
RNasin® Ribonuclease Inhibitor (40u/ $\mu\text{l}$ )	1 $\mu\text{l}$
Luciferase Control RNA, 1mg/ml	1 $\mu\text{l}$
Nuclease-Free Water	<u>10.6<math>\mu\text{l}</math></u>
final volume	50 $\mu\text{l}$

2. Incubate the translation reaction at  $30^{\circ}\text{C}$  for 60–90 minutes.

The luciferase control reactions can be stored at  $-20^{\circ}\text{C}$  for up to 2 months or at  $-70^{\circ}\text{C}$  for up to 6 months with little loss of luciferase activity.

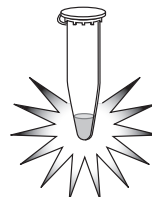
3. Detect the synthesis of functional luciferase using the standard luciferase assay (see Section 7 of #TB127).



Assemble reaction components. Gently mix. Return unused components to  $-70^{\circ}\text{C}$ .



Incubate at  $30^{\circ}\text{C}$  for 60–90 minutes.



Perform luciferase assay.

### Radioactive Luciferase Control Reaction

1. Assemble the following reaction:

Component	Volume
Flexi® Rabbit Reticulocyte Lysate	35 $\mu\text{l}$
Amino Acid Mixture, Minus Methionine, 1mM	1 $\mu\text{l}$
[ $^{35}\text{S}$ ]methionine (1,200Ci/mmol) at 10mCi/ml	2 $\mu\text{l}$
Potassium Chloride, 2.5M	1.4 $\mu\text{l}$
RNasin® Ribonuclease Inhibitor (40u/ $\mu\text{l}$ )	1 $\mu\text{l}$
Luciferase Control RNA, 1mg/ml	<u>1<math>\mu\text{l}</math></u>
Nuclease-Free Water to a final volume of	50 $\mu\text{l}$

2. Follow Steps 2–3 for the non-radioactive luciferase control reaction.

See additional protocol information in Technical Bulletin #TB127, available online at: [www.promega.com](http://www.promega.com)

3064MA09\_0A

### ORDERING/TECHNICAL INFORMATION:

[www.promega.com](http://www.promega.com) • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601



**Promega**

Printed in USA Revised 3/09.  
Part #9FB041

© 2000, 2009 Promega Corporation. All Rights Reserved.