



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 ***TDO Inhibitor Screening Assay Kit*** Catalog # 72023

**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. Tryptophan 2,3 dioxygenase (TDO) is upregulated in many tumors, providing cancer cells with an avenue for immune evasion.

**DESCRIPTION:** The *TDO Inhibitor Screening Assay Kit* is designed to measure TDO enzyme inhibition. The kit comes in a convenient format, with enough reaction solution and enzyme to perform a total of 100 reactions. The *TDO 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 measured by measuring the absorption of reaction product at  $\lambda = 320-325$  nm.

### COMPONENTS:

| Catalog # | Component                    | Amount     | Storage    |                                    |
|-----------|------------------------------|------------|------------|------------------------------------|
| 71195     | TDO, His-tag*                | 50 $\mu$ g | -80°C      | <b>(Avoid freeze/thaw cycles!)</b> |
| 73005     | TDO Reaction Solution        | 2 x 10 ml  | -80°C      |                                    |
| 73006     | 1x TDO Assay Buffer          | 1 ml       | -80°C      |                                    |
| 79965     | UV transparent 96-well plate | 1          | Room Temp. |                                    |

\*The concentration of TDO is lot-specific and will be indicated on the tube containing the enzyme.

### 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 TDO enzymology, screening inhibitors, and selectivity profiling.

**CONTRAINDICATIONS:** DMSO >0.5%, strong acids or bases, ionic detergents, and high salt

**STABILITY:** Stable at least 6 months from date of receipt, when stored as directed. Kit components require different storage conditions. Be sure to store each component at the proper temperature upon arrival.

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. Li, J.S., *et al. Arch Insect Biochem Physiol.* 2007 Feb; **64(2)**: 74–87

#### ASSAY PROTOCOL:

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

- 1) Thaw **TDO Reaction Solution** and aliquot 180  $\mu$ l into each well. *Note: **TDO Reaction Solution** may contain a precipitate after thawing. Please ensure the mixture is fully solubilized before aliquoting by mixing thoroughly. Do not vortex.*
- 2) Add 10  $\mu$ l of inhibitor solution (no more than 10% DMSO) to each well designated “Test Inhibitor.” For the “Positive Control” and “Blank,” add 10  $\mu$ l of the same solution without inhibitor (inhibitor buffer). *Note: Keep the final DMSO concentration below 0.5%.*
- 3) Thaw **TDO, His-tag** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Aliquot **TDO, His-tag** into single use aliquots. Store remaining undiluted enzyme in aliquots at  $-80^{\circ}\text{C}$ . *Note: **TDO, His-tag** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 4) Dilute **TDO, His-tag** in **1x TDO Assay Buffer** at 50 ng/ $\mu$ l. Keep diluted protein on ice until use. Discard any unused diluted protein after use.

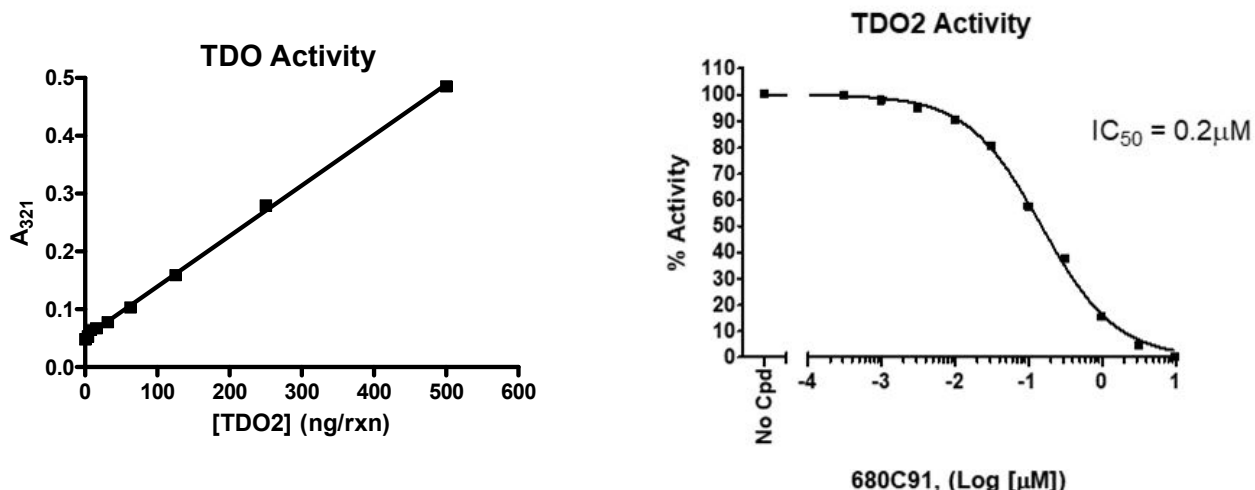
|                                 | Blank                        | Positive Control             | Test Inhibitor               |
|---------------------------------|------------------------------|------------------------------|------------------------------|
| Reaction Solution               | 180 $\mu$ l                  | 180 $\mu$ l                  | 180 $\mu$ l                  |
| Test Inhibitor                  | –                            | –                            | 10 $\mu$ l                   |
| Inhibitor buffer (no inhibitor) | 10 $\mu$ l                   | 10 $\mu$ l                   | –                            |
| 1x TDO Assay Buffer             | 10 $\mu$ l                   | –                            | –                            |
| TDO, His-tag (50 ng/ $\mu$ l)   | –                            | 10 $\mu$ l                   | 10 $\mu$ l                   |
| <b>Total</b>                    | <b>200 <math>\mu</math>l</b> | <b>200 <math>\mu</math>l</b> | <b>200 <math>\mu</math>l</b> |

- 5) Add 10  $\mu$ l of **1x TDO Assay Buffer** to the well designated “Blank.”
- 6) Initiate reaction by adding 10  $\mu$ l of diluted **TDO, His-tag** prepared as described above to the wells labeled “Positive Control,” and “Test Inhibitor.” Incubate at room temperature for 90 minutes.
- 7) Measure absorption at  $\lambda=320-325$  nm.

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)

**EXAMPLE OF ASSAY RESULTS:**



TDO activity (left) and inhibition (right), measured using the TDO Inhibitor Screening Assay Kit, BPS Bioscience, Catalog #72023. *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@bpsbioscience.com).*

**RELATED PRODUCTS:**

| <u>Product</u>  | <u>Catalog #</u> | <u>Size</u> |
|---|------------------|-------------|
| TDO, His-tag  | 71195            | 50 μg       |
| IDO1, His-tag   | 71182            | 50 μg       |
| IDO2, His-tag   | 71194            | 200 μg      |
| IDO1 Inhibitor Screening Assay Kit                            | 72021            | 96 rxns     |
| IDO2 Inhibitor Screening Assay Kit                            | 72022            | 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 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)