

Calcineurin Phosphatase Assay Kit

Catalog #: BML-AK804

A complete colorimetric assay kit for measuring calcineurin phosphatase activity

For the latest product information, including support documentation, visit us online:

www.enzolifesciences.com

Technical Support (US): 800-942-0430
Technical Support (EU): +41 61 926 8989



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 use only. Not for use in diagnostic procedures. 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 PROCEDURES.**

TABLE OF CONTENTS



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

BACKGROUND.....	2
MATERIALS SUPPLIED.....	3
STORAGE.....	4
ADDITIONAL MATERIALS NEEDED.....	4
SAFETY WARNINGS & PRECAUTIONS.....	5
PROCEDURE.....	5
PREPARING REAGENTS FOR ASSAY.....	5
PERFORMING THE ASSAY.....	5
DATA ANALYSIS.....	8
TROUBLESHOOTING.....	11
SELECTED REFERENCES FOR INHIBITORS SCREENED WITH ENZO'S CALCINEURIN KIT.....	12
REFERENCES.....	13

BACKGROUND

Calcineurin (CaN) is the neuronal form of the widely distributed Ca^{2+} /calmodulin-dependent Ser/Thr protein phosphatase 2B (PP-2B). CaN is a heterodimer consisting of a catalytic A subunit (57-61 kDa) and a regulatory B subunit (19 kDa). The catalytic A subunit is composed of four functional domains: the catalytic core with sequence homology to PP-1 and PP-2A (located between residues 71-235 in the rat brain $\alpha\delta$ isoform), binding sites for both calmodulin (residues 391-414) and CaN B-regulatory subunit, and a C-terminal (residues 457-482) autoinhibitory domain.

The *Calcineurin Phosphatase Assay Kit* is a complete colorimetric assay kit for measuring calcineurin phosphatase activity. It employs a convenient 96-well microtiter-plate format with all reagents necessary for measuring calcineurin (PP2B) phosphatase activity of purified enzyme. The RII phosphopeptide substrate, supplied with this kit, is the most efficient and outstanding peptide substrate known for calcineurin^{1,2}. The detection of free-phosphate released is based on the classic Malachite green assay^{3,4} and offers the following advantages:

- NON-RADIOACTIVE!
- CONVENIENT 1-STEP DETECTION -no mixing!
- MICROTITER-PLATE FORMAT -for high-throughput!

This new, improved version of the BML-AK804 kit incorporates human calcineurin A α (MW=60 kDa) + calcineurin B (MW=15 kDa) heterodimer expressed in an *E. coli* expression system. Both subunits were coexpressed in a construct with yeast myristoyl-CoA:protein N-myristoyltransferase. The resulting highly active calcineurin (protein phosphatase 2B) is N-myristoylated on the CaNB subunit, similar to the native protein⁵.

References

1. A. Enz *et al. Anal. Biochem.* 1994 **216** 147
2. A. Donella-Deana *et al. Eur. J. Biochem.* 1994 **219** 109
3. B. Martin *et al. J. Biol. Chem.* 1985 **260** 14932
4. K.W. Harder *et al. Biochem. J.* 1994 **298** 395
5. A. Mondragon *et al. Biochemistry* 1997 **36** 4934

MATERIALS SUPPLIED

BML-SE163-5000 CALCINEURIN ENZYME (human, recombinant)

FORM: 100 U/μl in 1X assay buffer (1:1 dilution of BML-KI128, below).

1 U=pmol/min @ 30°C.

STORAGE: -80°C; AVOID FREEZE/THAW CYCLES!

QUANTITY: 5000 U

BML-SE325-9090 CALMODULIN (human, recombinant)

FORM: 25 μM in dH₂O

STORAGE: -80°C

QUANTITY: 100μl

BML-P160-9090 SUBSTRATE (R11 phosphopeptide, sequence Asp-Leu-Asp-Val-Pro-Ile-Pro-Gly-Arg-Phe-Asp-Arg-Arg-Val-pSer-Val-Ala-Ala-Glu; MW=2192.0)

FORM: 1.5 mg net peptide/vial

STORAGE: -20°C

QUANTITY: 1 x 1.5 mg

BML-KI128-0020 2X ASSAY BUFFER

(100 mM Tris, pH 7.5, 200 mM NaCl, 12 mM MgCl₂, 1 mM DTT, 0.05% NP-40, 1 mM CaCl₂)

FORM: Liquid in screw-cap plastic bottle.

