

Data Sheet

IDO1 Fluorogenic Inhibitor Screening Assay Kit **(384 Well Format)** **Catalog #72047**

DESCRIPTION: The *IDO1 Fluorogenic Inhibitor Screening Assay Kit* is designed to measure enzyme inhibition of indoleamine 2,3-dioxygenase 1 (IDO1). The kit comes in a convenient format, with enough reaction solution and enzyme to perform a total of 400 reactions. The *IDO1 Fluorogenic Inhibitor Screening Assay Kit* is simple to use and detects fluorescence at long wavelengths, which minimizes potential errors due to compound interference. In the assay, the inhibitor and enzyme are added to a sample containing L-Trp substrate. After a 1 hour incubation at room temperature, the fluorescence solution is added and incubated at 37°C for four hours. Activity is measured by reading sample fluorescence at $\lambda=510$ nm following excitation of the reaction product at $\lambda=400$ 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	
71182	IDO1 His-Tag	2 x 80 μ g	-80°C	(Avoid freeze/thaw cycles!)
73009	IDO1 Fluorogenic Reaction Solution	3 x 10 ml	-80°C	
73002	IDO1 Assay Buffer	5 ml	-80°C	
	Fluorescence Solution	10 ml	-80°C	
	Black 384 Well Assay-Plate	1		
	Plate sealing film	1		

MATERIALS REQUIRED BUT NOT SUPPLIED:

Fluorimeter capable of excitation at 390-410 nm and detection at 500-520 nm
Adjustable micropipettor and sterile tips
Rotating or rocker platform

APPLICATIONS: Useful for the study of IDO1 enzymology, inhibitor screening, and selectivity profiling.

STABILITY: At least one year 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 **1.858.829.3082**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com

CONTRAINDICATIONS: DMSO >0.5%, strong acids or bases, ionic detergents, high salt.
Warning: the Fluorescence Solution contains a component that is known to be a skin and eye irritant. Use caution and appropriate personal protective equipment when handling this component.

REFERENCE(S):

1. 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 reaction solution and aliquot 50 µl into each well.
- 2) Add 5 µl of inhibitor solution (no more than 10% DMSO) to each well designated "Test Inhibitor." For the wells designated "Positive Control" and "Blank," add 5 µl of the same solution without inhibitor (inhibitor buffer). Note: Keep the final DMSO concentration below 0.5%.
- 3) Dilute **IDO1 His-Tag** in **IDO1 Assay Buffer** at 80 ng/µl. Keep diluted protein on ice until use. Discard any unused diluted protein after use.

	Blank	Positive Control	Test Inhibitor
IDO1 Fluorogenic Reaction Solution	50 µl	50 µl	50 µl
Test Inhibitor	–	–	5 µl
Inhibitor buffer (no inhibitor)	5 µl	5 µl	–
IDO1 Assay Buffer	5 µl	–	–
IDO1 His-Tag (80 ng/µl)	–	5 µl	5 µl
Total	60 µl	60 µl	60 µl

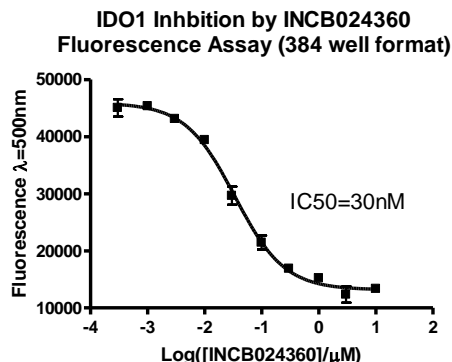
- 4) Add 5 µl of **IDO Assay Buffer** to the wells designated "Blank."
- 5) Initiate reaction by adding 5 µl of diluted **IDO1 His-Tag** prepared as described above to the wells labeled "Positive Control," and "Test Inhibitor." Incubate at room temperature for 1 hour.
- 6) Add 20 µl **Fluorescence Solution** to each well. Seal the plate and incubate at 37°C for four hours. Allow plate to cool for at least 10 minutes.
- 7) Unseal the plate. For best results, centrifuge plate to reduce condensation on the film. Inspect plate for wells with large changes in volume. Wells with significant changes should

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

be excluded from data analysis. Measure fluorescence in a fluorimeter capable of excitation at 400 nm and emission at 510 nm. Subtract "Blank" value from all other values.

EXAMPLE OF ASSAY RESULTS:



Inhibition of IDO1 activity by INCB024360 (Cat. #27338), measured using the *IDO1 Fluorogenic Inhibitor Screening Assay Kit (384 well format)*, BPS Bioscience, Catalog #72047. 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>
IDO1, His-tag	71182	50 μg
IDO2, His-tag	71194	50 μg
TDO, His-tag	71195	50 μg
IDO1 Inhibitor Screening Assay Kit	72021	96 rxns
IDO2 Inhibitor Screening Assay Kit	72022	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
IDO1-HEK293 Recombinant Cell line	60532	2 vials
IDO1 Cellular Activity QuickDetect™ Supplements	62000-1	100 rxns
N-formylkynurenine	73000	2 mg
NLG919	27337-1	10 mg
INCB024360	27338-1	10 mg
IDO1 Reaction Solution	73001	10 ml
IDO1 Assay Buffer	73002	1 ml
IDO2 Reaction Solution	73003	10 ml
IDO2 Assay Buffer	73004	1 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