

# **ab133026 – Cyclic GMP ELISA Kit**

## Instructions for Use

For quantitative detection of Cyclic GMP in saliva, plasma (Heparin, EDTA), urine, serum and tissue culture supernatants.

This product is for research use only and is not intended for diagnostic use.

# Table of Contents

---

## INTRODUCTION

- 1. BACKGROUND 2
- 2. ASSAY SUMMARY 3

## GENERAL INFORMATION

- 3. PRECAUTIONS 4
- 4. STORAGE AND STABILITY 5
- 5. MATERIALS SUPPLIED 5
- 6. MATERIALS REQUIRED, NOT SUPPLIED 6
- 7. LIMITATIONS 6
- 8. TECHNICAL HINTS 7

## ASSAY PREPARATION

- 9. REAGENT PREPARATION 8
- 10. STANDARD PREPARATIONS – NON-ACETYLATED FORMAT 9
- 11. STANDARD PREPARATIONS –ACETYLATED FORMAT 11
- 12. SAMPLE COLLECTION AND STORAGE 13
- 13. PLATE PREPARATION 14

## ASSAY PROCEDURE

- 14. ASSAY PROCEDURE 15

## DATA ANALYSIS

- 15. CALCULATIONS 17
- 16. TYPICAL DATA 18
- 17. TYPICAL SAMPLE VALUES 22
- 18. ASSAY SPECIFICITY 25

## RESOURCES

- 19. TROUBLESHOOTING 26
- 20. NOTES 27

## 1. BACKGROUND

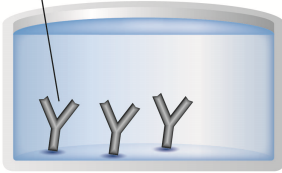
Abcam's Cyclic GMP (cGMP) *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Cyclic GMP in saliva, plasma (Heparin, EDTA), urine, and serum and tissue culture supernatants.

A goat anti-rabbit IgG antibody has been precoated onto 96-well plates. Standards or test samples are added to the wells, along with an alkaline phosphatase (AP) conjugated-cGMP antigen and a polyclonal rabbit antibody specific to cGMP. After incubation the excess reagents are washed away. pNpp substrate is added. After a short incubation the alkaline phosphatase enzyme reaction is stopped and the yellow color generated is read at 405 nm. The intensity of the yellow coloration is inversely proportional to the amount of cGMP captured in the plate.

Guanosine 3', 5'-cyclic monophosphate (cyclic GMP; cGMP) was identified in 1963. It has been shown to be present at levels typically 10-100 fold lower than cAMP in most tissues and is formed by the action of the enzyme guanylate cyclase on GTP. It is involved in a number of important biological reactions. Some hormones, such as acetylcholine, insulin and oxytocin, as well as certain other chemicals like serotonin and histamine, cause an increase in cGMP levels. Stimulators of guanylate cyclase such as the vasodilators nitroprusside, nitroglycerin, sodium nitrate and nitric oxide (NO) also stimulate cGMP levels. Peptides, such as atrial natriuretic peptide (ANP) that relaxes smooth muscle also increase cGMP concentrations. cGMP has been confirmed as a second messenger for ANP. NO can be synthesized from L-arginine and diffuse through cell membranes. The interaction of NO with guanylate cyclase allows cGMP to act as a third messenger in some cells.

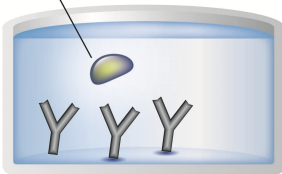
## 2. ASSAY SUMMARY

**Capture Antibody**



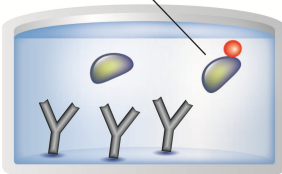
Prepare all reagents and samples as instructed.

**Sample**



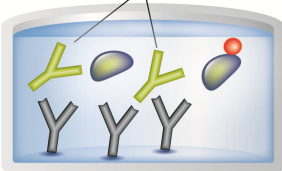
Add standards and samples to appropriate wells.

**Labeled AP-Conjugate**



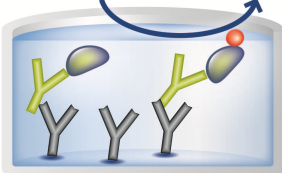
Add prepared labeled AP-conjugate to appropriate wells.

