



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 **JAK3 (Janus Kinase 3) Assay Kit** Catalog #79521

DESCRIPTION: Janus kinase (JAK) is a family of intracellular, nonreceptor tyrosine kinases that play an important role in cytokine-mediated signaling. This family includes four protein kinases; Jak1, Jak2, Jak3 and Tyk2. Various studies have suggested that regulation of Jak activity should be a promising approach to combat numerous diseases such as immune deficiencies and cancers. The *JAK3 (Janus Kinase 3) Assay Kit* is designed to measure JAK3 activity for screening and profiling applications using Kinase-Glo[®] MAX as a detection reagent. The *JAK3 (Janus Kinase 3) Assay Kit* comes in a convenient 96-well format, with enough purified recombinant JAK3 enzyme, JAK3 substrate peptide (Poly(Glu;Tyr, 4:1)), ATP and kinase assay buffer for 100 enzyme reactions.

COMPONENTS:

Catalog #	Reagent	Amount	Storage	
40452	JAK3	1 µg	-80°C	Avoid multiple freeze/thaw cycles!
79334	5x Kinase assay buffer 1	1.5 ml	-20°C	
79686	ATP (500 µM)	100 µl	-20°C	
40217	PTK substrate Poly (Glu:Tyr 4:1) (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:

Alicea, N.L. *et al.*, *Current Drug Targets* **12(4)**:546-555 (2011)

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

Please visit our website at: 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

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

- 1) Thaw **5x Kinase assay buffer**, **ATP** and **PTK Substrate Poly (Glu;Tyr, 4:1) (10 mg/ml)**.

(Please add DTT to 5x Kinase assay 1 buffer to make a 10 mM concentration; e.g. add 10 μ l of 1 M DTT to 1 ml 5x Kinase assay buffer 1. 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 μ l per well): N wells x (6 μ l **5x Kinase assay buffer 1** + 1 μ l **ATP (500 μ M)** + 1 μ l **PTK Substrate Poly (Glu;Tyr, 4:1) (10 mg/ml)** + 17 μ l distilled water). Add 25 μ l to every well.

	Positive Control	Test Inhibitor	Blank
5x Kinase assay buffer 1	6 μ l	6 μ l	6 μ l
ATP (500 μ M)	1 μ l	1 μ l	1 μ l
PTK Substrate (10 mg/ml)	1 μ l	1 μ l	1 μ l
Water	17 μ l	17 μ l	17 μ l
Test Inhibitor	-	5 μ l	-
Inhibitor Buffer (no inhibitor)	5 μ l	-	5 μ l
1x Kinase buffer 1	-	-	20 μ l
JAK3 (0.4 ng/ μ l)	20 μ l	20 μ l	-
Total	50 μ l	50 μ l	50 μ l

- 3) 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).
- 4) Prepare 3 ml of **1x Kinase assay buffer 1** by mixing 600 μ l of **5x Kinase assay buffer 1** with 2400 μ l water. 3 ml of **1x Kinase assay buffer 1** is sufficient for 100 reactions.
- 5) To the wells designated as "Blank", add 20 μ l of **1x Kinase assay buffer**.
- 6) Thaw **JAK3** enzyme on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Calculate the amount of **JAK3** required for the assay and dilute enzyme to ~0.4 ng/ μ l with **1x Kinase assay buffer**. Store remaining undiluted enzyme in aliquots at -80°C. *Note: JAK3 enzyme is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*

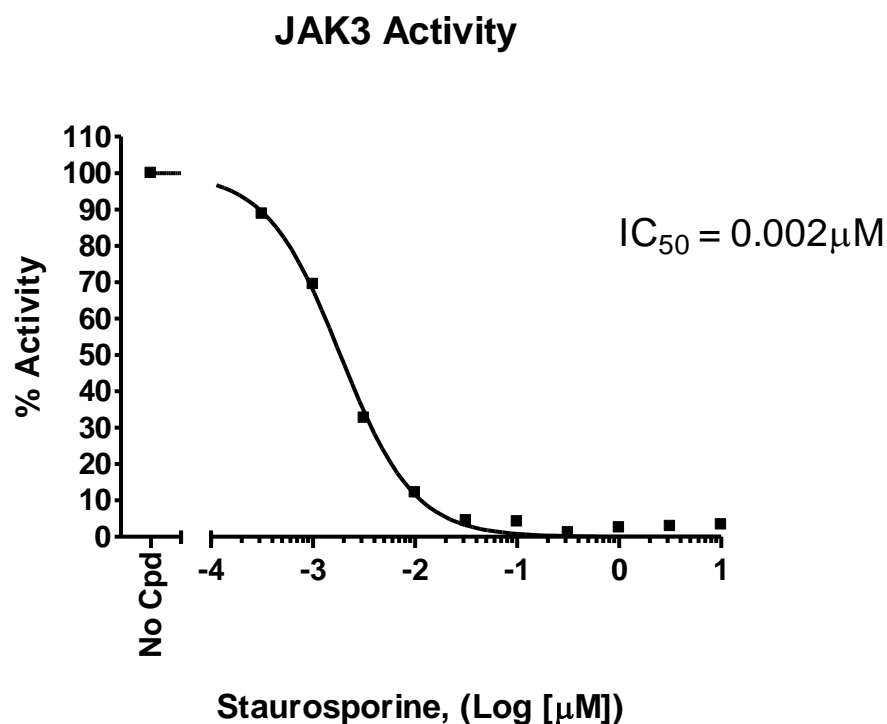
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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

Please visit our website at: www.bpsbioscience.com

- 7) Initiate reaction by adding 20 μ l of diluted **JAK3** enzyme to the wells designated "Positive Control" and "Test Inhibitor ". Incubate at 30°C for 45 minutes.
- 8) Thaw Kinase-Glo Max reagent.
- 9) After the 45 minute reaction, add 50 μ l of Kinase-Glo Max reagent to each well. Cover plate with aluminum foil and incubate the plate at room temperature for 15 minutes.
- 10) Measure luminescence using the microplate reader.

Example of Assay Results:

Inhibition of JAK3 enzyme by Staurosporine, measured using the *JAK3 (Janus Kinase 3) assay kit* (Cat. #79521). *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at support@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: support@bpsbioscience.com

Please visit our website at: 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

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
Jak1, GST-tag	40449	10 µg
Jak2 (JH1 domain), His-tag	40450	10 µg
Jak2 (JH1, JH2 domain), His-GST-tags	40451	10 µg
JAK2-His-Avi-Tag, Biotin-Labeled	79074	10 µg
Jak2 (V617F), GST/Avi-Tag, Biotin-Labeled	79115	10 µg
Jak3, His-tag	40452	10 µg
Tyk2, His-tag	40285	10 µg
ISRE Reporter Kit (JAK/STAT Pathway)	60613	500 rxns.
ISRE Reporter – HEK293 Cell Line	60510	2 vials
Transfection Collection™ : ISRE Transient Pack JAK/STAT Signaling Pathway	79264	500 rxns.
JAK1 (Janus Kinase 1) Assay Kit	79518	96 rxns.
JAK2 (Janus Kinase 2) Assay Kit	79520	96 rxns.
TYK2 (Tyrosine Kinase 2) Assay Kit	79521	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

Please visit our website at: www.bpsbioscience.com