



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
PIM1 Assay Kit
Catalog #79885
96 Reactions

Background: PIM kinases (PIM1, PIM2 and PIM3) are a family of serine/threonine protein kinases that play crucial roles in cell survival, proliferation, and drug resistance. PIM kinases are overexpressed in several tumors and promote growth and survival of malignant cells through cell cycle regulation and/or inhibition of apoptosis. Recently, PIM kinases were identified as a potential therapeutic target for precision medicine of advanced cancer.

DESCRIPTION: The *PIM1 Assay Kit* is designed to measure PIM1 activity for screening and profiling applications using Kinase-Glo[®] MAX as a detection reagent. The *PIM1 Assay Kit* comes in a convenient 96-well format, with enough purified recombinant PIM1 enzyme, PIM substrate (S6Ktide), ATP, and kinase assay buffer for 100 enzyme reactions.

COMPONENTS:

Catalog #	Reagent	Amount	Storage	
41107	PIM1	5 µg	-80°C	Avoid multiple freeze/ thaw cycles!
79334	5x Kinase assay buffer	1.5 ml	-20°C	
79686	ATP (500 µM)	100 µl	-20°C	
79884	PIM substrate (S6Ktide, 10 mg/ml)	100 µl	-20°C	
79696	96-well plate, white	1	Room Temp.	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Kinase-Glo MAX (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.

REFERENCE:

1. Jeyapal G. P., *et al.*, *Anticancer Agents Med Chem*, 2018, **18(8)**: 1100-1114
2. Asati V., *et al.*, *Eur J Med Chem*. 2019, **172**: 95-108

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

Please visit our website at: 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

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

- 1) Thaw **5x Kinase assay buffer**, **ATP (500 μ M)**, and **PIM Substrate (S6Ktide, 10 mg/ml)**.
(Optional: If desired, add DTT to **5x Kinase assay buffer** to make a 10 mM concentration; e.g. add 10 μ l of 1 M DTT to 1 ml **5x Kinase assay buffer**)
- 2) Prepare the master mixture (25 μ l per well): N wells x (6 μ l **5x Kinase assay buffer** + 1 μ l **ATP (500 μ M)** + 1 μ l **PIM Substrate (S6Ktide, 10 mg/ml)** + 17 μ l distilled water). Add 25 μ l to every well.

	Positive Control	Test Inhibitor	Blank
5x Kinase assay buffer	6 μ l	6 μ l	6 μ l
ATP (500 μ M)	1 μ l	1 μ l	1 μ l
PIM Substrate (S6Ktide, 10 mg/ml)	1 μ l	1 μ l	1 μ l
Distilled Water	17 μ l	17 μ l	17 μ l
Test Inhibitor	-	5 μ l	-
Inhibitor Buffer (e.g. 10% DMSO(aq))	5 μ l	-	5 μ l
1x Kinase buffer	-	-	20 μ l
PIM1 (2.5 ng/ μ l)	20 μ l	20 μ l	-
Total	50 μ l	50 μ l	50 μ l

- 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 μ M, dilute 1 mM inhibitor with water to make a 100 μ M inhibitor in 10% DMSO(aq). Then, add 5 μ l of the 100 μ M solution to the assay to make a 1% DMSO concentration in the final reaction mixture.*
- 4) Add 5 μ l of Inhibitor solution of each well labeled as "Test Inhibitor." For the "Positive Control" and "Blank," add 5 μ l of the same solution without inhibitor (Inhibitor buffer).
- 5) Prepare 3 ml of **1x Kinase assay buffer** by mixing 600 μ l of **5x Kinase assay buffer** with 2400 μ 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.
- 6) To the wells designated as "Blank," add 20 μ l of **1x Kinase assay buffer**.
- 7) Thaw **PIM1** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Calculate the amount of **PIM1** required for the assay and

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

Please visit our website at: www.bpsbioscience.com



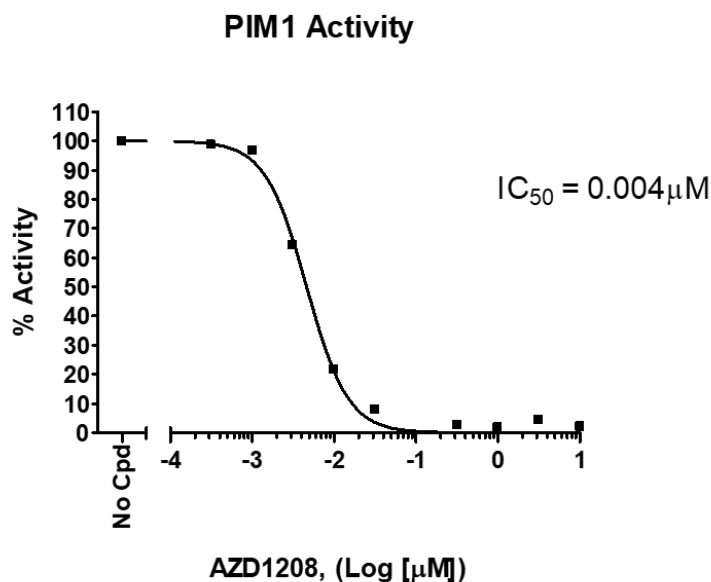
6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

dilute enzyme to ~2.5 ng/μl with **1x Kinase assay buffer**. Store remaining undiluted enzyme in aliquots at -80°C.

Note: PIM1 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 μl of diluted **PIM1** to the wells designated "Positive Control" and "Test Inhibitor." Incubate at 30°C for 45 minutes.
- 9) Thaw Kinase-Glo Max reagent.
- 10) After the 45 minutes reaction, add 50 μl 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 a microplate reader capable of reading chemiluminescence. "Blank" value is subtracted from all readings.

Example of Assay Results:



Inhibition of PIM1 by AZD1208 measured using the PIM1 assay kit (BPS Bioscience #79885). 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

Please visit our website at: 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

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
PIM1, Active, GST-tag	41107	10 µg
PIM2, Active, GST-tag	40153	10 µg
PIM3, Active, GST-tag	41108	10 µg
Kinase Buffer 1	79334	10 ml
ATP (500 µM)	79686	200 µl

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

Please visit our website at: www.bpsbioscience.com