



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

## **Data Sheet**

### **GSK3 $\beta$ Assay Kit**

**Catalog #79700**

**Background:** GSK3 (Glycogen synthase kinase) is a serine/threonine kinase which initially was found to phosphorylate and inactivate glycogen synthase in glycogen biosynthesis. It is an important regulator enzyme in many disease pathogeneses including cancer, immune disorders, metabolic disorders and neurological disorders like Alzheimer's disease. In fact, GSK3 $\beta$ , one of the isoforms of GSK3, was identified as a key regulator in tau-driven Alzheimer's disease.

**Description:** The *GSK3 $\beta$  Assay Kit* is designed to measure *GSK3 $\beta$*  activity for screening and profiling applications using Kinase-Glo<sup>®</sup> (Promega) as a detection reagent. The *GSK3 $\beta$  Assay Kit* comes in a convenient 96-well format, with enough purified recombinant *GSK3 $\beta$*  enzyme, *GSK3 $\beta$*  substrate (GSK substrate peptide), ATP, and kinase assay buffer for 100 enzyme reactions.

#### **COMPONENTS:**

Catalog #	Reagent	Amount	Storage	
40007	GSK3 $\beta$	1.5 $\mu$ 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	
79697	10x GSK substrate peptide	500 $\mu$ l	-20°C	
	96-well plate, white	1	Room Temp.	

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

Kinase-Glo<sup>®</sup> Max Assay (Promega #V6071)  
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.

**CONTRAINDICATION:** Keep final DMSO concentration below 1%

#### **REFERENCE:**

McCubrey JA., *et al.* *Oncotarget* **5(10)**: 2881-2911 (2014)

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)

## ASSAY PROTOCOL:

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

- 1) Thaw **5x Kinase assay buffer**, **ATP** and **10x GSK substrate peptide**. (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 (25  $\mu$ l per well): N wells x (5  $\mu$ l **5x Kinase assay buffer** + 1  $\mu$ l **ATP (500  $\mu$ M)** + 5  $\mu$ l **10x GSK substrate peptide** + 14  $\mu$ l distilled water). Add 25  $\mu$ l to every well.

	Positive Control	Test Inhibitor	Blank
5x Kinase assay buffer	5 $\mu$ l	5 $\mu$ l	5 $\mu$ l
ATP (500 $\mu$ M)	1 $\mu$ l	1 $\mu$ l	1 $\mu$ l
10x GSK substrate peptide	5 $\mu$ l	5 $\mu$ l	5 $\mu$ l
Water	14 $\mu$ l	14 $\mu$ l	14 $\mu$ l
Test Inhibitor	-	5 $\mu$ l	-
Inhibitor Buffer (e.g. 10% DMSO(aq))	5 $\mu$ l	-	5 $\mu$ l
1x Kinase buffer	-	-	20 $\mu$ l
GSK3 $\beta$ (~0.6 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>

- 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 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 5  $\mu$ l of Inhibitor solution of each well labeled as "Test Inhibitor". For the "Positive Control" and "Blank", add 5  $\mu$ l of the same solution without inhibitor (Inhibitor buffer, e.g. 10% DMSO(aq)).
- 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. Dilute only enough **5x kinase assay buffer** as required for the assay; discard any remaining **1x kinase assay buffer**.

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

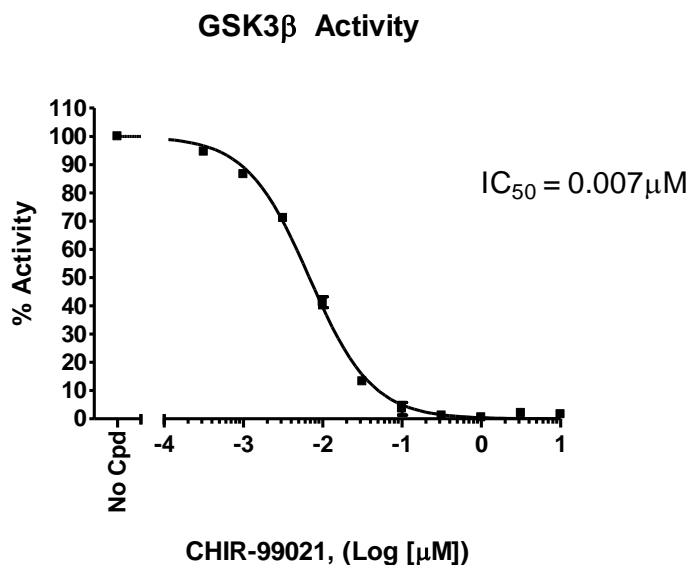
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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)

- 6) To the wells designated as "Blank", add 20  $\mu$ l of **1x Kinase assay buffer**.
- 7) Thaw **GSK3 $\beta$  enzyme** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Calculate the amount of **GSK3 $\beta$**  required for the assay and dilute enzyme to  $\sim$ 0.6 ng/ $\mu$ l with **1x Kinase assay buffer**. Store remaining undiluted enzyme in aliquots at  $-80^{\circ}\text{C}$ . *Note: GSK3 $\beta$  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 20  $\mu$ l of **diluted GSK3 $\beta$  enzyme** to the wells designated "Positive Control" and "Test Inhibitor Control". Carefully shake the plate well and incubate it at  $30^{\circ}\text{C}$  for 45 minutes.
- 9) Thaw Kinase-Glo Max reagent.
- 10) After the 42 minutes reaction, add 50  $\mu$ l of Kinase-Glo Max reagent to each well. Cover plate with aluminum foil and incubate the plate at room temperature for 15 minutes.
- 11) Measure luminescence using the microplate reader. "Blank" value should be subtracted from all wells.

### Example of Assay Results:



Inhibition of GSK3 $\beta$  enzyme by CHIR99021, measured using the *GSK3 $\beta$  kinase assay kit* (Cat. #79700). *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at 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:**

<b><u>Product Name</u></b>	<b><u>Catalog #</u></b>	<b><u>Size</u></b>
GSK3 $\alpha$	40006	10 $\mu$ g
GSK3 $\beta$	40007	10 $\mu$ g
TCF/LEF Reporter Kit (Wnt Signaling Pathway)	60500	500 rxns
CRE/CREB Reporter Kit (cAMP/PKA Signaling Pathway)	60611	500 rxns
CRE/CREB Luciferase Reporter Lentivirus	79580	500 $\mu$ l x 2

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)