STORAGE: -80°C

QUANTITY: 20 ml

BML-AK111-9090 BIOMOL[®] Green REAGENT

FORM: Liquid in screw-cap plastic bottle.

STORAGE: +4°C

QUANTITY: 20 ml

BML-KI132-0500 PHOSPHATE STANDARD

FORM: 80 μM in dH₂O

STORAGE: -80°C

QUANTITY: 0.5 ml

80-2404 1/2 VOLUME MICROPLATE

1 clear, 96-well

STORAGE: Ambient



Storage temp

STORAGE

Please note that all components, with the exception of Calcineurin (BML-SE163-5000) and Calmodulin (BML-SE325-9090) can be stored at the temperatures listed or at -80°C . The enzymes listed above must be stored at -80°C . *The calcineurin enzyme component BML-SE163 must be handled particularly carefully in order to retain maximal enzymatic activity. Thaw it quickly in a RT water bath or by rubbing between fingers, then immediately store on an ice bath. The remaining unused enzyme should be quickly refrozen by placing at -80°C . To minimize the number of freeze/thaw cycles, aliquot the calcineurin into separate tubes and store at -80°C .*

ADDITIONAL MATERIALS NEEDED

- Microplate reader capable of measuring A620 to ≥ 3 -decimal accuracy.
- Pipette capable of pipetting 5-100 μl accurately
- Multi-channel pipette capable of pipetting 100 μl (optional).
- Ice bucket to keep reagents cold until use.



Important/ Warning

SAFETY WARNINGS & PRECAUTIONS

1. Wear appropriate personnel protective apparel. Avoid contact with clothes and exposed skin. In case of accidental skin exposure, flush with water immediately. Consult a physician if required.
2. Use a safety pipetting device for all pipetting. Never pipet by mouth.
3. Interpretation of the results is the sole responsibility of the user.

PROCEDURE

NOTE ON HANDLING: *Hold all samples on ice bath until use, unless otherwise noted.*

PRECAUTIONS: *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 dH₂O or employ unused plasticware.*

PREPARING REAGENTS FOR ASSAY

1. Thaw all kit components and hold calcineurin, calmodulin and 2x assay buffer on an ice bath; equilibrate BIOMOL® Green reagent to room temperature.
2. Add calmodulin to the 2x assay buffer: Dilute calmodulin (BML-SE101-9090) 1/50 in 2X assay buffer (BML-KI128) to required quantity (25µl are required per assay well). For example, add 20µl to 980µl 2X assay buffer.
3. Reconstitute substrate (R11 phosphopeptide, BML-P160) with dH₂O to 0.75 mM (1.64 mg/ml): Add 915µl dH₂O per 1.5 mg vial (10µl are needed per assay well).

PERFORMING THE ASSAY

To prepare a standard curve:

4. Prepare 1 ml of 1X assay buffer (dilute 500µl of 2X assay buffer with 500µl dH₂O).
5. Perform 1:1 serial dilutions of phosphate standard and an assay buffer blank. Concentrations of 40, 20, 10, 5, 2.5, 1.25 and 0.625µM correspond to 2, 1, 0.5, 0.25, 0.125, 0.063 and 0.031 nmol PO₄ (see Table 1):

- a. Add 50 μ l of 2X assay buffer (BML-KI128) to each wells A1, and A2 (2 nmol PO₄ standards).
- b. Add 50 μ l 1X assay buffer (prepared in step 4 above) to wells B1-H1 and wells B2-H2 (remaining standard concentrations)
- c. Add 50 μ l of 80 μ M phosphate standard to well A1 and A2 of assay plate. Mix thoroughly by pipetting up and down several times.
- d. Remove 50 μ l from well A1 and add it to well B1. Mix thoroughly by pipetting up and down several times.
- e. Remove 50 μ l from well B1 and add it to well C1.
- f. Mix thoroughly and repeat for wells D1-G1. At well G1, remove 50 μ l and discard. DO NOT PROCEED TO WELL H1 (assay buffer blank). Final volume=50 μ l.
- g. Repeat serial dilution for the wells in column 2 (standard curve duplicates)

To prepare a time course/linearity assay:

6. Add 25 μ l 2X assay buffer (BML-KI128) w/ calmodulin (step 2) to microtiter plate wells designated for linearity assay (see Table 1).
7. Dilute the calcineurin (BML-SE163) to 8 U/ μ l, in 1X assay buffer, and add 5 μ l diluted calcineurin to wells. Final amount of calcineurin= 40 U per well.
8. Add 10 μ l dH₂O to each well.
9. Designate a reaction time to each well (e.g.: 60 min, 40 min, 30 min, 20 min, 10 min, 5 min, 2 min, 0 min).
10. Equilibrate microtiter plate to reaction temperature (e.g.: 30°C).
11. Start reaction by addition of 10 μ l phosphopeptide substrate (BML-P160; 0.75 mM from step 3) at appropriate time point. Make the addition in the reverse time order such that all incubations end at the same time (e.g.: Add 60 min time pt. at t=0; add 5 min at t=55 min, etc.). Final substrate concentration= 0.15 mM.

To prepare a test sample/inhibition assay:

12. Add 25 μ l assay buffer (BML-KI128) w/ calmodulin (step 2) to wells in microtiter plate. See Table 1.
13. Add 5 μ l diluted calcineurin (BML-SE163) to wells (step 7). Final amount of calcineurin= 40 U per well.
14. Add 10 μ l dH₂O to control wells.
15. Add 10 μ l of test sample/inhibitor (dissolved in dH₂O) to test sample wells.
16. Allow test sample/inhibitor to equilibrate to reaction temperature (e.g.: 30°C) for 10 minutes.
17. Start reaction by addition of 10 μ l phosphopeptide substrate (BML-P160; 0.75 mM from step 3). Final concentration= 0.15

mM. Allow reaction to proceed for a time period in which the reaction is linear (~10 min, see below).

To terminate reactions:

18. After incubating wells for desired duration, terminate reactions with 100µl BIOMOL® green reagent (BML-AK111-9090).
19. Allow color to develop 20-30 minutes, making sure all wells spend approximately the same time with the reagent before reading on microplate reader.
20. Read OD_{620nm} on microplate reader.
21. Perform data analysis (see below).

NOTE: Retain microplate for future use of unused wells!

Table 1. Example of Microtiter Plate Samples

Sample†	Std Curve	Time course	Test Samples
Well #	1,2	3,4	5,6
A	2 nmol PO ₄	60 min	Control
B	1	40	Inhibitor/test sample
C	0.5	30	
D	0.25	20	
E	0.125	10	
F	0.063	5	
G	0.031	2	
H	0	0	

† For highest accuracy, perform all samples in duplicate.

Table 2. Typical Assay Components

	2X Assay Buffer w/CaM	Calcineurin (40U)	H ₂ O	Test compound	Substrate (0.75 mM)
CONTROL	25µl	5µl	10µl	0	10µl
TEST SAMPLE	25µl	5µl	0	10µl	10µl

DATA ANALYSIS

Phosphate (PO₄) Standard Curve

1. Plot standard curve data as OD_{620nm} versus nmol PO₄ (Note that a background OD_{620nm} value for 0 nmol PO₄ has been subtracted from all data. See Figure 1. Data may also be plotted without subtracting the background. In that case, however, one should also not subtract background from experimental OD_{620nm} values before using the standard curve to convert them to nmol of PO₄).
2. Obtain a line-fit to the data using an appropriate routine.
3. Use the slope and Y-intercept to calculate amount of phosphate released for other experimental data (e.g., time course and experimental data).

