

# TRKC Kinase Assay Kit

## Description

The Tropomyosin receptor kinase C (TRKC) Kinase Assay Kit is designed to measure TRKC 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 TRKC kinase, kinase substrate, ATP, and kinase assay buffer for 100 enzyme reactions.

## Background

Tropomyosin receptor kinase C (TRKC), also known as neurotrophic tyrosine kinase receptor type 3, is a member of the receptor tyrosine kinase family that mediates neurotrophin-activated signaling. TRKC is a functional coordinator in the activation of a host of signaling cascades including the mitogen-activated protein kinase (MAPK) pathway that leads to cell differentiation. Most recently, it has been shown that TRKC levels in hippocampal neurons dictate TRKC-mediated synapse density. TRKC expression levels are known to play a role in tumorigenesis as well as neurodegenerative diseases. Studies involving TRKC may lead to further understanding of disease pathology and to therapeutic treatments.

## Application(s)

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

## Supplied Materials

Catalog #	Name	Amount	Storage
40282	TRKC, GST-tag Recombinant*	10 µg	-80°C
79334	Kinase assay buffer 1 (5x)	1.5 ml	-20°C
79686	ATP (500 µM)	100 µl	-20°C
40217	PTK substrate Poly (Glu:Tyr 4:1) (10 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
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 100 μM 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 **PTK substrate Poly (Glu:Tyr 4:1) substrate**.
2. Prepare 3 ml of **1x Kinase assay buffer 1** by mixing 600 μl of **5x Kinase assay buffer 1** with 2400 μl water. Three (3) ml of **1x Kinase assay buffer 1** is sufficient for 100 reactions.
3. Prepare the **Master Mix** (25 μl/well): N wells x (5 μl of **5x Kinase assay buffer 1** + 1 μl of **ATP (500 μM)** + 1 μl of PTK substrate Poly (Glu:Tyr 4:1) (10 mg/ml) + 18 μl of distilled water). Add 25 μl to every well.
4. 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.

If the Test Inhibitor is water-soluble (*without DMSO*):

- 4.1 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.
- 4.2 For the positive and negative controls, use 1x Kinase Assay Buffer 1 (Diluent Solution).

**Or**

If the Test Inhibitor is soluble in DMSO (*with DMSO*):

- 4.1 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%.
- 4.2 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.
- 4.3 For positive and negative controls, prepare 10% DMSO in water (vol/vol) so that all wells contain the same amount of DMSO (Diluent Solution).

*Note: The final concentration of DMSO should not exceed 1%.*

5. 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).
6. To the wells designated as "Blank," add 20 μl of **1x Kinase assay buffer 1**.

7. Thaw **TRKC kinase** on ice. Briefly spin the tube to recover its full contents. Dilute the protein kinase to 5 ng/μl using **1x Kinase assay buffer 1**.

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



*Note: 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.*

8. Initiate the reaction by adding 20 μl of diluted Kinase to the wells designated “Positive Control” and “Test Inhibitor”.

Fill in table as appropriate

<b>Component</b>	<b>Blank</b>	<b>Positive Control</b>	<b>Test Inhibitor</b>
Master Mix	25 μl	25 μl	25 μl
Test Inhibitor	-	-	5 μl
Diluent Solution	5 μl	5 μl	-
1x Kinase Buffer	20 μl	-	-
TRKC (5 ng/μl)	-	20 μl	20 μl
<b>Total</b>	<b>50 μl</b>	<b>50 μl</b>	<b>50 μl</b>

9. Incubate at 30°C for 45 minutes.
10. During the incubation, thaw the Kinase-Glo Max reagent. At the end of the 45-minute reaction, add 50 μl of Kinase-Glo Max reagent to each well. Cover the plate with aluminum foil and incubate the plate at room temperature for 15 minutes.
11. 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

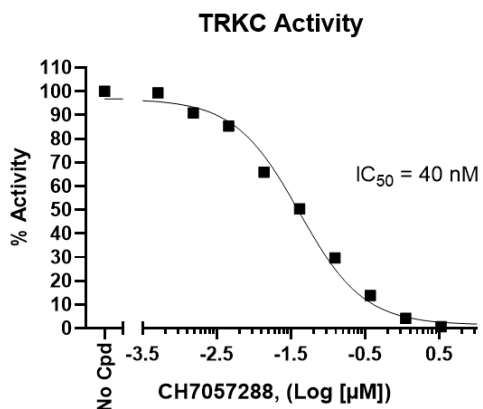


Figure 1: Inhibition of TRKC kinase Activity by CH7057288 (Selleckchem #S8788).

The inhibition of TRKC kinase activity was measured in the presence of increasing inhibitor concentrations using the TRKC Kinase Assay Kit (BPS Bioscience #78528). The Blank value was subtracted from all other values. Results are expressed as percent of control (kinase activity in the absence of inhibitor, set at 100%).

Data shown is representative, 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

1. Cocco, E., *et al.* Nature medicine, 25(9), 1422–1427.
2. Han, K. A., *et al.* Journal of Neuroscience 27 April 2016, 36 (17) 4816-4831;
3. Jin W. Roles of TrkC Signaling in the Regulation of Tumorigenicity and Metastasis of Cancer. *Cancers (Basel)*. 2020;12(1):147. Published 2020 Jan 8. doi:10.3390/cancers12010147

## Related Products

Products	Catalog #	Size
TRKC (G623R) Mutant, Active) Recombinant	40203	10 µg
TRKC (G623E) Mutant, Active) Recombinant	40204	10 µg
TRKC (G623R L686M) Mutant, Active Recombinant	40215	10 µg
TRKC (L686M), Mutant, Active Recombinant	40216	10 µg