



Product Manual

BIOMOL[®] GREEN

Reagent for phosphate detection

Catalog #: BML-AK111



Product Manual

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON- INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURE**

TABLE OF CONTENTS



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

Introduction	2
Materials Supplied.....	2
Storage	2
Stability	2
Additional Materials Needed	3
Precautions.....	3
Reagent Preparation.....	3
I. Microplate Assay	3
II. Cuvette Assay	5
Data Analysis	6
Reagent Stability.....	7
Notes.....	8
Contact Information.....	10

INTRODUCTION

The *BIOMOL[®] GREEN Reagent* is a convenient 1-step reagent for measuring free-phosphate released during enzymatic phosphatase assays. It is a modification of the classic Malachite green reagent^{1, 2} and it offers the following advantages:

- Non-Radioactive
- Convenient 1-Step Reagent -No Mixing!
- Long Shelf-Life > 6 Months at 4°C
- Excellent Sensitivity (See Below)
- Suitable For Microplate Assays (100 µL Volume) or Cuvette Assays (1 mL Volume)

Each *BIOMOL[®] GREEN Reagent* bottle includes 1 vial of phosphate standard (1 mL) suitable for several phosphate standard curves.

References:

1. B. Martin *et al. J. Biol. Chem.* 1985 **260** 14932
2. K.W. Harder *et al. Biochem. J.* 1994 **298** 395

MATERIALS SUPPLIED

<i>BIOMOL[®] GREEN</i> Reagent (BML-AK111-1000) or (BML-AK111-0250)	1000 mL or 250 mL
Phosphate Standard in dH ₂ O (BML-KI102) (800 µM in distilled water)	1 mL

STORAGE

Room Temperature. Long-term at +4°C. *Please note that BIOMOL[®] GREEN can be stored frozen without deleterious effects.*

***Note: Due to potential expansion during freezing, it is recommended that the contents be divided into smaller volumes prior to freezing.**

STABILITY

When used and stored as directed, the kit is stable until the expiration date indicated on the box.

ADDITIONAL MATERIALS NEEDED

- Microplate reader or visible wavelength spectrophotometer capable of measuring absorbance at 600 to 680 nm to ≥ 3 -decimal accuracy. Although OD_{620nm} is used in the protocols and examples below, the colored phosphate complex formed in the reaction has a broad absorbance peak with a maximum at ~ 650 nm. With appropriate calibration with the Phosphate Standard, any wavelength in the 600 to 680 nm range may be used in the assay.
- $\frac{1}{2}$ -Volume or standard 96-well microplate (ENZO catalog no. ADI-80-2404) or 1 mL cuvette (plastic/glass/quartz)
- Pipetman capable of pipetting 5 μ L -1 mL accurately
- Multi-channel pipettor capable of pipetting 100 μ L (optional).



PRECAUTIONS

Important/ Warning

The BIOMOL[®] GREEN reagent is a highly sensitive phosphate detection solution. Free phosphate present on labware and in reagent solutions will greatly increase the background absorbance of the assay. This is detected visually as a change in color from yellow to green. Detergents used to clean labware may contain high levels of phosphate. Use caution by either rinsing labware with dH_2O or employ unused plasticware.

REAGENT PREPARATION

I. Microplate Assay (Linear Range: 0.03 nmol to 2 nmol):

1. Prepare 1:1 serial dilutions of phosphate standard (BML-KI102) plus a blank (50 μ L per well). Dilutions of 40, 20, ... 0.625 μ M correspond to 2, 1, ...0.031 nmol PO_4 (see Table 1):
 - a. Dilute 800 μ M phosphate standard (BML-KI102) 1/20 into the experimental assay buffer (e.g., 20 μ L phosphate + 380 μ L assay buffer).
 - b. Add 100 μ L of the 40 μ M phosphate standard to sample #1 of microtiter-plate (1/2 volume or standard).
 - c. Add 50 μ L assay buffer to samples #2 through #8.
 - d. Remove 50 μ L from sample #1 and add it to sample #2.
 - e. Mix thoroughly by pipetting up and down several times.
 - f. Remove 50 μ L from sample #2 and add it to sample #3.
 - g. Mix thoroughly and repeat for samples #4-7. At sample #7, remove 50 μ L and discard. DO NOT PROCEED TO sample #8 (blank).
2. Prepare experimental phosphatase samples in a total volume of 50 μ L.

