

## Description

The HIPK3 Kinase Assay Kit is designed to measure HIPK3 kinase activity for screening and profiling applications using Kinase-Glo<sup>®</sup> MAX as a detection reagent. The assay kit comes in a convenient 96-well format, with enough purified recombinant HIPK3 kinase, kinase substrate, ATP, and kinase assay buffer for 100 enzyme reactions.

## Background

HIPK3 (Homeodomain Interacting Protein Kinase 3) is a serine/threonine protein kinase involved in the regulation of gene transcription induced by steroid hormones such as testosterone. Also known as ANPK (Androgen receptor-interacting nuclear protein kinase), it amplifies androgen receptor-mediated transcription. The kinase phosphorylates transcription factors JUN and RNUX2 and may decrease apoptosis through phosphorylation of FADD (FAS-associated death domain protein).

## Applications

Study enzyme kinetics and screen small molecular inhibitors for drug discovery and high throughput (HTS) applications.

## Supplied Materials

Catalog #	Name	Amount	Storage
40110	HIPK3*	8 µg	-80°C
79334	Kinase assay buffer 1 (5x)	1.5 ml	-20°C
79686	ATP (500 µM)	100 µl	-20°C
40535	Myelin basic protein (MBP), 5 mg/ml	100 µl	-20°C
79696	White 96-well plate	1	Room Temperature

\*The concentration of the protein is lot-specific and will be indicated on the tube.

## Materials Required but Not Supplied

Name	Catalog #
Kinase-Glo MAX	Promega #V6071
DTT (Dithiothreitol), 1 M, optional	
Microplate reader capable of reading luminescence	
Adjustable micropipettor and sterile tips	
30°C incubator	

## Storage Conditions



This assay kit will perform optimally for up to **6 months** from date of receipt when the materials are stored as directed.

**Safety**

This product is for research purposes only and not for human or therapeutic use. This product should be considered hazardous and is harmful by inhalation, in contact with skin, eyes, clothing, and if swallowed. If contact occurs, wash thoroughly.

**Assay Principle**

Kinase activity is measured using Kinase-Glo™ Max (Promega, #V6071). The addition of the reagent results in the generation of a luminescent signal that correlates with the amount of ATP. The reagent is linear to 100µM ATP.

**Contraindications**

The final concentration of DMSO in the assay should not exceed 1%.

**Assay Protocol**

All samples and controls should be tested in duplicate.

1. Thaw **5x Kinase assay buffer**, **ATP**, and **Myelin basic protein (MBP), 5 mg/ml substrate**.  
Optional: If desired, add DTT to **5x Kinase assay buffer** to make a 10 mM DTT concentration (for example, add 10 µl of 1 M DTT to 1 ml of **5x Kinase assay buffer**).
2. Prepare 3 ml of **1x Kinase assay buffer** by mixing 600 µl of **5x Kinase assay buffer** with 2400 µl water. Three (3) ml of **1x Kinase assay buffer** is sufficient for 100 reactions.
3. Prepare the Test Inhibitor (5 µl/well): for a titration, prepare serial dilutions at concentrations 10-fold higher than the desired final concentrations. The final volume of the reaction is 50 µl.
  - a. If the Test Inhibitor is water-soluble, prepare serial dilutions in the 1x Kinase Assay Buffer, 10-fold more concentrated than the desired final concentrations. For the positive and negative controls, use 1x Kinase Assay Buffer (Diluent Solution).
  - b. If the Test inhibitor is soluble in DMSO, prepare the test inhibitor at 100-fold the highest desired concentration in DMSO, then dilute the inhibitor 10-fold in 1x Kinase Assay Buffer to prepare the highest concentration of the 10-fold intermediate dilutions. The concentration of DMSO is now 10%.  
Prepare serial dilutions of the Test Inhibitor at 10-fold the desired final concentrations using 10% DMSO in 1x Kinase Assay Buffer to keep the concentration of DMSO constant.

For positive and negative controls, prepare 10% DMSO in water (vol/vol) so that all wells contain the same amount of DMSO (Diluent Solution).

The final concentration of DMSO should not exceed 1%.

4. Add 5 µl of Test Inhibitor to each well labeled "Test Inhibitor." For the "Positive Control" and "Blank," add Diluent Solution (either distilled water or 10% DMSO in water, as described above).
5. To the wells designated as "Blank", add 20 µl of **1x Kinase assay buffer**.

6. Thaw **HIPK3 kinase** on ice. Briefly spin the tube to recover its full contents. Dilute the protein kinase to 4 ng/ $\mu$ l using **1x Kinase assay buffer**.

Notes: the concentration of protein is lot-specific and is indicated on the tube. Verify the initial concentration and dilute accordingly.

This kinase is particularly sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use the thawed protein and do not re-use the diluted kinase.

