



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

ALK2 Kinase Assay Kit

Catalog #79605

BACKGROUND:

ALK2, also known as activin receptor-like kinase 2 (ACVRLK2) is a receptor from the TGF- β family of transmembrane serine/threonine kinases that plays an important role in the bone morphogenic protein (BMP) pathway. Mutation of ALK2 is seen to play a role in cancer and fibrodysplasia ossificans progressiva (FOP).

DESCRIPTION:

The *ALK2 Kinase Assay Kit* is designed to measure ALK2 kinase activity for screening and profiling applications using ADP-Glo[®] Kinase Assay as a detection reagent. The *ALK2 Kinase Assay Kit* comes in a convenient 96-well format, with enough purified recombinant ALK2 enzyme, ALK2 substrate, ATP and kinase assay buffer for 100 enzyme reactions.

COMPONENTS:

Catalog #	Reagent	Amount	Storage	
40019	ALK2	10 μ g	-80°C	<i>Avoid multiple freeze/thaw cycles!</i>
79334	5x Kinase assay buffer	1.5 ml	-20°C	
	ATP (500 μ M)	100 μ l	-20°C	
79550	Casein (10 mg/ml)	100 μ l	-20°C	
	96-well plate, white	1	Room Temp.	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

ADP-Glo[®] 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 **1.858.829.3082** 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: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

REFERENCES:

1. Cunha, Sara I., and Kristian Pietras. "ALK2 as an Emerging Target for Antiangiogenic Therapy of Cancer." *Blood* **117(26)**: 6999–7006 (2011).
2. Minsung Kim, *et al.* Identification of novel ALK2 inhibitors and their effect on cancer cells, *Biochemical and Biophysical Research Communications*, **492(1)**:121-127 (2017).

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

- 1) Thaw **5x Kinase assay buffer**, **ATP (500 μ M)**, and **Casein (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 (12.5 μ l per well): N wells x (3 μ l **5x Kinase assay buffer** + 0.5 μ l **ATP (500 μ M)** + 0.5 μ l **Casein (10 mg/ml)** + 8.5 μ l distilled water). Add 12.5 μ l to every well.

	Positive Control	Test Inhibitor	Blank
5x Kinase assay buffer	3 μ l	3 μ l	3 μ l
ATP (500 μ M)	0.5 μ l	0.5 μ l	0.5 μ l
Casein 10 mg/ml	0.5 μ l	0.5 μ l	0.5 μ l
Water	8.5 μ l	8.5 μ l	8.5 μ l
Test Inhibitor	–	2.5 μ l	–
Inhibitor Buffer (no inhibitor)	2.5 μ l	–	2.5 μ l
1x Kinase buffer	–	–	10 μ l
ALK2 (5 ng/ μ l)	10 μ l	10 μ l	–
Total	25 μ l	25 μ l	25 μ l

- 3) Add 2.5 μ l of Inhibitor solution of each well labeled as "Test Inhibitor." For the "Positive Control" and "Blank," add 2.5 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 10 μ 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 **1.858.829.3082** 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: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 6) Thaw **ALK2** enzyme on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Calculate the amount of **ALK2** required for the assay and dilute enzyme to 5 ng/μl with **1x Kinase assay buffer**. Store remaining undiluted enzyme in aliquots at -80°C. *Note: ALK2 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 **ALK2** 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).

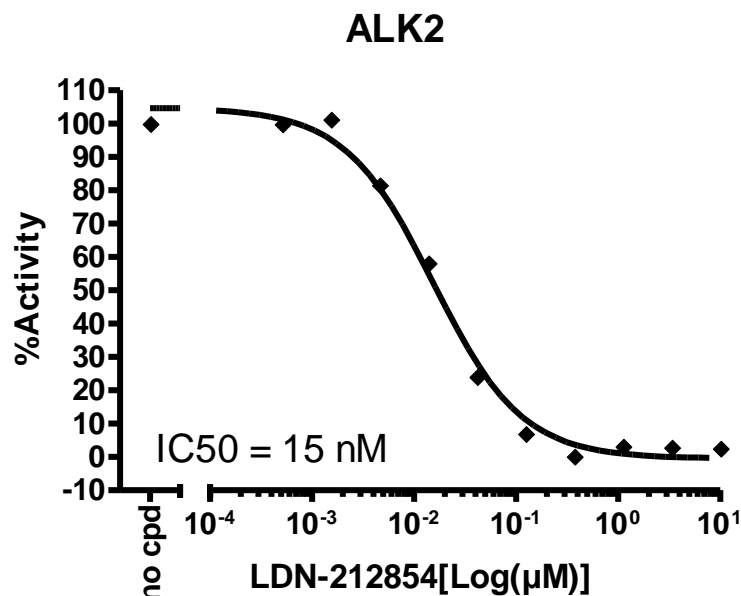
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com

Example of Assay Results:



Inhibition of ALK2 enzyme by LDN-212854, measured using the *ALK2 kinase assay kit* (Cat. #79605). *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com*

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
ALK1, GST-tag	40018	10 μg
ALK2, GST-tag	40019	10 μg
ALK4, GST-tag	40020	10 μg
ALK1 Kinase Assay Kit	79549	96 rxns.
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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com