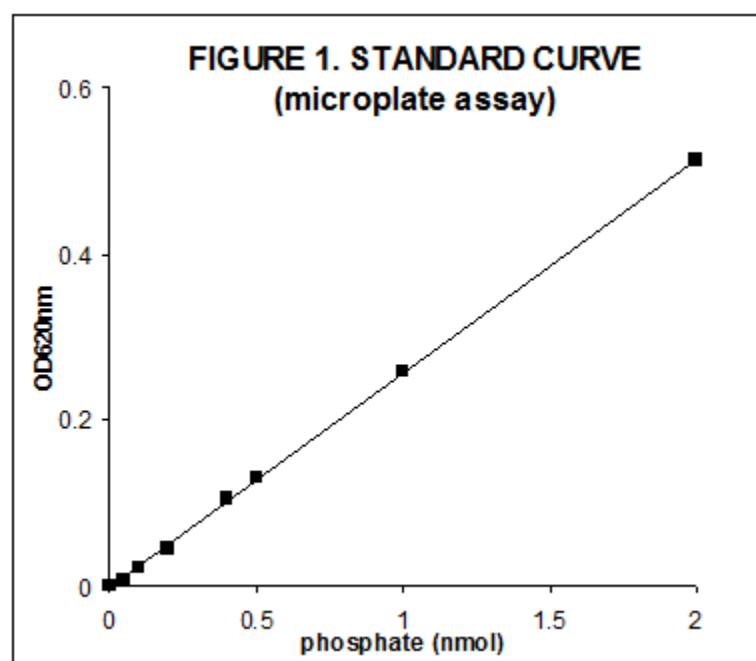
Sample #	†Well #	Phosphate (nmol)
1	A1,2	2
2	B1,2	1
3	C1,2	0.5
4	D1,2	0.25
5	E1,2	0.125
6	F1,2	0.063
7	G1,2	0.031
8	H1,2	0

† For highest accuracy, perform all samples in duplicate.

Table 1. Example of Microtiter Plate Standard Curve Samples

To Terminate Reaction:

1. Add 100 μ L *BIOMOL*[®] *GREEN* Reagent (BML-AK111) to each well of the microtiter plate (Hint: Use a multichannel pipetman). The experiment should be performed such that all wells will be terminated at approximately the same time.
2. Incubate plate at room temperature for 20-30 minutes to allow development of the green color.
3. Measure OD_{620nm} on a microtiter-plate reader (see Figure 1, 1/2 volume microtiter-plate, below).
4. Perform data analysis (see below).

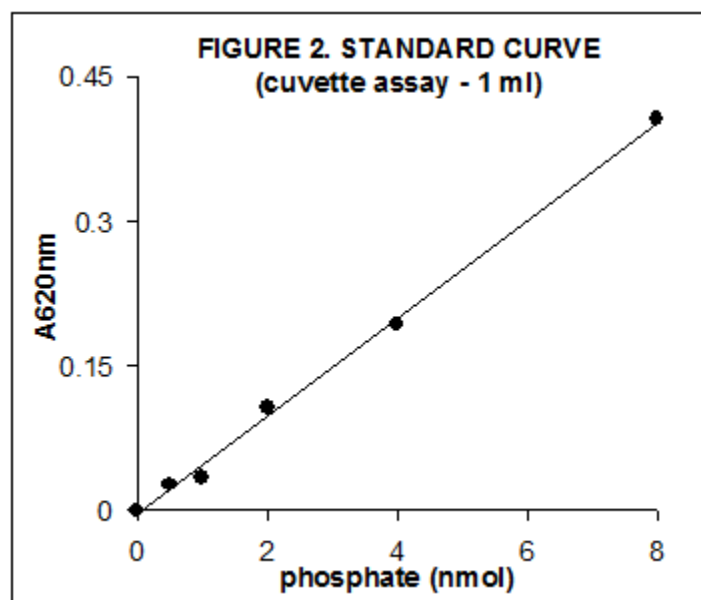


II. Cuvette Assay (Linear Range: 0.05 nmol to 8 nmol):

1. Prepare in tubes 1:1 serial dilutions of phosphate standard (BML-KI102) plus a blank (100 μ L per tube). Dilutions of 80, 40, ...5 μ M correspond to 8, 4, ...0.5 nmol PO_4 :
 - a. Dilute 800 μ M phosphate standard (BML-KI102) 1/10 into the experimental assay buffer (e.g., 50 μ L phosphate + 450 μ L assay buffer).
 - b. Add 200 μ L of the 80 μ M phosphate standard to sample #1.
 - c. Add 100 μ L assay buffer to samples #2 through #6.
 - d. Remove 100 μ L from sample #1 and add it to sample #2.
 - e. Vortex thoroughly.
 - f. Remove 100 μ L from sample #2 and add it to sample #3.
 - g. Vortex thoroughly and repeat for samples #4-5. At sample #5, remove 100 μ L and discard. DO NOT PROCEED TO SAMPLE #6 (blank).
2. Prepare experimental phosphatase samples in a total volume of 100 μ L.