**Target Specific Antibody**



Add Cyclic GMP antibody to appropriate wells. Incubate at room temperature.

**Substrate**      **Colored Product**



Add pNpp substrate to each well. Incubate at room temperature. Add Stop Solution to each well. Read immediately.

### 3. PRECAUTIONS

**Please read these instructions carefully prior to beginning the assay.**

- All kit components have been formulated and quality control tested to function successfully as a kit. Modifications to the kit components or procedures may result in loss of performance
- Some kit components contain azide, which may react with lead or copper plumbing. When disposing of reagents always flush with large volumes of water to prevent azide build-up
- Some Solutions supplied in this kit are caustic; care should be taken with their use
- The activity of the alkaline phosphatase conjugate is dependent on the presence of  $Mg^{2+}$  and  $Zn^{2+}$  ions. The activity of the conjugate is affected by concentrations of chelators (>10 mM) such as EDTA and EGTA
- This kit's performance is tested with a variety of samples, however it is possible that high levels of interfering substances may cause variation in assay results
- The Cyclic GMP Complete Standard provided is supplied in ethanolic buffer at a pH optimized to maintain Cyclic GMP integrity. Care should be taken handling this material because of the known and unknown effects of Cyclic GMP

## 4. STORAGE AND STABILITY

Store kit at +4°C immediately upon receipt, apart from the AP Conjugate & cGMP Standard, which should be stored at -20°C. Avoid multiple freeze-thaw cycles. Components are stable until expiry date.

Refer to list of materials supplied for storage conditions of individual components.

## 5. MATERIALS SUPPLIED

Item	Amount	Storage Condition (Before Preparation)
Goat anti-Rabbit IgG Microplate (12 x 8 wells)	96 Wells	+4°C
Cyclic GMP Complete Alkaline Phosphatase Conjugate	5 mL	-20°C
Cyclic GMP Complete Antibody	5 mL	+4°C
Cyclic GMP Complete Standard	500 µL	-20°C
20X Wash Buffer Concentrate	27 mL	+4°C
pNpp Substrate	20 mL	+4°C
Stop Solution	5 mL	+4°C
Acetylation kit - Triethylamine	2 mL	+4°C
Acetylation kit - Acetic Anhydride	1 mL	+4°C
Assay Buffer 2	27 mL	+4°C

### **6. MATERIALS REQUIRED, NOT SUPPLIED**

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Standard microplate reader - capable of reading at 405 nm, preferably with correction between 570 and 590 nm.
- Automated plate washer (optional)
- Adjustable pipettes and pipette tips. Multichannel pipettes are recommended when large sample sets are being analyzed
- Eppendorf tubes
- Microplate Shaker
- Absorbent paper for blotting
- Triton X-100 (optional for sample preparation)
- Optional (for tissue samples): Liquid nitrogen, mortar & pestle, and concentrated HCl

### **7. LIMITATIONS**

- Assay kit intended for research use only. Not for use in diagnostic procedures
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted

### 8. TECHNICAL HINTS

- Standards can be made up in either glass or plastic tubes.
- Pre-rinse the pipette tip with the reagent, use fresh pipette tips for each sample, standard and reagent.
- Pipette standards and samples to the bottom of the wells.
- Add the reagents to the side of the well to avoid contamination.
- This kit uses break-apart microtiter strips, which allow the user to measure as many samples as desired. Unused wells must be kept desiccated at 4°C in the sealed bag provided. The wells should be used in the frame provided.
- Care must be taken to minimize contamination by endogenous alkaline phosphatase. Contaminating alkaline phosphatase activity, especially in the substrate solution, may lead to high blanks. Care should be taken not to touch pipet tips and other items that are used in the assay with bare hands.
- Prior to addition of substrate, ensure that there is no residual wash buffer in the wells. Any remaining wash buffer may cause variation in assay results.
- **This kit is sold based on number of tests. A ‘test’ simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions**

## 9. REAGENT PREPARATION

Equilibrate all reagents and samples to room temperature (18 - 25°C) prior to use.

### 9.1 **Cyclic GMP Complete Alkaline Phosphatase Conjugate**

Allow the Cyclic GMP Complete Alkaline Phosphatase Conjugate to warm to room temperature. Any unused conjugate should be aliquoted and re-frozen at or below -20°C.

