



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

BTK Assay Kit

Catalog #79568

DESCRIPTION: Bruton's tyrosine kinase or BTK, is an enzyme that plays a role in the functionality and maturation of B cells. The BTK pathway has implications for a number of autoimmune disorders including isolated growth hormone deficiency type III and rheumatoid arthritis. The *BTK Assay Kit* is designed to measure BTK activity for screening and profiling applications using Kinase-Glo[®] MAX as a detection reagent. The *BTK Assay Kit* comes in a convenient 96-well format, with enough purified recombinant BTK enzyme, PolyGluTyr peptide, ATP, and kinase assay buffer for 100 enzyme reactions.

COMPONENTS:

Catalog #	Reagent	Amount	Storage	
40405	BTK	20 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	
40217	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:

Rawlings, D., Saffran, D., Tsukada, S., Largaespada, D., Grimaldi, J., Cohen, L., et al. (1993). Mutation of unique region of Bruton's tyrosine kinase in immunodeficient XID mice. *Science* **261(5119)**: 358–361.

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 (500 μ M)**, and **Poly(Glu,Tyr) (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 **PolyGluTyr (10 mg/ml)**.+ 17 μ l 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
Poly(Glu,Tyr) (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	-	-	20 μ l
BTK (10 ng/ μ l)	20 μ l	20 μ l	-
Total	50 μl	50 μl	50 μl

- 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). *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 μ l of **5x Kinase assay buffer** with 2400 μ l water. 3 ml of **1x Kinase assay buffer** is sufficient for 100 reactions.
- 5) To the wells designated as "Blank," add 20 μ l of **1x Kinase assay buffer**.
- 6) Thaw **BTK enzyme** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Calculate the amount of **BTK** required for the assay and dilute enzyme to 10 ng/ μ l with **1x Kinase assay buffer**. Store remaining undiluted enzyme in aliquots at -80°C. *Note: BTK 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



6405 Mira Mesa Blvd Ste. 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

- 7) Initiate reaction by adding 20 μ l of **diluted BTK enzyme** to the wells designated "Positive Control" and "Test Inhibitor Control." Incubate at 30°C for 60 minutes.
- 8) Thaw Kinase-Glo Max reagent.
- 9) After the 60 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) 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.

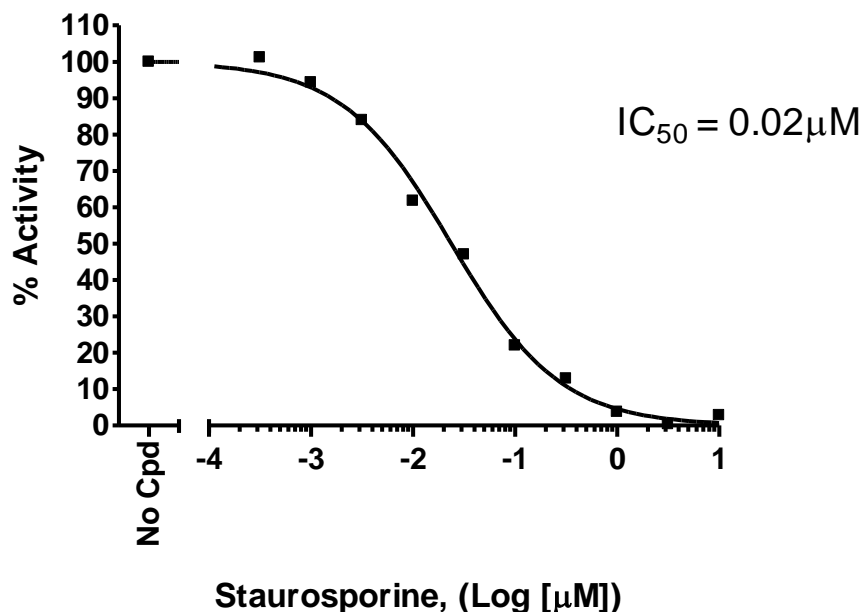
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

Example of Assay Results:

BTK Activity



Inhibition of BTK enzyme by Staurosporine (BPS Bioscience #27002), measured using the BTK assay kit (BPS Bioscience #79568). *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>
BTK, GST-tag	40405	10 µg
poly-(Glu,Tyr 4:1)	40217	1 mg
Staurosporine	27002	10 mg
5x Kinase assay buffer	79334	10 ml

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