7. Add 20  $\mu$ l of diluted HIPK3 to the wells designated "Positive Control" and "Test Inhibitor."
8. Preincubate the test inhibitor and HIPK3 for 30 minutes at room temperature.
9. Prepare the Master Mix (25  $\mu$ l/well): N wells x (5  $\mu$ l of **5x Kinase assay buffer** + 1  $\mu$ l of **ATP (500  $\mu$ M)** + 1  $\mu$ l of **Myelin basic protein (MBP), 5 mg/ml** + 18  $\mu$ l of distilled water). Add 25  $\mu$ l to every well.
10. Initiate the reaction by adding 25  $\mu$ l to the wells designated "Blank," "Positive Control," and "Test Inhibitor."

Component	Blank	Positive Control	Test Inhibitor
Master Mix	25 $\mu$ l	25 $\mu$ l	25 $\mu$ l
Test Inhibitor	-	-	5 $\mu$ l
Diluent Solution	5 $\mu$ l	5 $\mu$ l	-
1x Kinase Buffer	20 $\mu$ l	-	-
HIPK3 (4 ng/ $\mu$ l)	-	20 $\mu$ l	20 $\mu$ l
<b>Total</b>	<b>50 <math>\mu</math>l</b>	<b>50 <math>\mu</math>l</b>	<b>50 <math>\mu</math>l</b>

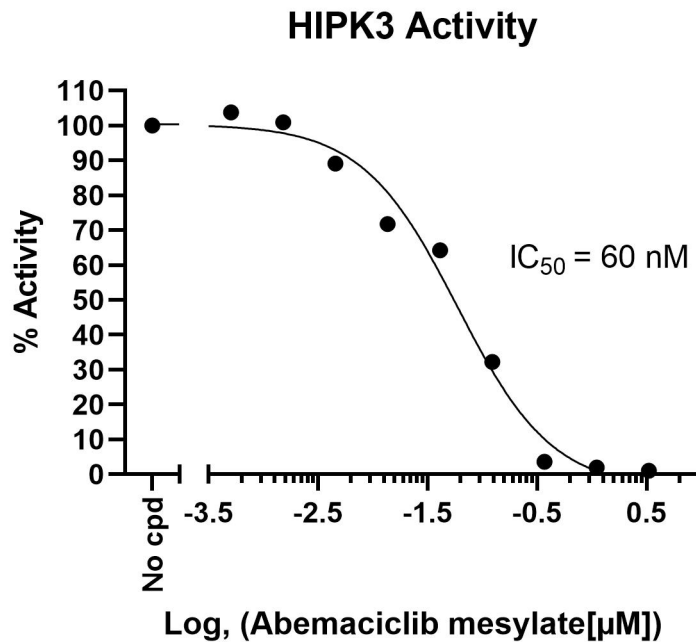
11. Incubate at 30°C for 45 minutes.
12. During the incubation, thaw the Kinase-Glo Max reagent. At the end of the 45-minute reaction, add 50  $\mu$ l of Kinase-Glo Max reagent to each well. Cover the plate with aluminum foil and incubate the plate at room temperature for 15 minutes.
13. Immediately read in a luminometer or a microplate reader capable of reading luminescence. The "Blank" value is subtracted from all other readings.

### Reading Luminescence

Luminescence is the emission of light resulting from a chemical reaction. The detection of luminescence requires no wavelength selection because the method used is emission photometry and not emission spectrophotometry.

To properly read luminescence, make sure the plate reader is set for LUMINESCENCE mode. Typical integration time is 1 second, delay after plate movement is 100 msec. Do not use a filter when measuring light emission. Typical settings for the Synergy 2 BioTek plate reader: use the "hole" position on the filter wheel; Optics position: Top; Read type: endpoint. Sensitivity may be adjusted based on the luminescence of a control assay without enzyme (typically we set this value as 100).

## Example Results



*Figure 1: Inhibition of HIPK3 kinase activity by Abemaciclib mesylate.* The inhibition of HIPK3 kinase activity was measured in the presence of increasing inhibitor concentrations using the HIPK3 Kinase Assay Kit (BPS Bioscience #78382). Results are expressed as percent of control (kinase activity in the absence of inhibitor, set at 100%).

For lot-specific information, please contact BPS Bioscience, Inc. at [support@bpsbioscience.com](mailto:support@bpsbioscience.com).

### Troubleshooting Guide

Visit [bpsbioscience.com/assay-kits-faq](https://bpsbioscience.com/assay-kits-faq) for detailed troubleshooting instructions. For all further questions, please email [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

### References

- Li, Yawei, *et al.* "Circ HIPK 3 sponges miR-558 to suppress heparanase expression in bladder cancer cells." *EMBO reports* 2017; **18.9**: 1646-1659.
- Kai, Ding, *et al.* "Circular RNA HIPK3 promotes gallbladder cancer cell growth by sponging microRNA-124." *Biochemical and biophysical research communications* 2018; **503.2**: 863-869.

### Related Products

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
HIPK3, His-Tag	40110	10 µg
HIPK1, GST-Tag	40179	10 µg
HIPK4, GST-Tag	40111	10 µg