### 9.2 **1X Wash Buffer**

Prepare the 1X Wash Buffer by diluting 5 mL of the 20X Wash Buffer Concentrate in 95 mL of deionized water. Mix thoroughly and gently.

### 9.3 **Acetylation Reagent (optional)**

Prepare the Acetylating Reagent by adding 0.5 mL of Acetic Anhydride to 1 mL of Triethylamine. Note that this volume is sufficient to add to 30 mL of diluted standards and samples. Use the prepared reagent within 60 minutes of preparation.

Discard any unused portion of the Acetylating Reagent.

## 10. STANDARD PREPARATIONS – NON-ACETYLATED FORMAT

Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of standards for every use. Diluted standards should be used within 60 minutes of preparation.

- 10.1 Allow the 5,000 pmol/mL Cyclic GMP **Stock Standard** solution to equilibrate to room temperature. The standard solution should be stored at -20°C. Avoid repeated freeze-thaw cycles. Dilute with either Assay Buffer 2 or Tissue Culture Media as appropriate.
- 10.2 Label six tubes with numbers #1 – #6.
- 10.3 Add 900 µL of appropriate diluent (Assay Buffer or Tissue Culture Media) to tube #1.
- 10.4 Add 800 µL appropriate diluent to tubes #2 through #6.
- 10.5 Prepare a 500 pmol/mL **Standard 1** by adding 100 µL of 5,000 pmol/mL Stock Standard to tube #1. Vortex thoroughly.
- 10.6 Add 800 µL of the appropriate diluent into tubes #2 – #6.
- 10.7 Prepare **Standard 2** by transferring 200 µL from **Standard 1** to tube 2. Mix thoroughly and gently.
- 10.8 Prepare **Standard 3** by transferring 200 µL from **Standard 2** to tube 3. Mix thoroughly and gently.
- 10.9 Using the table below as a guide, repeat for tubes #4 and #6.

## ASSAY PREPARATION

Standard	Sample to Dilute	Volume to Dilute ( $\mu\text{L}$ )	Volume of Diluent ( $\mu\text{L}$ )	Starting Conc. ( $\mu\text{mol/mL}$ )	Final Conc. ( $\mu\text{mol/mL}$ )
1	Standard	100	900	5,000	500
2	Standard 1	200	800	500	100
3	Standard 2	200	800	100	20
4	Standard 3	200	800	20	4
5	Standard 4	200	800	4	0.8
6	Standard 5	200	800	0.8	0.16



## 11. STANDARD PREPARATIONS –ACETYLATED FORMAT (optional)

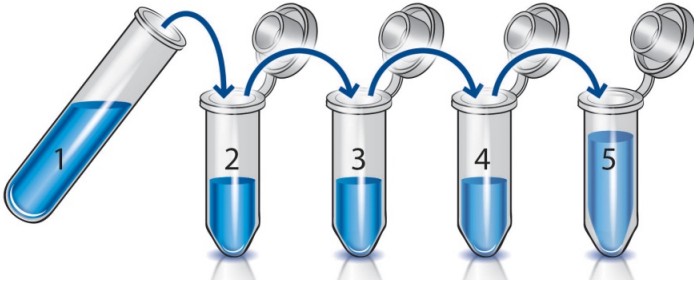
Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of standards for every use Diluted acetylated standards should be used within 30 minutes of preparation.

- 11.1 Allow the 5,000 pmol/mL Cyclic GMP **Stock Standard** solution to equilibrate to room temperature. The standard solution should be stored at -20°C. Avoid repeated freeze-thaw cycles.
- 11.2 Label five tubes with numbers #1 – #5.
- 11.3 Add 980 µL of appropriate diluent (Assay Buffer or Tissue Culture Media) to tube #1.
- 11.4 Add 800 µL appropriate diluent to tubes #2 through #5
- 11.5 Prepare a 500 pmol/mL **Standard 1** by adding 20 µL of 5,000 pmol/mL Stock Standard to tube 1. Vortex thoroughly.
- 11.6 Prepare **Standard 2** by transferring 200 µL from **Standard 1** to tube #2. Vortex thoroughly.
- 11.7 Prepare **Standard 3** by transferring 200 µL from **Standard 2** to tube #3. Vortex thoroughly.
- 11.8 Using the table below as a guide, repeat for tubes #4 through #5.
- 11.9 Acetylate all standards by adding 10 µL of the Acetylating Reagent (see section 9.3) for each 200 µL of the standard. Add the Acetylating Reagent directly to the diluted standard or sample and vortex immediately after the addition of the Acetylating Reagent.

