



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Data Sheet**

### **IDO1 - HEK293 Recombinant Cell Line**

#### **Cat #: 60532**

**Description:** Stable Recombinant HEK293 cell line expressing tetracycline-inducible human indoleamine 2,3-dioxygenase (IDO1) Genbank accession number NM\_002164.

**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.

#### **Application**

- Monitor IDO1 pathway activity
- Screen for activators or inhibitors of IDO1 in a cellular context

#### **Format**

Each vial contains  $\sim 1.5 \times 10^6$  cells in 1 ml of 10% DMSO

#### **Storage**

Immediately upon receipt, store in liquid nitrogen.

#### **Mycoplasma testing**

The cell line has been screened using the PCR-based Venor<sup>®</sup>GeM Mycoplasma Detection kit (Sigma-Aldrich) to confirm the absence of Mycoplasma species.

#### **Materials Required but Not Supplied**

- Thaw Medium 1 (BPS Cat. #60187)
- Growth Medium 1D (BPS Cat. #79536)
- Geneticin (Sigma, #G8168) and Blasticidin S (Sigma (#15205)
- Doxycycline (MP Biomedicals #0219504401)
- PBS
- 0.05% Trypsin EDTA
- 6.1 N Trichloroacetic acid (Sigma #T0699)
- Acetic acid (Sigma #320099)
- IDO1 Cellular Activity QuickDetect<sup>™</sup> Supplements (BPS Cat. #62000). *Note: other formulations can be used, but significant optimization may be required.*

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](mailto:info@bpsbioscience.com)

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



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

### General Culture Conditions

**Thaw Medium 1 (BPS Cat. #60187):** MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS (Invitrogen #26140-079), 1% non-essential amino acids (Hyclone #SH30238.01), 1 mM Na pyruvate (Hyclone #SH30239.01), 1% Penicillin/Streptomycin (Hyclone SV30010.01).

**Growth Medium 1D (BPS Cat. #79536):** Thaw Medium 1 (BPS Cat. #60187) plus 600 µg/ml of Geneticin (Life Technologies #11811031), and 5 µg/ml of Blastidicin (Life Technologies # R210-01) to ensure the recombinant expression plasmid is maintained.

Cells should be grown at 37°C with 5% CO<sub>2</sub> Growth Medium 1D. hIDO1-HEK293 cells should exhibit a typical cell division time of ~50 hours.

**To thaw the cells,** it is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath, transfer to a tube containing 10 ml of Thaw Medium 1 (**no Geneticin or Blastidicin**). Spin down cells, resuspend cells in pre-warmed Thaw Medium 1 (**no Geneticin or Blastidicin**) transfer the resuspended cells to a T25 flask and culture in 37°C in a CO<sub>2</sub> incubator overnight. The next day, replace the medium with fresh Thaw Medium 1 (**no Geneticin or Blastidicin**) and continue growing culture in a CO<sub>2</sub> incubator at 37°C until the cells are ready to be split. Cells should reach ~80% confluence two days after being thawed. Cells should be split before they reach complete confluence. At first passage, switch to Growth Medium 1D (**contains Geneticin and Blastidicin**).

**To passage the cells,** rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with 0.05% Trypsin/EDTA, add Growth Medium 1D (**contains Geneticin and Blastidicin**) and transfer to a tube. Spin down cells, resuspend cells and seed appropriate aliquots of cell suspension into new culture vessels. Subcultivation ratio: 1:5 to 1:10 weekly or twice a week.

### Induction of the target protein expression

Induce cells in MEM medium, 10% FBS, 1% non-essential amino acids, 1 mM Na pyruvate, 1% Penicillin/Streptomycin, 0.2 µg/ml Doxycycline for 24 hours before cell harvesting or assay.

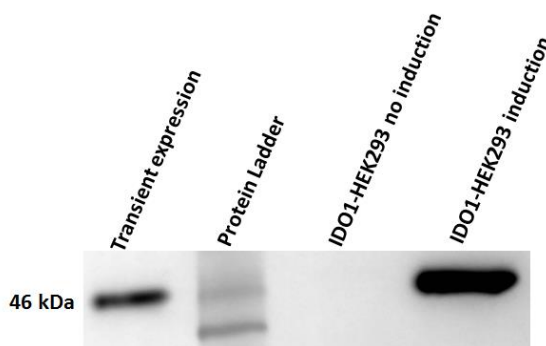
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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](mailto:info@bpsbioscience.com)

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

**Figure 1 Western blot of hIDO1 in hIDO1-HEK293 cells.**



Data shows the tetracycline-inducible expression of hIDO1 in HEK293 cells. Western blot was probed with anti-FLAG antibody (Sigma # F7425).

### Functional validation

N<sup>-</sup>-terminal FLAG tagged human IDO1 has been stably integrated into HEK293 cells and its expression can be induced by tetracycline (doxycycline). The tetracycline-inducible expression of IDO1 was confirmed by Western blotting. IDO1 activity was confirmed by an absorbance-based assay measuring the catalyzed production of kynurenine in cell culture medium.

When IDO1 is expressed in hIDO1-HEK293 cells, it catalyzes L-Trp conversion to Kyn, which gets released in the assay medium and can be easily detected by a reaction with Ehrlich's reagent, producing a yellow color. The hIDO1 enzymatic activity in hIDO1-HEK293 cells can be blocked by a known hIDO1 inhibitor, INCB024360, as shown by the drop in the absorbance signal to the basal level found in the un-induced HEK293 cells.

