



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Data Sheet**

### ***TNKS2 Histone Ribosylation Assay Kit (Antibody Detection)*** **Catalog # 80576**

**DESCRIPTION:** The *TNKS2 Histone Ribosylation Assay Kit (Antibody Detection)* is designed to measure Tankyrase-2 activity for screening and profiling applications. The *TNKS2 Histone Ribosylation Assay Kit (Antibody Detection)* comes in a convenient format, with 8-well strips, histone mixture substrate, antibody against poly (ADP-ribose) modified histone, the secondary HRP-labeled antibody, NAD<sup>+</sup>, PARP assay buffer, and purified TNKS2 enzyme for 100 enzyme reactions. The key to the *TNKS2 Histone Ribosylation Assay Kit (Antibody Detection)* is a highly specific antibody that recognizes poly (ADP-ribose)-modified histone. After coating the plate with the histone mixture, only three simple steps are required. First, NAD<sup>+</sup> is incubated with a sample containing assay buffer and enzyme for one hour. Next, primary antibody is added. Finally, the plate is treated with an HRP-labeled secondary antibody followed by addition of the HRP substrate to produce chemiluminescence that can be measured using a chemiluminescence reader.

*Note: This kit is also suitable for use with cell extracts, but non-specific ribosylation activity may be detected.*

#### **COMPONENTS:**

| Catalog # | Reagent                            | Amount                          | Storage |                                    |
|-----------|------------------------------------|---------------------------------|---------|------------------------------------|
| 80515     | TNKS2                              | 2 µg                            | -80°C   | <b>(Avoid freeze/thaw cycles!)</b> |
| 52029     | 5x histone mixture                 | 1 ml                            | -80°C   |                                    |
|           | NAD <sup>+</sup>                   | 250 µl                          | -80°C   |                                    |
| 80602     | 10x PARP Assay Buffer              | 1 ml                            | -20°C   |                                    |
| 52140K    | Primary antibody 11                | 12.5 µl                         | -80°C   |                                    |
| 52130H    | Secondary HRP-labeled antibody 1   | 10 µl                           | -80°C   |                                    |
| 79743     | Blocking buffer 3                  | 2 x 40 ml                       | +4°C    |                                    |
| 79670     | ELISA ECL substrate (2 components) | 6 ml each                       | +4°C    |                                    |
|           | 8-well strip plate module          | 1 plate<br>(12 x 8-well strips) | +4°C    |                                    |

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

PBS buffer  
PBST buffer (1x PBS, containing 0.05% Tween-20)  
Luminometer or fluorescent microplate reader capable of reading chemiluminescence  
Adjustable micropipettor and sterile tips  
Rotating or rocker platform  
Paper towels

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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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

**CONTRAINDICATIONS:** DMSO >1%, strong acids or bases, ionic detergents, high salt

**STABILITY:** 6 months from date of receipt when stored as directed.

**REFERENCE:** Dillon SC, Zhang X, Trievel RC, Cheng X. *Genome Biology* 2005; **6**:227.

**ASSAY PROTOCOL:**

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

Coating the plate with the histone mixture:

- 1) Dilute 5x histone mixture 1:5 in PBS.
- 2) Add 50 µl diluted histone mixture to each well and incubate for overnight at 4°C.
- 3) Wash the plate 3 times with 200 µl PBST buffer.
- 4) Block the wells by adding 150 µl of Blocking buffer 3 to every well. Incubate for 30 min at room temperature.
- 5) Wash the plate 3 times with 200 µl PBST buffer

**Step 1:**

- 1) Prepare the master mixture: N wells x (5 µl **10x PARP assay buffer** + 2.5 µl **NAD<sup>+</sup>** + 17.5 µl **H<sub>2</sub>O**)
- 2) Thaw **TNKS2 enzyme** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **TNKS2 enzyme** into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C immediately. *Note: TNKS2 enzyme is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 3) Dilute **TNKS2 enzyme** in 1X PARP assay buffer at 1 ng/µl (20 ng/20 µl). Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.
- 4) Add 25 µl of master mixture to each well designated for the "Positive Control", "Test Inhibitor", and "Blank". For the "Substrate Control", add 5 µl **10x PARP assay buffer** + 20 µl **H<sub>2</sub>O**.