*Note: If acetylating standards, then samples must be acetylated in the same format by adding 20 µL of the Acetylating Reagent for each 100 µL of the sample.*

# ASSAY PREPARATION

Standard	Sample to Dilute	Volume to Dilute ( $\mu\text{L}$ )	Volume of Diluent ( $\mu\text{L}$ )	Starting Conc. ( $\text{pmol/mL}$ )	Final Conc. ( $\text{pmol/mL}$ )
1	Standard	20	980	5,000	100
2	Standard 1	200	800	100	20
3	Standard 2	200	800	20	4.0
4	Standard 3	200	800	4.0	0.8
5	Standard 4	200	800	0.8	0.16



## 12. SAMPLE COLLECTION AND STORAGE

- Store samples to be assayed within 24 hours at 2-8°C. For long-term storage, aliquot and freeze samples at -20°C. Avoid repeated freeze-thaw cycles.
- Samples containing rabbit IgG will interfere with the assay. EDTA plasma may precipitate during acetylation.
- Samples in the majority of tissue culture media, including those containing fetal bovine serum, can also be read in the assay, provided the standards have been diluted into the tissue culture media instead of Assay Buffer 2.
- Urine samples may be used in the assay directly after dilution in Assay Buffer 2. Plasma samples should be drawn in tubes containing EDTA. EDTA chelates calcium and will stop phosphodiesterase activity. The plasma collected should be assayed immediately or frozen below -20°C.
- If samples cannot be diluted, the following extraction method may be used. (Add 2 mL of 95% ethanol to 1 mL of the collected plasma. Vortex for 15 seconds and let sit at room temperature for 5 minutes. Centrifuge for 10 minutes at 600 x g at room temperature. Decant the supernatant into a clean tube. These samples should be dried down, reconstituted in Assay Buffer 2 and then used directly in the assay.)
- Tissue samples should be rapidly frozen in liquid nitrogen. Grind the frozen tissue to a fine powder under liquid nitrogen in a stainless steel mortar. Weigh the frozen tissue and homogenize in 10 volumes of cold 5% TCA in a glass-Teflon tissue grinder. Centrifuge at 600 x g for 10 minutes. Extract the supernatants with 3 volumes of water-saturated ether. Dry the aqueous extracts and run the reconstituted samples directly in the assay.

## 13. PLATE PREPARATION

- The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents
- Unused well strips should be returned to the plate packet and stored at 4°C
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates)
- Well effects have not been observed with this assay. Contents of each well can be recorded on the template sheet included in the Resources section.

	1	2	3	4
A	B <sub>s</sub>	Std 1	Std 5	etc
B	B <sub>s</sub>	Std 1	Std 5	etc
C	TA	Std 2	Std 6	
D	TA	Std 2	Std 6	
E	NSB	Std 3	Sample 1	
F	NSB	Std 3	Sample 1	
G	B <sub>0</sub>	Std 4	Sample 2	
H	B <sub>0</sub>	Std 4	Sample 2	

Plate layout shows controls, blanks and standards required for each assay. Use additional strips of wells to assay all your samples.

### Key:

**B<sub>s</sub>** = Blank; contains substrate only.

**TA** = Total Activity; contains conjugate (5 µL) and substrate.

**NSB** = Non-specific binding; contains standard diluent, assay buffer, conjugate and substrate.

**B<sub>0</sub>** = 0 pg/mL standard; contains standard diluent, conjugate, antibody and substrate

## **14. ASSAY PROCEDURE**

- **Equilibrate all materials and prepared reagents to room temperature prior to use**
- **It is recommended to assay all standards, controls and samples in duplicate**
- **Refer to the recommended plate layout in Section 12 before proceeding with the assay**
- **If Acetylated Version of the kit is to be run, acetylate all standards as described in Section 11. Acetylate all samples by adding 10  $\mu\text{L}$  of the Acetylating Reagent (see step 9.3) for each 200  $\mu\text{L}$  of sample. Add the reagent directly to the samples and vortex for 2 seconds. Use the acetylated samples within 30 minutes.**

14.1 Add 150  $\mu\text{L}$  appropriate diluent\* into the NSB (non-specific binding) wells. (\*Use the same diluent used to prepare standards in section 10, either Assay Buffer or Tissue Culture Media).