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.*

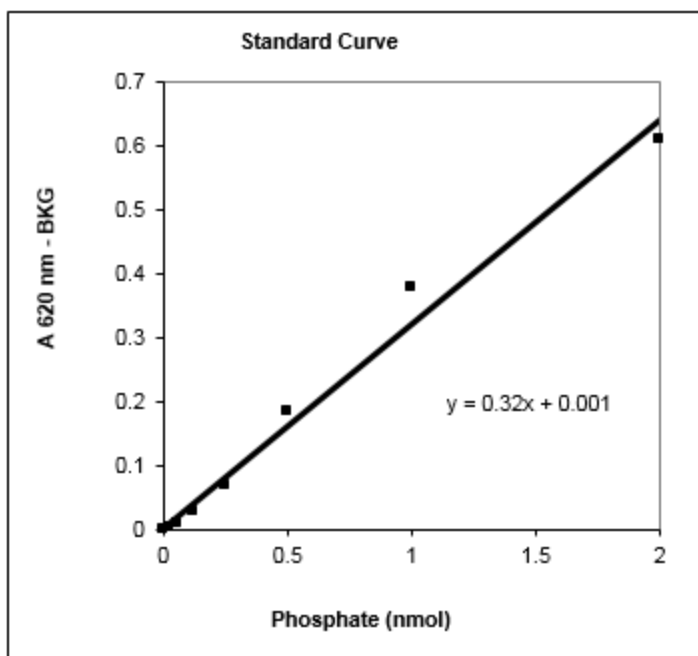


Figure 1. BIOMOL[®] Green Phosphate Standard Curve

Conversion of OD_{620nm} to Amount Phosphate Released

1. 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} = 0.001 OD_{620nm}

Sample OD_{620nm} = 0.4

$$\text{Phosphate released} = (0.4 - 0.001) / 0.3 = 1.33 \text{ nmol}$$

Time Course/Linearity Curve

1. If the 0 time (Table 1, wells H3,4) has a significant value, subtract this number from all samples. This is background phosphate in the samples.
2. Plot OD_{620nm} versus reaction time. See Figure 2. Alternatively, the OD_{620nm} can be converted to phosphate released, as above.
3. Determine the reaction time range in which the amount of phosphate released is linear. In Figure 2, this range is from 0-60 min. This value is variable depending on reaction conditions and storage/handling of the calcineurin. The time range can be lengthened by decreasing the amount of calcineurin in the assay and lowering the assay temperature. For accurate results, it is important to perform inhibitor/agonist assays under linear assay conditions.

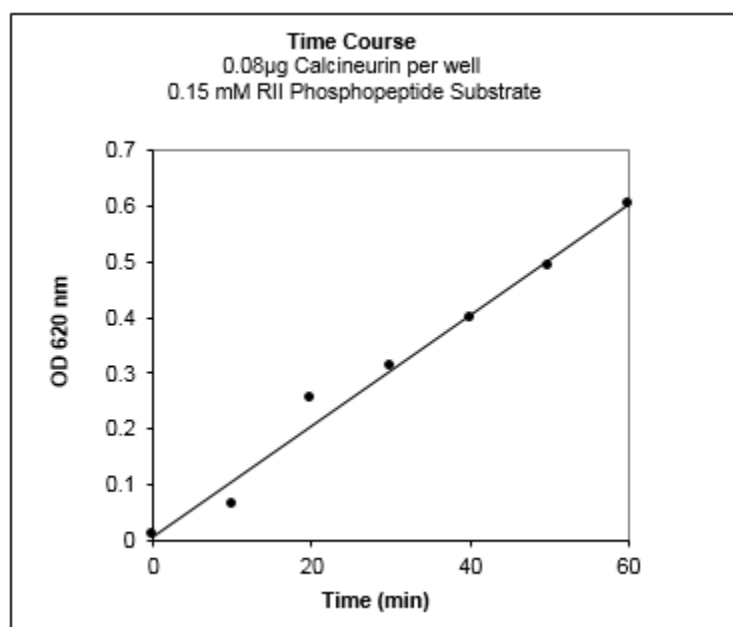


Figure 2. Time Course of Phosphate Released by Calcineurin

Calmodulin Activation of Calcineurin Activity

1. Figure 3 illustrates the activation of calcineurin's phosphatase activity by calmodulin. In the presence (+) of calmodulin, calcineurin's activity is high. In the absence (-) of calmodulin, calcineurin activity is relatively low.

