



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**  
***GLS2 Inhibitor Screening Assay Kit***  
Catalog #79925  
Size: 96 reactions

**BACKGROUND:** Liver-type glutaminase (GLS2) is a phosphate-activated amidohydrolase that catalyzes the hydrolysis of L-glutamine to L-glutamate and ammonia. GLS2 is related to tumor progression and cancer, including breast and hepatocellular carcinoma.

**DESCRIPTION:** The *GLS2 Inhibitor Screening Assay Kit* is designed to measure the hydrolase activity of GLS2 for screening and profiling applications. The GLS2 assay kit comes in a convenient 96-well format, with purified GLS2, its substrates, the Coupling reagent, and buffer for 100 enzyme reactions.

**COMPONENTS:**

Catalog #	Component	Amount	Storage	
71242	GLS2, His-tag	30 µg	-80 °C	Avoid multiple freeze/thaw cycles!
	L-Glutamine (100 mM)	100 µl	-20 °C	
	NAD <sup>+</sup> (20 mM)	1 ml	-20 °C	
	Coupling reagent	10 µl	-20 °C	
	4X GLS assay buffer	2.5 ml	-20 °C	
79685	96-well black microplate	1	Room Temp	

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

0.5 M dithiothreitol (DTT) in aqueous solution  
Adjustable micropipettor and sterile tips  
Fluorescent microplate reader capable of reading  $\lambda_{excitation} = 340 \text{ nm}$ ;  $\lambda_{emission} = 460 \text{ nm}$

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

**STABILITY:** Up to 6 months from date of receipt, when stored as recommended.

**REFERENCES:**

1. Xiang, Lisha, *et al.* (2013) "Knock-down of glutaminase 2 expression decreases glutathione, NADH, and sensitizes cervical cancer to ionizing radiation." *Biochimica et Biophysica Acta (BBA)-Molecular Cell Research* **1833.12**: 2996-3005.
2. Zhang, Jianbin, *et al.* (2013) "Epigenetic silencing of glutaminase 2 in human liver and colon cancers." *BMC cancer* **13.1**: 601-609.

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)



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

#### ASSAY PROTOCOL:

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

- 1) Add 20  $\mu$ l of 0.5 M DTT (not supplied) to 2.5 ml 4X GLS2 assay buffer. Prepare **1X GLS buffer** by diluting **4X GLS buffer** with DTT 4-fold into water. For example, to prepare 10 ml, add 2.5 ml of **4X GLS buffer** and 20  $\mu$ l of 0.5 M DTT to 7.5 ml of water. (Final assay concentration of DTT is 1 mM).
- 2) Add 20  $\mu$ l **1X GLS buffer** to each well designated "Blank."
- 3) Thaw **GLS2** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Prepare **Enzyme solution (15 ng/ $\mu$ l GLS2)** by diluting **GLS2** in **1X GLS buffer**. Store remaining undiluted enzyme in aliquots at  $-80^{\circ}\text{C}$ . Note: **GLS2** is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.
- 4) Add 20  $\mu$ l **Enzyme solution (15 ng/ $\mu$ l GLS2)** to each well designated "Positive Control," and "Test Inhibitor."
- 5) Prepare 100x concentration of test compound in DMSO. Dilute test compound 1:10 in water, and add 5  $\mu$ l to each well designated "Test Inhibitor." For the wells labeled "Positive Control" and "Blank," add 5  $\mu$ l of 10% DMSO in water (Inhibitor buffer).
- 6) Prepare **Substrate solution** by diluting **L-Glutamine (100 mM)** 62.5-fold, **NAD<sup>+</sup> (20 mM)** 5-fold and **Coupling reagent** 600-fold in 1X GLS buffer. For example, to prepare 1000  $\mu$ l, add 16  $\mu$ l L-Glutamine (100 mM), 200  $\mu$ l NAD<sup>+</sup> (20 mM) and 1.66  $\mu$ l coupling reagent to 782  $\mu$ l 1X GLS buffer. Do not re-use Substrate solution.
- 7) Add 25  $\mu$ l **Substrate solution** to all wells. Read fluorescence intensity of the samples ( $\lambda_{\text{excitation}} = 340 \text{ nm}$ ;  $\lambda_{\text{emission}} = 460 \text{ nm}$ ) in an appropriate microplate reader ("t = 0 min reading"). Incubate reaction for 30 minutes at room temperature.

	Positive Control	Test Inhibitor	Blank
Enzyme solution (15 ng/ $\mu$ l GLS2)	20 $\mu$ l	20 $\mu$ l	-
1X GLS buffer	-	-	20 $\mu$ l
Test Inhibitor	-	5 $\mu$ l	-
10% DMSO in water (Inhibitor buffer)	5 $\mu$ l	-	5 $\mu$ l
Substrate solution	25 $\mu$ l	25 $\mu$ l	25 $\mu$ l
Total	50 $\mu$ l	50 $\mu$ l	50 $\mu$ l

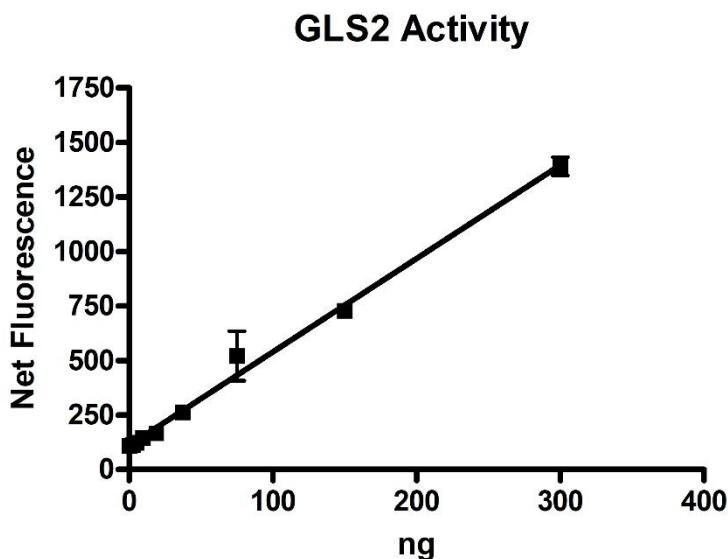
- 8) Read fluorescence intensity of the samples ( $\lambda_{\text{excitation}} = 340 \text{ nm}$ ;  $\lambda_{\text{emission}} = 460 \text{ nm}$ ) ("t = 30 min reading"). Subtract background fluorescence intensity values to get net fluorescence intensity: "t = 60 min reading" - "t = 0 min reading."

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)

Example of assay results:



GLS2 titration measured using the *GLS2 Inhibitor Screening Assay Kit*, BPS Bioscience, #79925. Fluorescence was measured using a Bio-Tek microplate reader. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**RELATED PRODUCTS:**

<u>Product</u>	<u>Catalog#</u>	<u>Size</u>
GLS1, His-tag	71102	20 µg
GLS2, His-tag	71242	20 µg
GLS1 Inhibitor Screening Assay Kit	79596	96 rxns.

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)