14.2 Add 100  $\mu\text{L}$  appropriate diluent (Assay Buffer or tissue culture media) into the  $B_0$  (0 pg/mL standard) wells.

14.3 Add 100  $\mu\text{L}$  of prepared standards and 100  $\mu\text{L}$  diluted samples to appropriate wells.

14.4 Add 50  $\mu\text{L}$  of Cyclic GMP Complete Alkaline Phosphatase Conjugate (blue) into NSB,  $B_0$ , standard and sample wells, i.e. not the Total Activity (TA) and  $B_s$  wells.

14.5 Add 50  $\mu\text{L}$  of Cyclic GMP (yellow) into  $B_0$ , standard and sample wells, i.e. not  $B_s$ , TA and NSB wells.

*Note:* Every well used should be Green in color except the NSB wells which should be Blue. The Blank and TA wells are empty at this point and have no color.

14.6 Incubate the plate at room temperature on a plate shaker for 2 hours at  $\sim 500$  rpm. The plate may be covered with the plate sealer provided.

If using the acetylated format incubate for 18-24 hours at 4°C.

- 14.7 Empty the contents of the wells and wash by adding 400  $\mu$ L of 1X Wash Buffer to every well. Repeat the wash 2 more times for a total of 3 washes. After the final wash, empty or aspirate the wells, and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
- 14.8 Add 5  $\mu$ L of the Cyclic GMP Complete Alkaline Phosphatase Conjugate to the TA wells only.
- 14.9 Add 200  $\mu$ L of the pNpp Substrate solution to every well. Incubate at room temperature for 1 hour without shaking.
- 14.10 Add 50  $\mu$ L Stop Solution into each well. The plate should be read immediately.
- 14.11 Blank the plate reader against the blank wells, read the O.D. absorbance at 405 nm, preferably with correction between 570 and 590 nm. If the plate reader is not able to be blanked against the blank wells, manually subtract the mean optical density of the blank wells from all readings.

## 15. CALCULATIONS

- 15.1 Calculate the average net absorbance measurement (Average Net OD) for each standard and sample by subtracting the average NSB absorbance measurement from the average absorbance measurement (Average OD) for each standard and sample.

$$\text{Average Net OD} = \text{Average Bound OD} - \text{Average NSB OD}$$

- 15.2 Calculate the binding of each pair of standard wells as a percentage of the maximum binding wells ( $B_0$ ), using the following formula

$$\text{Percent Bound} = \frac{\text{Average Net OD}}{\text{Average Net } B_0 \text{ OD}} \times 100$$

- 15.3 Plot the Percent Bound ( $B/B_0$ ) and the net OD versus concentration of Cyclic GMP for the standards. The concentration of Cyclic GMP in the unknowns can be determined by interpolation of net OD values.

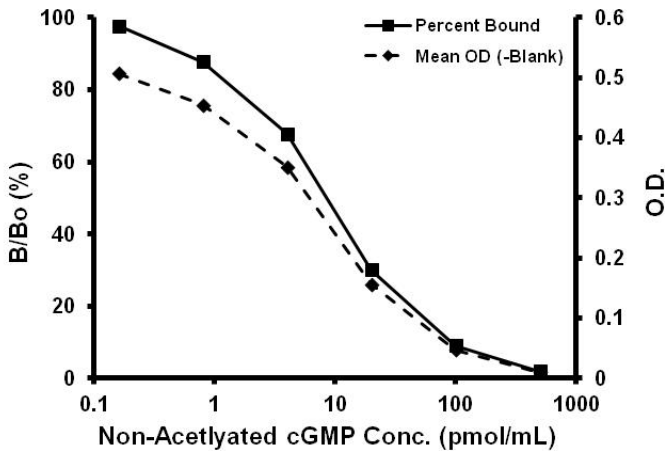
A four parameter algorithm (4PL) provides the best fit, though other equations can be examined to see which provides the most accurate (e.g. linear, semi-log, log/log, 4 parameter logistic). Interpolate protein concentrations for unknown samples from the standard curve plotted.

Samples producing signals greater than that of the highest standard should be further diluted and reanalyzed, then multiplying the concentration found by the appropriate dilution factor

## 16. TYPICAL DATA

**NOTE** – This data is provided for **demonstration purposes only**. A new standard curve must be generated for each assay performed.