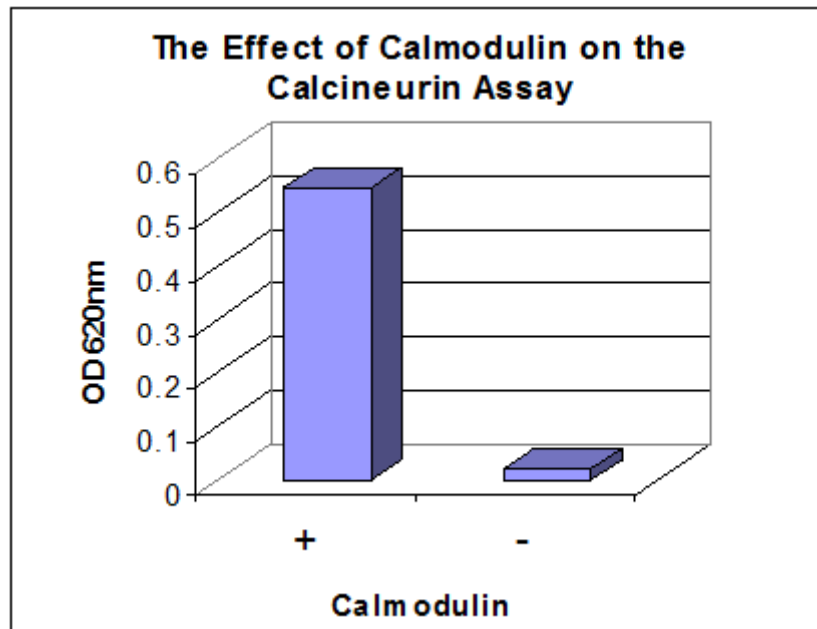


Figure 3. Calmodulin Activation of Calcineurin Phosphatase Activity

TROUBLESHOOTING

Problem	Potential Cause	Solution
High background (High signal with no added calcineurin)	<p>Interfering substance present.</p> <p>Phosphatase contamination of substrate stock solution.</p>	<p>Remove trace amounts of phosphate from assay buffers/reagents.</p> <p>Only use 18 MΩ deionized water, such as Milli-Q™ water, in preparation of all buffers.</p> <p>Soaps and detergents may cause high background. Any container coming into contact with any solutions used in the assay should be triple washed with deionized water prior to use.</p> <p><i>Milli-Q is a trademark of Millipore Corporation.</i></p>
Expected calcineurin activity is not detected	<p>Calcineurin has been inactivated.</p> <p>Activators are not present.</p>	<p>Use positive control phosphatase or standard curve to check assay performance.</p> <p>Be sure necessary co-factors, such as calcium and calmodulin, are in reaction mix.</p>
Cloudy precipitate	Interfering substances present.	Identify and remove incompatible metals, phosphate or detergents. Certain divalent cations (Magnesium, Copper, Zinc) and detergents (SDS and deoxycholate) should be avoided.
Weak signal	Calcineurin is too dilute.	Increase amount of calcineurin used or increase the assay time.
Tested calcineurin inhibitor fails to demonstrate expected activity	Inhibitor concentration employed in assay was too low.	Perform assay using a broader range of inhibitor concentrations. Verify inhibition with other well-characterized inhibitors (e.g. Cyclosporin A, FK506, Cypermethrin, Deltamethrin, or Fenvalerate)

SELECTED REFERENCES FOR INHIBITORS SCREENED WITH ENZO'S CALCINEURIN KIT

	Inhibitor	Reference
1	FK506 Inhibition of CN activity	Toxicological Sciences (2007) 100: 474–485
2	Targeted inhibition of calcineurin with cain or an adenovirus expressing calcineurin inhibitory domain	PNAS (2000) 97: 1196–1201
3	RCAN1 inhibition of calcineurin	Arch. Biochem and Biophys (2007) 467: 185–192
4	cyclosporin A, FK506	Pharmacology, Biochemistry and Behavior (2008) 90: 763–768
5	Minimal RCAN derived sequence (part of the RCAN CIC motif) that inhibits Cn-NFAT signaling in vivo.	J. Biol Chem (2009) 284: 9394–9401
6	Inhibition of calcineurin by caffeoyl phenylethanoid glycosides	J. of Ethnopharmacology (2011) 137: 1306– 1310
7	Jasminum humile leaf and root ethanolic extracts inhibit CN-dependent gene expression in a yeast model system	J. of Ethnopharmacology (2012) 140: 293– 297
8	Effects of CsA and FK506 on aquaporin-2	Pflugers Arch - Eur J Physiol (2011) 462: 611–622
9	Multiple Domains of MCIP1 Contribute to Inhibition of Calcineurin Activity	J Biol Chem. (2002) 277: 30401-7.
10	Tacrolimus (TRL) calcium-dependent protein phosphatase calcineurin CN	Analytical Biochemistry (2006) 358: 104–110

REFERENCES

1. B. Mehul *et al.* *J. Biol. Chem.* 2000 **275** 12841
2. T. Taigen *et al.* *Proc. Natl. Acad. Sci.* 2000 **97** 1196
3. G. Mallert *et al.* *Cell* 2001 **104** 675
4. M. Ichida and T. Finkel *J. Biol. Chem.* 2001 **276** 3524
5. O. Bueno *et al.* *Proc. Natl. Acad. Sci.* 2002 **99** 4586



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