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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

|                                 | Positive Control | Test Inhibitor | Substrate Control | Blank   |
|---------------------------------|------------------|----------------|-------------------|---------|
| 10X PARP Assay Buffer           | 5 µl             | 5 µl           | 5 µl              | 5 µl    |
| NAD <sup>+</sup>                | 2.5 µl           | 2.5 µl         | –                 | 2.5 µl  |
| H <sub>2</sub> O                | 17.5 µl          | 17.5 µl        | 20 µl             | 17.5 µl |
| Test Inhibitor                  | –                | 5 µl           | –                 | –       |
| Inhibitor buffer (no inhibitor) | 5 µl             |                | 5 µl              | 5 µl    |
| 1 x PARP buffer                 | -                | -              | -                 | 20 µl   |
| TNKS2 (1 ng/µl)                 | 20 µl            | 20 µl          | 20 µl             | –       |
| Total                           | 50 µl            | 50 µl          | 50 µl             | 50 µl   |

- 5) Add 5 µl of inhibitor solution to each well designated "Test Inhibitor". For the "Positive Control", "Substrate Control", and "Blank", add 5 µl of the same solution without inhibitor (inhibitor buffer).
- 6) Add 20 µl of 1 x PARP buffer to the well designated "Blank".
- 7) Initiate TNKS reactions by adding 20 µl of diluted TNKS2 prepared above. Incubate the reactions for 1 hour at room temperature.
- 8) Wash the plate 3 times with 200 µl PBST.
- 9) Add 100 µl of Blocking buffer 3 to every well. Shake on a rotating platform for 10 min. Remove supernatant as above.

## Step 2:

- 1) Dilute **Primary antibody 11** 800-fold with Blocking buffer 3.
- 2) Add 100 µl per well. Incubate 1 hour at room temperature with slow shaking.
- 3) Wash strip plate with PBST buffer and Blocking buffer 3 as in steps 1-8 and 1-9.

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

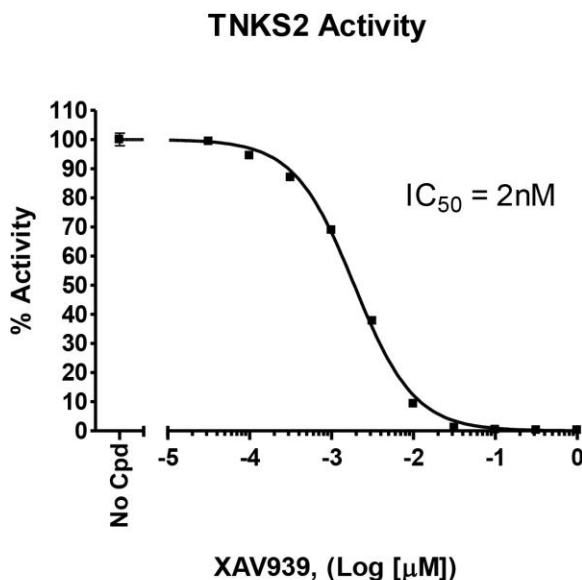


6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

### Step 3:

- 1) Dilute **Secondary HRP-labeled antibody 1** 1,000-fold with Blocking buffer 3.
- 2) Add 100  $\mu$ l per well. Incubate for 30 min. at room temperature with slow shaking.
- 3) Wash strip plate with PBST buffer and Blocking buffer 3 as in steps 1-8 and 1-9.
- 4) Just before use, mix on ice 50  $\mu$ l **ELISA ECL substrate A** and 50  $\mu$ l **ELISA ECL substrate B** and add 100  $\mu$ l per well. Discard any unused chemiluminescent reagent after use.
- 5) Immediately read sample in a luminometer or microtiter-plate capable of reading chemiluminescence.