**NON-ACETYLATED:**



Sample	Average Net OD	Percent Bound	cGMP (pmol/mL)
B <sub>s</sub>	(0.176)		
TA	0.296		
NSB	0	0	
Standard 1	0.01	1.9	500
Standard 2	0.047	9.1	100
Standard 3	0.156	30.1	20
Standard 4	0.352	67.8	4
Standard 5	0.455	87.7	0.8
Standard 6	0.507	97.7	0.16
B <sub>0</sub>	0.519	100	0
Unknown 1	0.3	58.9	6.05
Unknown 2	0.098	18.9	39.3

## Typical Quality Control Parameters – Non-Acetylated

Total Activity Added =  $0.296 \times 10 = 2.96$

%B<sub>0</sub>/TA = 17.6%

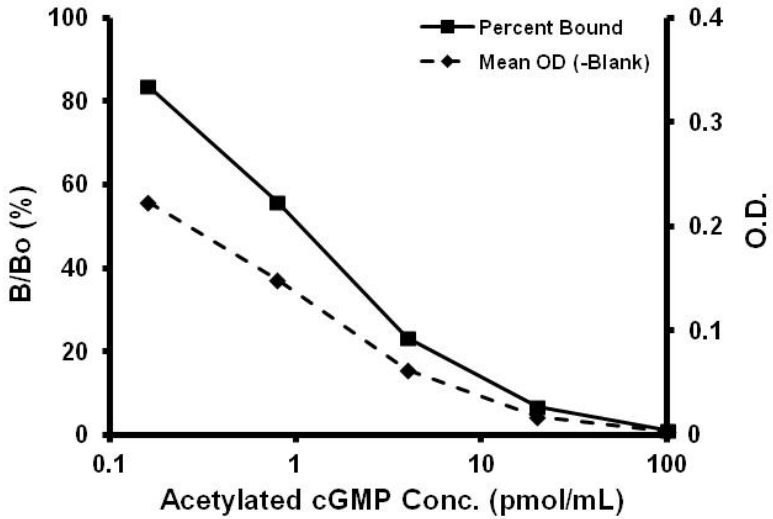
Quality of Fit = 0.9999

20% Intercept = 36.5 pmol/mL

50% Intercept = 8.4 pmol/mL

80% Intercept = 1.9 pmol/mL

## ACETYLATED:



Sample	Average Net OD	Percent Bound	cGMP (pmol/mL)
B <sub>s</sub>	(0.146)		
TA	0.32		
NSB	0.002	0	
Standard 1	0.003	1.1	100
Standard 2	0.018	6.8	20
Standard 3	0.062	23.3	4
Standard 4	0.149	56	0.8
Standard 5	0.223	83.8	0.16
B <sub>0</sub>	0.266	100	0
Unknown 1	0.13	48.9	1.12
Unknown 2	0.01	3.8	33.9

## Typical Quality Control Parameters

Total Activity Added =  $0.320 \times 10 = 3.20$

%B<sub>0</sub>/TA = 8.3%

Quality of Fit = 1.0000

20% Intercept = 5.1 pmol/mL

50% Intercept = 1.1 pmol/mL

80% Intercept = 0.2 pmol/mL

**17. TYPICAL SAMPLE VALUES**

**SENSITIVITY –**

The sensitivity, minimum detectable dose of Cyclic GMP using this Abcam ELISA kit was measured at 2 standard deviations from the mean of 16 zeros along the standard curve, and was determined to be 0.37 pmol/mL in the non-acetylated assay format and 0.088 pmol/mL in the acetylated assay format.

**SAMPLE RECOVERY –**

**Non-Acetylated Format:**

Sample Type	Average % Recovery	Recommended Dilution
Tissue Culture Media	101.7	1:100
Human Serum	102.9	1:10
Human Saliva	101.3	1:10-1:100
Human Heparin Plasma	104.4	1:10
Human EDTA Plasma	115.0	1:10-1:100
Human Urine	97.7	1:100-1:1000

**Acetylated Format:**

Sample Type	Average % Recovery	Recommended Dilution
Tissue Culture Media	95.8	Undiluted
Human EDTA Plasma	93.6	≥ 1:2

### **LINEARITY OF DILUTION –**

#### **Non-Acetylated Version**

A buffer sample containing 800 pmol/mL cGMP was serially diluted 8 times 1:2 in the kit Assay Buffer 2 and measured in the assay. The data was plotted graphically as actual cGMP concentration versus measured cGMP concentration.

