



6042 Cornerstone Ct. West, Ste. B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: 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 μ g	-80°C	(Avoid freeze/thaw cycles!)
73005	TDO Reaction Solution	2 x 10 ml	-80°C	
73006	1x TDO Assay Buffer	1 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 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.

REFERENCE(S):

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

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
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Ct. West, Ste. B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

ASSAY PROTOCOL:

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

Step 1:

- 1) Thaw **TDO Reaction Solution** and aliquot 180 μ 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 μ l of inhibitor solution (no more than 10% DMSO) to each well designated "Test Inhibitor." For the "Positive Control" and "Blank," add 10 μ 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°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/ μ l. Keep diluted protein on ice until use. Discard any unused diluted protein after use.

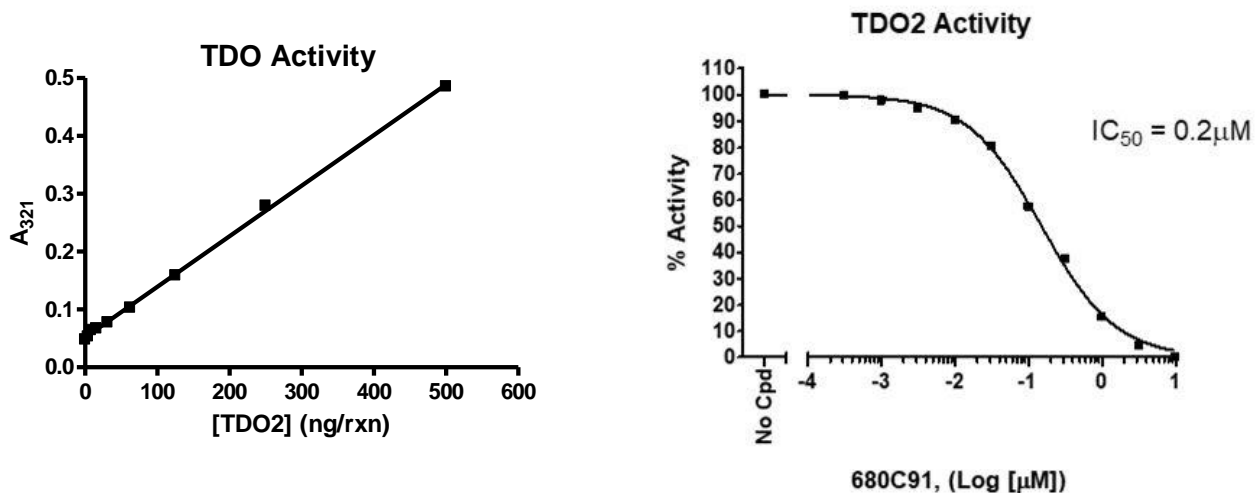
	Blank	Positive Control	Test Inhibitor
Reaction Solution	180 μ l	180 μ l	180 μ l
Test Inhibitor	-	-	10 μ l
Inhibitor buffer (no inhibitor)	10 μ l	10 μ l	-
1x TDO Assay Buffer	10 μ l	-	-
TDO, His-tag (50 ng/ μ l)	-	10 μ l	10 μ l
Total	200 μl	200 μl	200 μl

- 5) Add 10 μ l of **1x TDO Assay Buffer** to the well designated "Blank."
- 6) Initiate reaction by adding 10 μ 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
Please visit our website at: 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.

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
 Please visit our website at: www.bpsbioscience.com