



6405 Mira Mesa Blvd, Ste 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

## Data Sheet

### ***RIPK1 Kinase Assay Kit (96-well)***

**Catalog #79560**

**Background:** RIPK1 is a member of the receptor-interacting protein kinase (RIP) family. It plays an important role in cell survival as well as cell death, including apoptosis and necroptosis. RIPK1 is highly expressed in Alzheimer's disease (AD) brains. Importantly, it was shown that inhibition of RIPK1 results in decreased amyloid and inflammation levels as well as accelerated amyloid beta peptide (A $\beta$ ) degradation by microglia *in vitro*.

**Description:** The *RIPK1 Kinase Assay Kit* is designed to measure RIPK1 kinase activity for screening and profiling applications using ADP-Glo<sup>®</sup> Kinase Assay as a detection reagent. The *RIPK1 Kinase Assay Kit* comes in a convenient 96-well format, with enough purified recombinant RIPK1 enzyme, RIPK1 substrate, ATP and kinase assay buffer for 100 enzyme reactions.

MBP (Myelin Basic Protein) is a non-specific protein substrate that is used as a "universal substrate" for many in-vitro kinase activity assays. This protein is targeted by many serine/threonine kinases at conserved amino acids. We use the dephosphorylated version of the MBP substrate in our assays to determine the kinase-mediated phosphorylation of MBP. Our assays are not suitable for studying autophosphorylation of the kinase due to the presence of the MBP substrate.

#### **COMPONENTS:**

| Catalog # | Reagent                | Amount      | Storage    |   |
|-----------|------------------------|-------------|------------|---|
| 40371     | RIPK1                  | 10 $\mu$ g  | -80°C      | <b>Avoid multiple freeze/thaw cycles!</b> |
| 79334     | 5x Kinase assay buffer | 1.5 ml      | -20°C      |   |
| 79686     | ATP (500 $\mu$ M)      | 100 $\mu$ l | -20°C      |   |
| 78514     | MBP (5 mg/ml)          | 100 $\mu$ l | -20°C      |   |
| 79696     | 96-well plate, white   | 1           | Room Temp. |   |

#### **MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:**

ADP-Glo<sup>®</sup> Kinase Assay (Promega #V6930)  
Dithiothreitol (DTT, 1 M; optional)  
Microplate reader capable of reading luminescence  
Adjustable micropipettor and sterile tips  
30°C incubator

**APPLICATIONS:** Useful for studying enzyme kinetics and screening small molecular inhibitors for drug discovery and HTS applications.

**STABILITY:** Up to 6 months when stored as recommended.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd, Ste 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: support@bpsbioscience.com

**CONTRAINDICATION:** Avoid >1% DMSO

**REFERENCE:**

Ofengeim D., *et al. Proc. Natl. Acad. Sci. USA* **114(41)**: E8788-E8797 (2017)

**ASSAY PROTOCOL:**

*All samples and controls should be tested in duplicate.*

- 1) Thaw **5x Kinase assay buffer**, **ATP** and **RIPK1 substrate (MBP)**. (Optional: If desired, add DTT to **5x Kinase assay buffer** to make a 10 mM concentration; e.g. add 10  $\mu$ l of 1 M DTT to 1 ml **5x Kinase assay buffer**. Prepare only enough 5x Kinase assay buffer with DTT as required for the assay, as any excess 5x kinase buffer/DTT cannot be stored and should be discarded.)
- 2) Prepare the master mixture (12.5  $\mu$ l per well): N wells x (3  $\mu$ l **5x Kinase assay buffer** + 0.5  $\mu$ l **ATP (500  $\mu$ M)** + 1  $\mu$ l **RIPK1 substrate (MBP)** + 8  $\mu$ l distilled water). Add 12.5  $\mu$ l to every well.

|                                      | Positive Control            | Test Inhibitor              | Blank                       |
|--------------------------------------|-----------------------------|-----------------------------|-----------------------------|
| 5x Kinase assay buffer               | 3 $\mu$ l                   | 3 $\mu$ l                   | 3 $\mu$ l                   |
| ATP (500 $\mu$ M)                    | 0.5 $\mu$ l                 | 0.5 $\mu$ l                 | 0.5 $\mu$ l                 |
| RIPK1 substrate (MBP)                | 1 $\mu$ l                   | 1 $\mu$ l                   | 1 $\mu$ l                   |
| Water                                | 8 $\mu$ l                   | 8 $\mu$ l                   | 8 $\mu$ l                   |
| Test Inhibitor                       | -                           | 2.5 $\mu$ l                 | -                           |
| Inhibitor Buffer (e.g. 10% DMSO(aq)) | 2.5 $\mu$ l                 | -                           | 2.5 $\mu$ l                 |
| 1x Kinase buffer                     | -                           | -                           | 10 $\mu$ l                  |
| RIPK1 (~10 ng/ $\mu$ l)              | 10 $\mu$ l                  | 10 $\mu$ l                  | -                           |
| <b>Total</b>                         | <b>25 <math>\mu</math>l</b> | <b>25 <math>\mu</math>l</b> | <b>25 <math>\mu</math>l</b> |

