

# RayBio<sup>®</sup> Magnesium Assay Kit

User Manual Version 1.0  
January 15, 2015

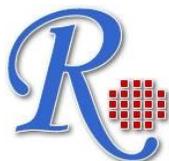
RayBio<sup>®</sup> Magnesium Assay  
Kit Protocol

(Cat#:68-Mag-S100)



We Provide You With Excellent  
Support And Service

Tel:(Toll Free)1-888-494-8555 or 770-729-2992; Fax:770-206-2393;  
Web: [www.raybiotech.com](http://www.raybiotech.com) Email: [info@raybiotech.com](mailto:info@raybiotech.com)



RayBiotech, Inc.

---

## RayBio<sup>®</sup> Magnesium Assay Kit

### TABLE OF CONTENTS

I. Introduction.....	1
II. Kit Contents.....	2
III. Storage and Handling.....	2
IV. Reagent Preparation .....	2
V. Magnesium Assay Protocol.....	2

## I. INTRODUCTION

Magnesium is the 11th most abundant element by mass in the human body.  $Mg^{+2}$  is essential to all living cells where it plays an important role in facilitating the processing of biological polyphosphates like ATP, DNA, RNA and enzyme functions.  $Mg^{+2}$  is the metallic ion at the center of chlorophyll, and a common additive to fertilizers.  $Mg^{+2}$  compounds are used as laxatives, antacids, and used to stabilize abnormal nerve excitation and blood vessel spasm i.e., eclampsia. The RayBiotech Magnesium Assay Kit provides a simple sensitive means of quantitating magnesium in a variety of biological samples. The kit takes advantage of the specific requirement of glycerol kinase for  $Mg^{+2}$ . An enzyme linked reaction leads to formation of an intensely colored ( $\lambda_{max} = 450nm$ ) product whose formation is proportional to  $Mg^{+2}$  concentration. The linear range of the assay is 2-15 nmoles with detection sensitivity~ 40  $\mu M$ .

## II. KIT CONTENTS

Store Kit at -20 °C

Components	Size	Cap Code	Part Number
Magnesium Assay Buffer	25 ml	WM	Item A
Magnesium Developer	Lyophilized	Red	Item B
Magnesium Enzyme Mix	Lyophilized	Green	Item C
Magnesium Standard (150 nmol/μl)	0.1 ml	Yellow	Item D

## III. STORAGE AND HANDLING

Store kit at -20°C, protect from light. Warm buffer to room temperature before use. Briefly centrifuge all small vials prior to opening.

## IV. REAGENT PREPARATION

**Developer:** Dissolve with 1.1 ml dH<sub>2</sub>O. Stable for two months at 4°C.

**Magnesium Enzyme Mix:** Dissolve in 550 μl Assay Buffer. Aliquot and store at -20 °C. Use within two months.

**Magnesium Standard:** Ready to use as supplied. 150 nmol/μl of Mg<sup>+2</sup> Standard stock solution. Store at -20 °C. Mix before each use.

## VII. MAGNESIUM ASSAY PROTOCOL

- 1) Standard Curve Preparations:** Dilute the standard to 1.5 nmol/μl by adding 10 μl of the 150 nmol/μl Magnesium Standard to 990 μl of distilled water, mix well. Add 0, 2, 4, 6, 8, 10 μl into a series of wells. Adjust volume to 50 μl/well with distilled water to generate 0, 3, 6, 9, 12, 15 nmol/well of Magnesium Standard.
- 2) Sample Preparation:** Tissue or cells can be extracted with 4 volume of Magnesium Assay Buffer, spin 16000g for 10 min to get clear extract. Add 1-50 μl of liquid sample into 96 well plate, bring total volume to 50 μl with water.

Normal serum contains  $Mg^{2+}$  0.7-1.05 mM (1.65-2.55 mg/dL), use 5  $\mu$ l serum for testing. Urine should be diluted 10X. For unknown samples, we suggest testing different amount of samples to ensure OD is in the linear range.

**3) Magnesium Reaction Mix:** Mix enough reagent for the number of samples and standards to be performed: For each well, prepare a total 50  $\mu$ l Reaction Mix containing:

35  $\mu$ l Magnesium Assay Buffer  
10  $\mu$ l Developer  
5  $\mu$ l Magnesium Enzyme Mix

**4)** Add 50  $\mu$ l of the Reaction Mix to each well containing the Magnesium Standard and test samples. For best results, use a multichannel pipettor to initiate reaction in all samples at the same time. Mix well.

**5)** Incubate at 37°C for 10 min. Read the plate OD<sub>450nm</sub> to get  $A_0$  for each standard or sample.

*Notes:*

- Since enzyme kinetics are sensitive to temperature variation, the reaction rate will increase as the temperature rises. The reaction takes ~ 10 minutes to reach a linear reaction rate.
- NAD(P)H etc. in samples may generate background, the 10 min waiting time can correct these nonspecific background.
- $Mn^{2+}$ ,  $Zn^{2+}$ ,  $Ni^{2+}$ ,  $Fe^{2+}$ ,  $Cu^{2+}$ ,  $Co^{2+}$ ,  $Ca^{2+}$  do not interfere with the assay.

**6)** Incubate the reaction for additional 10-30 min, read the OD again to get reading A. We recommend monitor the reaction kinetics to ensure the readings are in linear range when read the plate for the additional 10-30 minutes. All readings should not exceed 1.5 OD.

**7) Calculation:** Subtract  $A_0$  from standard and sample readings to get  $\Delta OD = A - A_0$ . Plot Magnesium standard curve. Apply sample  $\Delta OD$  to the standard curve to get  $Mg^{2+}$  amount B (nmol) in the reaction well.  $Mg^{2+}$  concentration:

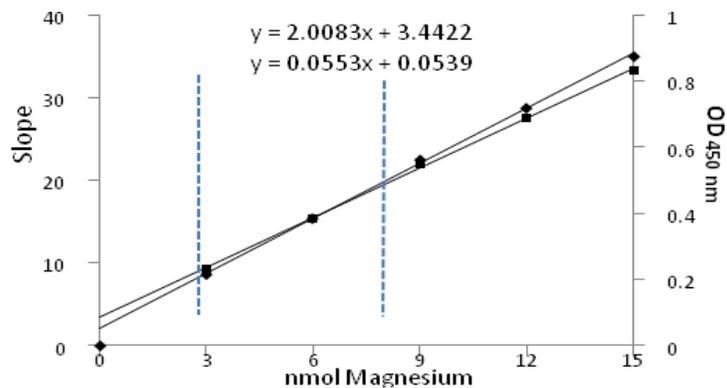
$$C = B/V \quad (\text{nmol/ml or } \mu\text{M})$$

Where: **B** is  $Mg^{2+}$  amount in reaction well (in nmol).

**V** is the sample volume added into the reaction well (in ml).

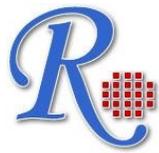
Magnesium molecular weight: 24.3 g/mol, 1mM=2.43 mg/dL.

The assay may also be calculated by monitoring reaction slopes in the standards and samples reactions.



**Magnesium standard curve:** Assay is performed according to kit protocol. Vertical dotted lines indicate the lower and upper limits of normal serum  $Mg^{2+}$  concentrations.

This product is for research use only.



*©2004 RayBiotech, Inc.*