### Example of Assay Results:



Inhibition of TNKS2 by XAV939, measured using the *TNKS2 Histone Ribosylation Assay Kit (Antibody Detection)*, BPS Bioscience #80576. Luminescence was measured using a Bio-Tek fluorescent microplate reader.

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

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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## RELATED PRODUCTS

| <u>Product Name</u>  | <u>Catalog #</u> | <u>Size</u> |
|--|------------------|-------------|
| TNKS1 Histone Ribosylation Assay Kit<br>(Biotin-labeled NAD <sup>+</sup> ) | 80574            | 96 rxns.    |
| PARP1 Chemiluminescent Assay Kit   | 80551            | 96 rxns.    |
| PARP2 Chemiluminescent Assay Kit   | 80552            | 96 rxns.    |
| PARP3 Chemiluminescent Assay Kit   | 80553            | 96 rxns.    |
| PARP6 Chemiluminescent Assay Kit   | 80556            | 32 rxns.    |
| PARP7 Chemiluminescent Assay Kit   | 80557            | 96 rxns.    |
| PARP11 Chemiluminescent Assay Kit  | 80561            | 96 rxns.    |
| PARP1 Enzyme   | 80501            | 10 µg       |
| PARP2 Enzyme   | 80502            | 10 µg       |
| PARP3 Enzyme   | 80503            | 10 µg       |
| PARP6 Enzyme   | 80506            | 10 µg       |
| TNKS1 (PARP5A) Enzyme  | 80504            | 10 µg       |
| TNKS2 (PARP5B), (667-end) Enzyme   | 80505            | 10 µg       |
| TNKS2 (PARP5B), (849-end) Enzyme   | 80515            | 10 µg       |
| PARP7 Enzyme   | 80507            | 10 µg       |
| PARP9 Enzyme   | 80509            | 10 µg       |
| PARP11 Enzyme  | 80511            | 10 µg       |
| PARP12 Enzyme  | 80512            | 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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



6042 Cornerstone Court W, Ste B  
 San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## TROUBLESHOOTING GUIDE

| Problem  | Possible Cause                                    | Solution   |
|--|---|--|
| Luminescence signal of positive control reaction is weak   | TNKS2 enzyme has lost activity                    | Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh enzyme (TNKS2, BPS Bioscience #80515). Store enzyme in single-use aliquots. Increase time of enzyme incubation. Increase enzyme concentration. |
|  | Antibody reaction is insufficient                 | Increase time for primary antibody incubation. Avoid freeze/thaw cycles of antibodies.   |
|  | Incorrect settings on instruments                 | Record light signals at 5 second intervals. Refer to instrument instructions for settings to increase sensitivity of light detection.  |
|  | Chemiluminescent reagents mixed too soon          | Chemiluminescent solution should be used within 15 minutes of mixing. Ensure both reagents are properly mixed.   |
| Luminescent signal is erratic or varies widely among wells | Inaccurate pipetting/technique                    | Run duplicates of all reactions. Use a multichannel pipettor. Use master mixes to minimize errors.   |
|  | Bubbles in wells                                  | Pipette slowly to avoid bubble formation. Tap plate lightly to disperse bubbles; be careful not to splash between wells.   |
| Background (signal to noise ratio) is high                 | Insufficient washes                               | Increase number of washes. Increase wash volume. Increase Tween-20 concentration to 0.1% in PBST.  |
|  | Sample solvent is inhibiting the enzyme           | Run negative control assay including solvent. Maintain DMSO level at <1%. Increase time of enzyme incubation.  |
|  | Results are outside the linear range of the assay | Use different concentrations of enzyme (TNKS2, BPS Bioscience #80515) to create a standard curve.  |

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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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