

Calcineurin Phosphatase Activity Assay Kit (Colorimetric) ab139461

10 References 3 Images

Overview

Product name	Calcineurin Phosphatase Activity Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Purified protein, Inhibitor compounds
Assay type	Enzyme activity
Product overview	Abcam Calcineurin Phosphatase Activity Assay Kit (Colorimetric) (ab139461) is a complete non-radioactive colorimetric assay kit designed for measuring calcineurin phosphatase activity through kinetic assay or for inhibitor screening. The activity is determined using the RII phosphopeptide substrate, the most efficient and selective peptide known for calcineurin, and detection of free phosphate released is based on the classic malachite green assay.
Notes	Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.
Platform	Microplate reader

Properties

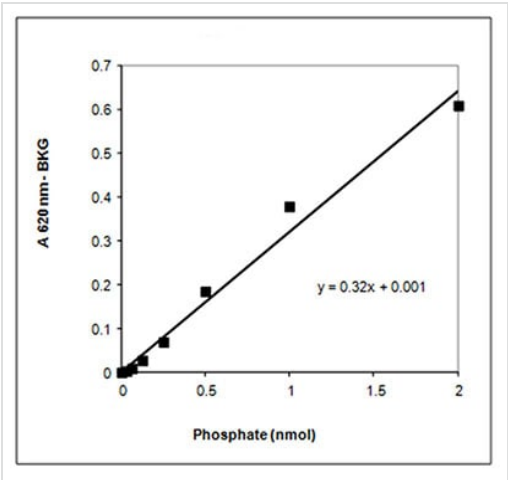
Storage instructions Store at -80°C. Please refer to protocols.

Components	1 x 96 tests
96-well Clear Microplate (1/2 Volume)	1 unit
Active Calcineurin Enzyme	1 x 50µl
Calcineurin Assay Buffer	1 x 20ml
Calcineurin Substrate	1 x 1.5mg
Calmodulin	1 x 100µl

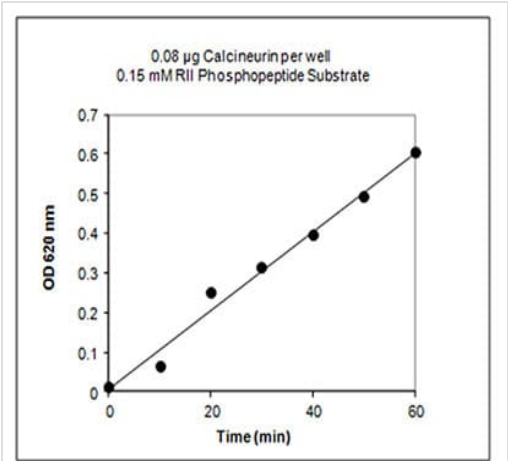
Components	1 x 96 tests
Green Assay Reagent	1 x 20ml
Phosphate Standard	1 x 0.5ml

Function	Calcium-dependent, calmodulin-stimulated protein phosphatase. This subunit may have a role in the calmodulin activation of calcineurin. Dephosphorylates DNM1L, HSPB1 and SSH1.
Sequence similarities	Belongs to the PPP phosphatase family. PP-2B subfamily.
Cellular localization	Nucleus. Colocalizes with ACTN1 and MYOZ2 at the Z line in heart and skeletal muscle.

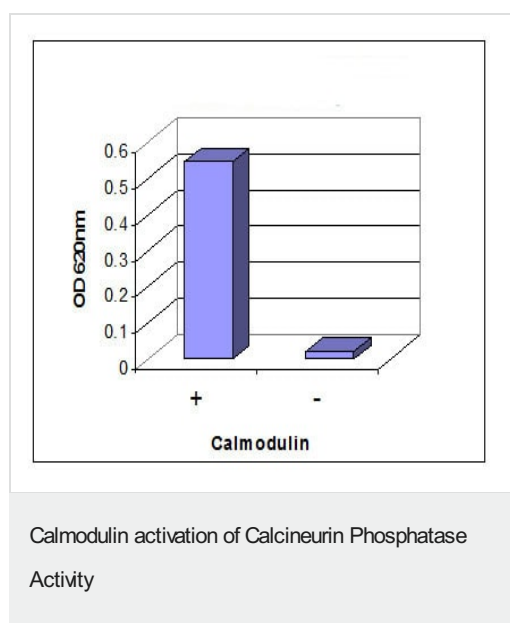
Images



Standard Curve



Time Course



In the presence (+) of Calmodulin, Calcineurin's activity is high. In the absence (-) of Calmodulin, Calcineurin activity is relatively low.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors