



6405 Mira Mesa Blvd. Suite 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet ***IDO2 Inhibitor Screening Assay Kit*** Catalog # 72022

**DESCRIPTION:** The *IDO2 Inhibitor Screening Assay Kit* is designed to measure IDO2 enzyme inhibition. The kit comes in a convenient format with enough reaction solution and enzyme to perform a total of 100 reactions. The *IDO2 Inhibitor Screening Assay Kit* is simple to use. Inhibitor and enzyme are added to a sample containing L-Trp substrate. After a room temperature incubation, activity is determined by measuring the absorption of reaction product at  $\lambda=320 - 325$  nm.

**BACKGROUND:** L-tryptophan (L-Trp) is an essential amino acid necessary for protein synthesis in mammalian cells and the L-Trp to kynurenine (Kyn) pathway is firmly established as a key regulator of innate and adaptive immunity. Catabolism of L-Trp to Kyn maintains an immunosuppressive microenvironment by starving immune cells of L-Trp and releasing degradation products of L-Trp that have immunosuppressive functions. Indoleamine 2,3-dioxygenases (IDO1 & IDO2), two of the rate limiting enzymes in this pathway, are upregulated in many tumors, providing cancer cells with an avenue for immune evasion.

### COMPONENTS:

Catalog #	Component	Amount	Storage	
71194	IDO2 His-Tag	2 x 500 $\mu$ g	-80°C	<b>(Avoid freeze/thaw cycles!)</b>
	IDO2 Reaction Solution component 1	2 x 10 ml	-80°C	
	IDO2 Reaction Solution component 2	2 x 100 $\mu$ l	-80°C	
	IDO2 Substrate	2 x 1 ml	-80°C	
	1x IDO2 Assay Buffer	5 ml	-80°C	
	UV transparent 96-well plate	1	Room Temp.	

### MATERIALS REQUIRED BUT NOT SUPPLIED:

Spectrophotometer capable of measuring absorbance at  $\lambda=320 - 325$  nm.  
Adjustable micropipettor and sterile tips

**APPLICATIONS:** Useful for the study of IDO2 enzymology, screening inhibitors, and selectivity profiling.

### CONTRAINDICATIONS:

DMSO >0.5%, strong acids or bases, ionic detergents, high salt

**STABILITY:** At least 6 months from date of receipt when stored as directed.

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](mailto:info@bpsbioscience.com)

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



6405 Mira Mesa Blvd. Suite 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**REFERENCE(S):**

1. Liu, X., *et al.*, *Blood*. 2010; **115(17)**: 3520-3530.
2. Seegers, N., *et al.* *J. Biomol. Screen.* 2014; **19(9)**: 1266-74.

**ASSAY PROTOCOL:**

All samples and controls should be tested in duplicate. Use slow shaking for all incubations.

**Step 1:**

- 1) Thaw **IDO2 reaction solution component 1** and **component 2**. Prepare **complete IDO2 reaction solution** by mixing 50  $\mu$ l of **component 2** with 4.95 ml of **component 1**. Only prepare enough complete reaction solution needed for assay. *Protect from light, and do not re-use freeze-thawed complete IDO2 reaction solution.*
- 2) Add 50  $\mu$ l of **complete IDO2 reaction solution** to each well.
- 3) Add 5  $\mu$ l of inhibitor solution (containing no more than 10% DMSO in 1x IDO2 assay buffer) to each well designated "Test Inhibitor." For the wells labeled "Positive Control" and "Blank," add 5  $\mu$ l of the same solution without inhibitor (inhibitor buffer). Note: Keep the final DMSO concentration below 0.5%.
- 4) Thaw **IDO2 His-Tag** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Aliquot **IDO2 His-Tag** into single use aliquots. Store remaining undiluted enzyme in aliquots at  $-80^{\circ}\text{C}$ . Note: **IDO2 His-Tag** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.
- 5) Dilute **IDO2 His-Tag** in **1x IDO2 Assay Buffer** to 400 ng/ $\mu$ l. Keep diluted protein on ice until use. Discard any unused diluted protein after use.

	<b>Blank</b>	<b>Positive Control</b>	<b>Test Inhibitor</b>
Complete IDO2 Reaction Solution	50 $\mu$ l	50 $\mu$ l	50 $\mu$ l
Test Inhibitor	–	–	5 $\mu$ l
Inhibitor buffer (no inhibitor)	5 $\mu$ l	5 $\mu$ l	–
1x IDO2 Buffer	25 $\mu$ l	–	–
IDO2 (400 ng/ $\mu$ l)	–	25 $\mu$ l	25 $\mu$ l
IDO2 Substrate	20 $\mu$ l	20 $\mu$ l	20 $\mu$ l
<b>Total</b>	<b>100 <math>\mu</math>l</b>	<b>100 <math>\mu</math>l</b>	<b>100 <math>\mu</math>l</b>

- 6) Add 25  $\mu$ l of **1x IDO2 Assay Buffer** to the well designated "Blank".

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

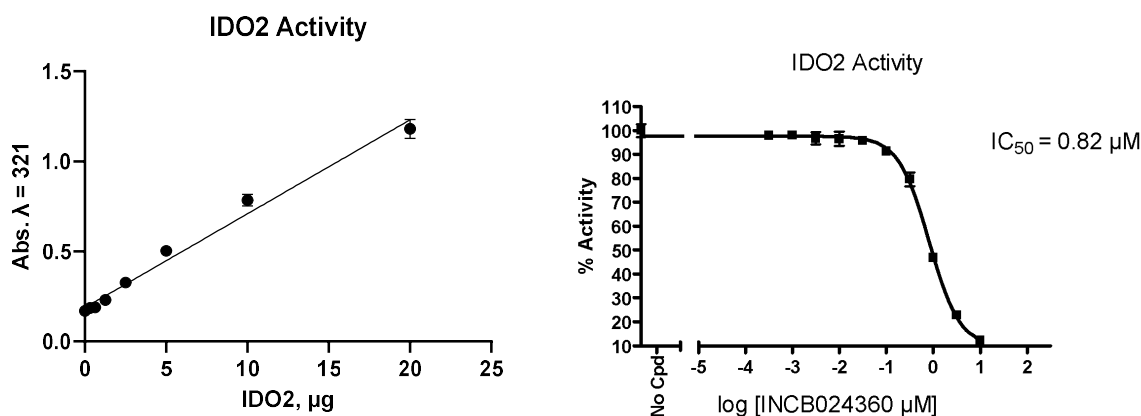
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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

- 7) Add 25  $\mu$ l of diluted **IDO2 His-Tag** prepared as described above to the wells labeled "Positive Control," and "Test Inhibitor." Cover plate with foil and pre-incubate for 30 minutes at room temperature with slow shaking.
- 8) Initiate reaction by adding 20  $\mu$ l of IDO2 substrate. Cover plate with foil and incubate for 2 hours at 30 °C.
- 9) Measure absorption at  $\lambda=320 - 325$  nm.

**EXAMPLE OF ASSAY RESULTS:**



IDO2 activity measured using the IDO2 protocol for the Universal Inhibitor Screening Assay Kit, BPS Bioscience, Catalog #72022. *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto: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](mailto:info@bpsbioscience.com)

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



6405 Mira Mesa Blvd. Suite 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**RELATED PRODUCTS:**

<b><u>Product</u></b>	<b><u>Catalog #</u></b>	<b><u>Size</u></b>
IDO1, His-tag	71182	50 µg
IDO2, His-tag	71194	200 µg
TDO, His-tag	71195	50 µg
IDO1 Inhibitor Screening Assay Kit	72021	96 rxns
TDO Inhibitor Screening Assay Kit	72023	96 rxns
IDO1 Cell-Based Assay Kit	72031	100 rxns
TDO Cell-Based Assay Kit	72033	100 rxns
PD-1[Biotinylated]:PD-L2 Inhibitor Screening Colorimetric Assay Kit	72019	96 rxns
PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit	72003	96 rxns
PD-1:PD-L2[Biotinylated] Inhibitor Screening Assay Kit	72004	96 rxns
PD-1[Biotinylated]:PD-L1 Inhibitor Screening Assay Kit	72005	96 rxns
PD-1[Biotinylated]:PD-L2 Inhibitor Screening Assay Kit	72006	96 rxns
CD28:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72007	96 rxns
BTLA:HVEM[Biotinylated] Inhibitor Screening Assay Kit	72008	96 rxns
CTLA4:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72009	96 rxns
NLG919	27337-1	10 mg
NLG919	27337-2	50 mg
INCB024360	27338-1	10 mg

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](mailto:info@bpsbioscience.com)

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