To Terminate Reaction:

1. Add 1 mL *BIOMOL*[®] *GREEN* Reagent (BML-AK111) to each sample. The experiment should be performed such that all wells will be terminated at approximately the same time.
2. Incubate samples at room temperature for 20-30 minutes to allow development of the green color.
3. Measure $\text{OD}_{620\text{nm}}$ using 1 mL cuvettes (plastic/glass/quartz) on a visible-wavelength spectrophotometer (see Figure 2, below).
4. Perform data analysis (see below).



DATA ANALYSIS

Phosphate (PO₄) Standard Curve:

NOTE: For highest accuracy, a standard curve must be performed for each new set of assay data. This will normalize for variations in free phosphate in samples, time of incubation with the BIOMOL® GREEN reagent, and other experimental factors.

1. Plot standard curve data as OD_{620nm} versus nmol PO₄ (see Figure 1 and 2).
2. Obtain a line-fit to the data using an appropriate routine.
3. Convert OD_{620nm} data into the amount of phosphate released using the standard curve line-fit data, from above:

$$\text{Phosphate released} = (\text{OD}_{620\text{nm}} - Y_{\text{int}})/\text{slope}$$

EXAMPLE:

Std curve slope=0.3 OD_{620nm}/nmol phosphate

Std curve Y_{int}=0.003 OD_{620nm}

Sample OD_{620nm}=0.4

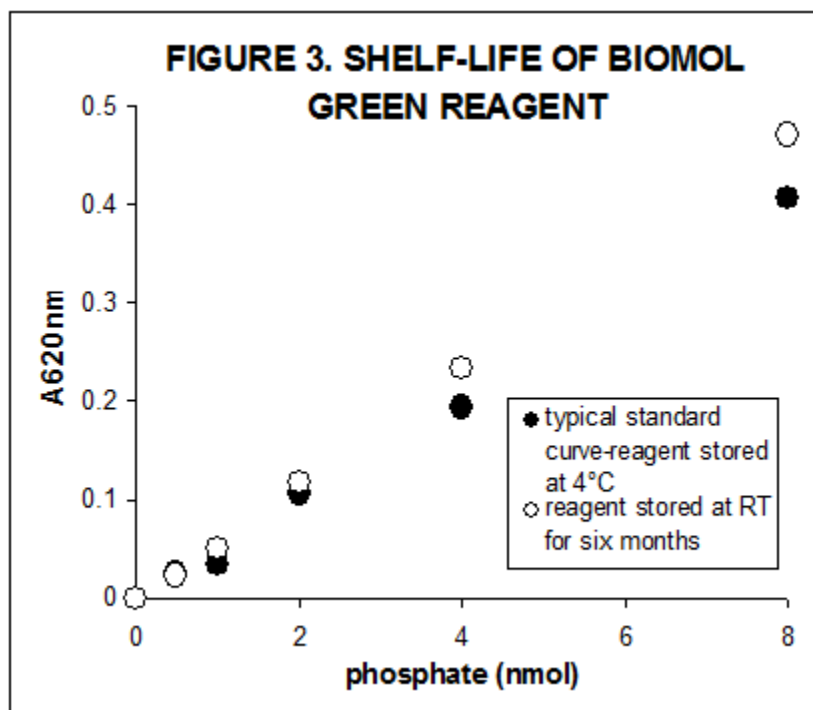
Phosphate released= (0.4-0.003)/0.3 = 1.32 nmol

Other Considerations:

1. The range and background of phosphate standard curves using the BIOMOL® GREEN Reagent may be affected by the composition of assay buffers employed. Therefore, phosphate standard curves should always be performed in a buffer identical to that of the experimental samples. The presence of bovine serum albumin (BSA) in an assay buffer can diminish the signal and/or delay color development. If BSA must be used, a concentration ≤0.1 mg/mL and a 60 min. incubation following addition of BIOMOL® GREEN are recommended. The detergents NP-40 (0.05%) and Brij 35 (0.01%) are compatible with the assay. Tween detergents can suppress color development, especially at lower phosphate concentrations, and are therefore not recommended.
2. Cell extracts or other samples containing extraneous free phosphate and/or interfering substances may be used if desalted prior to analysis. This can be accomplished by dialysis or by gel-filtration (Hint: Use spin-columns).
3. Suitable substrates for phosphatase assays include phosphopeptides, phosphorylated proteins and phosphoamino acids. The substrate must be matched to the target phosphatase of interest. The minimal amount of substrate used in the assay must consider the usable range defined by the phosphate standard curve (e.g., Figure 1 and 2). Typically, <10% of total substrate should be hydrolyzed during the assay to allow linear

kinetics. Thus, a general minimal concentration of substrate is 200 μ M for the microplate (50 μ L assay volume) or cuvette assay (100 μ L assay volume).

REAGENT STABILITY





Product Manual

NOTES



Product Manual



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:

www.enzolifesciences.com