- 3) Prepare 10X concentrated inhibitor in an aqueous-based solution. *Note: Final DMSO concentration must be  $\leq$ 1%. Higher DMSO levels can significantly decrease the enzyme activity. For example, to test an inhibitor dissolved in 100% DMSO at 10  $\mu$ M, dilute 1 mM inhibitor with water to make a 100  $\mu$ M inhibitor in 10% DMSO(aq). Then, add 2.5  $\mu$ l of the 100  $\mu$ M solution to the assay to make a 1% DMSO concentration in the final reaction mixture.*
- 4) Add 2.5  $\mu$ l of Inhibitor solution of each well labeled as "Test Inhibitor". For the "Positive Control" and "Blank", add 2.5  $\mu$ l of the same solution without inhibitor (Inhibitor buffer, e.g. 10% DMSO(aq)).

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd, Ste 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

- 5) Prepare 3 ml of **1x Kinase assay buffer** by mixing 600  $\mu$ l of **5x Kinase assay buffer** with 2400  $\mu$ l water. 3 ml of **1x Kinase assay buffer** is sufficient for 100 reactions.
- 6) To the wells designated as "Blank", add 10  $\mu$ l of **1x Kinase assay buffer**.
- 7) Thaw **RIPK1 enzyme** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Calculate the amount of **RIPK1** required for the assay and dilute enzyme to 10 ng/ $\mu$ l with **1x Kinase assay buffer**. Store remaining undiluted enzyme in aliquots at -80°C. *Note: RIPK1 enzyme is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 8) Initiate reaction by adding 10  $\mu$ l of **diluted RIPK1 enzyme** to the wells designated "Positive Control" and "Test Inhibitor Control". Carefully shake the plate well and incubate it at 30°C for 50 minutes.
- 9) Thaw ADP-Glo reagent.
- 10) After the 50 minutes reaction, add 25  $\mu$ l of ADP-Glo reagent to each well. Cover plate with aluminum foil and incubate the plate at room temperature for 45 minutes.
- 11) Thaw Kinase Detection reagent.
- 12) After the 45 minutes incubation, add 50  $\mu$ l of Kinase Detection reagent to each well. Cover plate with aluminum foil and incubate the plate at room temperature for another 45 minutes.
- 13) Measure luminescence using the microplate reader. "Blank" value should be subtracted from all wells.

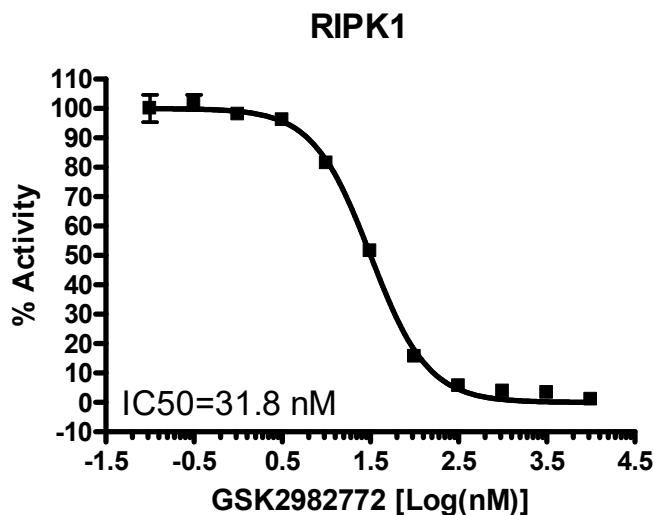
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

**Example of Assay Results:**



Inhibition of RIPK1 enzyme by GSK2982772, measured using the *RIPK1 kinase assay kit* (Cat. #79560). *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com*

**RELATED PRODUCTS:**

| <u>Product Name</u>     | <u>Catalog #</u> | <u>Size</u> |
|-------------------------|------------------|-------------|
| RIPK1, GST-Th-Tag       | 40371            | 10 µg       |
| RIPK2, GST-tag          | 40173            | 10 µg       |
| RIPK5, GST-tag          | 40174            | 10 µg       |
| LRRK2 (RIPK7), GST-Tag  | 40842            | 10 µg       |
| LRRK2 (Y1699C), GST-Tag | 40838            | 10 µg       |
| LRRK2 (Y1699G), GST-Tag | 40839            | 10 µg       |
| LRRK2 (R1441H), GST-Tag | 40837            | 10 µg       |
| LRRK2 (R1441G), GST-Tag | 40836            | 10 µg       |
| LRRK2 (R1441C), GST-Tag | 40835            | 10 µg       |
| LRRK2 (G2385R), GST-Tag | 40833            | 10 µg       |
| LRRK2 (G2019S), GST-Tag | 40832            | 10 µg       |
| LRRK2 (D1994A), GST-Tag | 40831            | 10 µg       |
| LRRK2 (I2020T), GST-Tag | 40834            | 10 µg       |

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)