The line obtained had a slope of 1.060 with a correlation coefficient of 0.998.

#### **Acetylated Version**

A buffer sample containing 5.79 pmol/mL cGMP was serially diluted 3 times 1:2 in the kit Assay Buffer 2 and measured in the Acetylated version of the assay. The data was plotted graphically as actual cGMP concentration versus measured cGMP concentration.

The line obtained had a slope of 0.903 with a correlation coefficient of 0.999.

## PRECISION –

Intra-assay precision was determined by taking samples containing low, medium and high concentrations of Cyclic GMP and running these samples multiple times (n=16) in the same assay.

Inter-assay precision was determined by measuring three samples with low, medium and high concentrations of Cyclic GMP in multiple assays (n=16).

The precision numbers listed below represent the percent coefficient of variation for the concentrations of cGMP determined in these assays as calculated by a 4 parameter logistic curve fitting program.

### Non-Acetylated Format:

#### Intra-Assay

	pmol/mL	%CV
Low	1.5	5.2
Medium	16.6	4.0
High	481	7.6

#### Inter-Assay

	pmol/mL	%CV
Low	1.8	13.7
Medium	16.9	3.5
High	359	5.0

## Acetylated Format:

### Intra-Assay

	<b>pmol/mL</b>	<b>%CV</b>
Low	0.54	6.5
Medium	1.50	4.6
High	6.81	4.5

### Inter-Assay

	<b>pmol/mL</b>	<b>%CV</b>
Low	0.70	5.9
Medium	1.96	6.2
High	8.61	6.8

## 18. ASSAY SPECIFICITY

### CROSS REACTIVITY –

The cross reaction of the antibody calculated at 50% is:

Cyclic GMP	100 %
GMP	<0.001 %
GTP	<0.001 %
cAMP	<0.001 %
AMP	<0.001 %
ATP	<0.001 %
cUMP	<0.001 %
CTP	<0.001 %

## 19. TROUBLESHOOTING

<b>Problem</b>	<b>Cause</b>	<b>Solution</b>
Poor standard curve	Inaccurate pipetting	Check pipettes
	Improper standards dilution	Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing
Low Signal	Incubation times too brief	Ensure sufficient incubation times; change to overnight standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
Samples give higher value than the highest standard	Starting sample concentration is too high.	Dilute the specimens and repeat the assay
Large CV	Plate is insufficiently washed	Review manual for proper wash technique. If using a plate washer, check all ports for obstructions
	Contaminated wash buffer	Prepare fresh wash buffer
Low sensitivity	Improper storage of the kit	Store the all components as directed.

20. NOTES







**UK, EU and ROW**

Email: [technical@abcam.com](mailto:technical@abcam.com) | Tel: +44-(0)1223-696000

**Austria**

Email: [wissenschaftlicherdienst@abcam.com](mailto:wissenschaftlicherdienst@abcam.com) | Tel: 019-288-259

**France**

Email: [supportscientifique@abcam.com](mailto:supportscientifique@abcam.com) | Tel: 01-46-94-62-96

**Germany**

Email: [wissenschaftlicherdienst@abcam.com](mailto:wissenschaftlicherdienst@abcam.com) | Tel: 030-896-779-154

**Spain**

Email: [soportecientifico@abcam.com](mailto:soportecientifico@abcam.com) | Tel: 911-146-554

**Switzerland**

Email: [technical@abcam.com](mailto:technical@abcam.com)

Tel (Deutsch): 0435-016-424 | Tel (Français): 0615-000-530

**US and Latin America**

Email: [us.technical@abcam.com](mailto:us.technical@abcam.com) | Tel: 888-77-ABCAM (22226)

**Canada**

Email: [ca.technical@abcam.com](mailto:ca.technical@abcam.com) | Tel: 877-749-8807

**China and Asia Pacific**

Email: [hk.technical@abcam.com](mailto:hk.technical@abcam.com) | Tel: 400 921 0189 / +86 21 2070 0500

**Japan**

Email: [technical@abcam.co.jp](mailto:technical@abcam.co.jp) | Tel: +81-(0)3-6231-0940

[www.abcam.com](http://www.abcam.com) | [www.abcam.cn](http://www.abcam.cn) | [www.abcam.co.jp](http://www.abcam.co.jp)