



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**

### ***AXL Kinase Assay Kit***

**Catalog #79711**

**Background:** AXL (TYRO7) is a receptor tyrosine kinase that is overexpressed in various forms of cancer, including breast, prostate, colorectal, and non-small cell cancers. AXL overexpression has been shown to encourage tumor progression and metastasis and to decrease the anti-tumor immune response, making AXL an attractive candidate for anticancer therapies.

**Description:** The *AXL Kinase Assay Kit* is designed to measure AXL kinase activity for screening and profiling applications using ADP-Glo<sup>®</sup> Kinase Assay as a detection reagent. The *AXL Kinase Assay Kit* comes in a convenient 96-well format, with enough purified recombinant AXL enzyme, AXL substrate (IRS1-tide), ATP and kinase assay buffer for 100 enzyme reactions.

#### **COMPONENTS:**

Catalog #	Reagent	Amount	Storage	
40180	AXL*	10 µg	-80°C	<b>Avoid multiple freeze/thaw cycles!</b>
79334	5x Kinase assay buffer	1.5 ml	-20°C	
79686	ATP (500 µM)	100 µl	-20°C	
79519	10x IRS1-tide (1 mg/ml)	500 µl	-20°C	
79696	96-well plate, white	1	Room Temp.	

\*The concentration of AXL is lot-specific and will be indicated on the tube containing the enzyme.

#### **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)

**REFERENCE:**

1. Wu, X., Liu, X., Koul, S., Lee, C.Y., Zhang, Z., Halmos, B. AXL kinase as a novel target for cancer therapy. *Oncotarget*. 2014; **5(20)**:9546-63.
2. Rankin, E.B., Giaccia, A.J. The Receptor Tyrosine Kinase AXL in Cancer Progression. *Cancers (Basel)*. 2016; **8(11)**:103.

**ASSAY PROTOCOL:**

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

- 1) Thaw **5x Kinase assay buffer**, **ATP (500  $\mu$ M)**, and **10x IRS1tide (1 mg/ml)**. (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**).
- 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)** + 5  $\mu$ l **10x IRS1tide (1 mg/ml)** + 4  $\mu$ l 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
IRS1tide 1 mg/ml	5 $\mu$ l	5 $\mu$ l	5 $\mu$ l
Water	4 $\mu$ l	4 $\mu$ l	4 $\mu$ l
Test Inhibitor	-	2.5 $\mu$ l	-
Inhibitor Buffer (no inhibitor)	2.5 $\mu$ l	-	2.5 $\mu$ l
1x Kinase buffer	-	-	10 $\mu$ l
AXL (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) 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). *Note: Keep DMSO concentration of the Test Inhibitor at  $\leq$ 10%, as final DMSO concentration in the reaction should be  $\leq$ 1%.*
- 4) 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.
- 5) To the wells designated as "Blank," add 10  $\mu$ l of **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: [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)

- 6) Thaw **AXL** enzyme on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Calculate the amount of **AXL** required for the assay and dilute enzyme to 10 ng/μl with **1x Kinase assay buffer**. Store remaining undiluted enzyme in aliquots at -80°C. *Note: AXL enzyme is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 7) Initiate reaction by adding 10 μl of diluted **AXL** enzyme to the wells designated "Positive Control" and "Test Inhibitor Control." Incubate at 30°C for 45 minutes.
- 8) Thaw ADP-Glo reagent.
- 9) After the 45 minutes reaction, add 25 μl of ADP-Glo reagent to each well. Cover plate with aluminum foil and incubate the plate at room temperature for 45 minutes.
- 10) Thaw Kinase Detection reagent.
- 11) After the 45 minutes incubation, add 50 μl of Kinase Detection reagent to each well. Cover plate with aluminum foil and incubate the plate at room temperature for another 45 minutes.
- 12) Immediately read sample in a luminometer or microtiter-plate capable of reading chemiluminescence. "Blank" value is subtracted from all readings.

### **Reading Chemiluminescence:**

Chemiluminescence is the emission of light (luminescence) which results from a chemical reaction. The detection of chemiluminescence requires no wavelength selection because the method used is emission photometry and is not emission spectrophotometry.

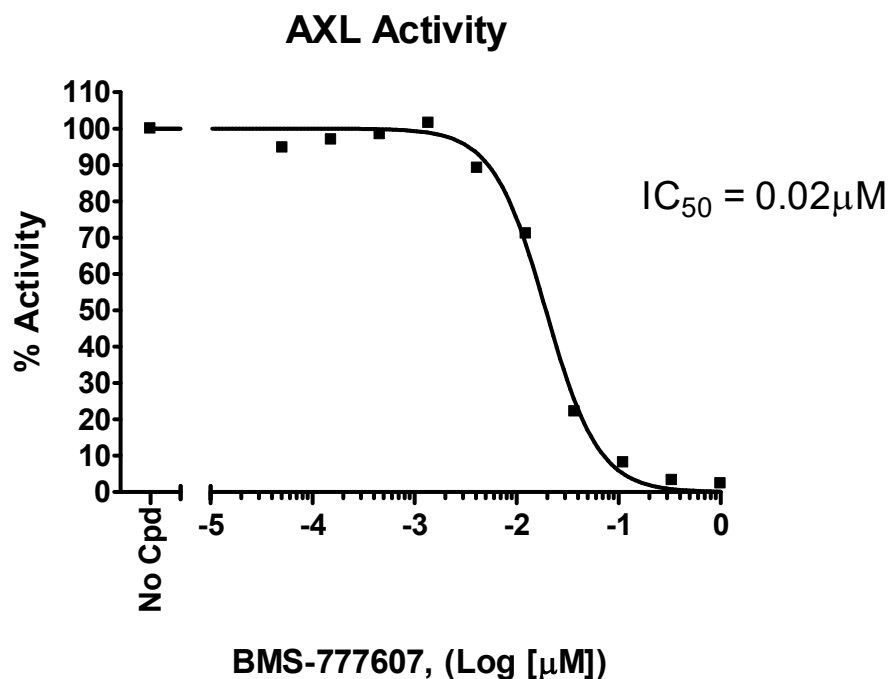
To properly read chemiluminescence, make sure the plate reader is set for LUMINESCENCE mode. Typical integration time is 1 second, delay after plate movement is 100 msec. Do not use a filter when measuring light emission. Typical settings for the Synergy 2 BioTek plate reader are: use the "hole" position 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).

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 AXL enzyme by BMS-777607, measured using the *AXL kinase assay kit* (Cat. #79711). *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at support@bpsbioscience.com*

**RELATED PRODUCTS:**

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
AXL (TYRO7), GST-tag	40180	10 µg
TYRO3, GST-tag	40293	10 µg
EPHA1, GST-tag	40191	10 µg
EPHA2, GST-tag	40190	10 µg
EPHA3 (TYRO4), GST-tag	40192	10 µg
EPHA4 (TYRO1), GST-tag	40193	10 µg
EPHB2 (TYRO5), His-tag	40200	10 µg
EPHB3 (TYRO6), His-tag	40186	10 µg
DDR2 (TYRO10), His-tag	40185	10 µg
5X Kinase assay buffer	79334	10 ml
TYRO3 Kinase Assay Kit	79593	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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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