### Sample protocol to determine the effect of reference inhibitor INCB024360 on doxycycline-induced hIDO1 in hIDO1-HEK293 cells:

*Note: We recommend each treatment be set up in at least triplicate.*

- 1) On day 0, seed hIDO-HEK293 cells at a density of 30,000 cells in 100 µl of Thaw Medium 1 (**no Geneticin or Blastidin**) into each well of a tissue culture-treated 96-well plate. Incubate cells at 37°C in a CO<sub>2</sub> incubator overnight. Leave a couple wells empty for use as a background control.
- 2) The next day (Day 1), treat the cells with 0.2 µg/ml of Doxycycline in Thaw Medium 1 without antibiotics to induce IDO1 expression.
- 3) On the next day (Day 2), prepare *Assay Medium* according to the protocol provided with BPS Cat. #62000. Briefly, after thawing, dilute Assay Supplement 1 1:100 and Assay Supplement 2 1:100 into cell culture medium without antibiotics.

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](mailto:info@bpsbioscience.com)

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



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

- 4) Remove culture medium and treat cells with the test inhibitor in *Assay Medium*, in this case, we used 1  $\mu$ M INCB024360 in 200  $\mu$ l of freshly prepared *Assay Medium*. Add 200  $\mu$ l of *Assay Medium* containing DMSO to cell-free control wells (for determining background absorbance) and un-induced cell control wells (as an optional negative control on any basal level or leaking expression of IDO1 from hIDO1-HEK293 cells). Incubate cells overnight at 37°C in a CO<sub>2</sub> incubator. *Note: The final DMSO concentration should not exceed 0.3%.*
- 5) On day 3, remove 140  $\mu$ l of medium from each well of the cell culture and transfer into a fresh 96-well plate. Add 10  $\mu$ l of 6.1 N trichloroacetic acid to each well. Incubate the plate at 50°C for 30 min. Centrifuge the plate at 2500 rpm for 10 minutes to remove any sediment. If a plate centrifuge is not available, the liquid can be transferred to a microcentrifuge tube and spun briefly to pellet any solids.
- 6) During the incubation, prepare *Detection Reagent Solution* by dissolving Detection Reagent (Provided in BPS Cat. #62000) at a 50-fold dilution in acetic acid, e.g. 200 mg in 10 ml undiluted acetic acid. Prepare only enough reagent required for the assay.
- 7) Transfer 100  $\mu$ l of supernatant to a transparent 96-well plate and mix with 100  $\mu$ l of freshly prepared *Detection Reagent Solution*. Incubate the plate at room temperature for 10 minutes, then measure absorbance at 480 nm using a microplate reader.
- 8) Data analysis: in the absence of the reference inhibitor the absorbance (At) in each should be set to 100%. The absorbance of cell-free control wells (Ab) in each data set should be defined as 0%. The percent absorbance in the presence of reference inhibitor compound is calculated according to the following equation: % Absorbance = (A-Ab)/(At-Ab), where A= the absorbance in the presence of the compound.

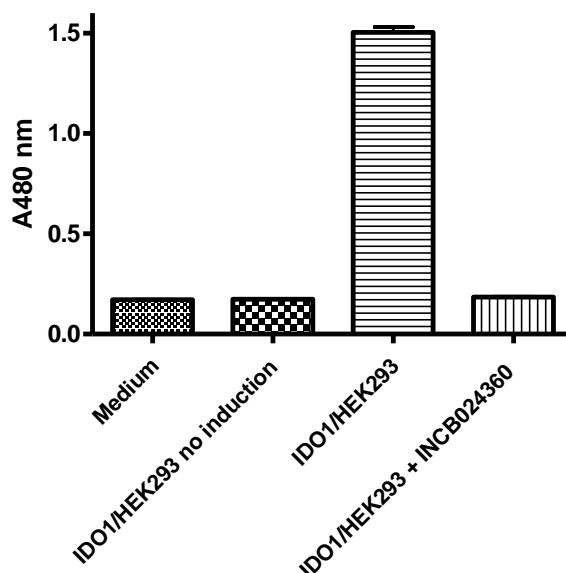
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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](mailto:info@bpsbioscience.com)

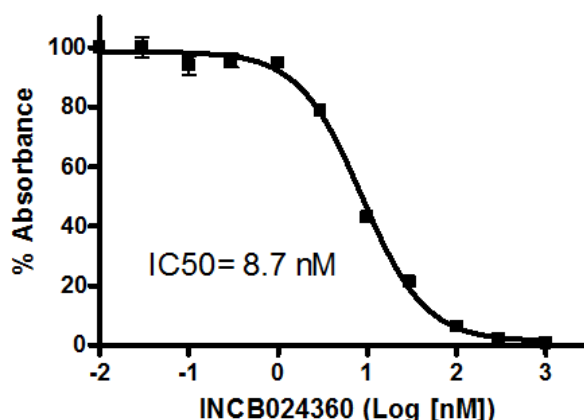
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

**Figure 1. hIDO1-catalyzed Kyn production from L-Trp in hIDO1 – HEK293 Recombinant Cell Line (BPS Cat. #60532).**



INCB024360 completely blocks hIDO1 enzyme activity at a concentration of 1  $\mu$ M. The results are shown as raw absorbance data at 480 nm. Conditions from left to right: medium only (no cells), hIDO1 – HEK293 Cells with no induction plus all assay components, hIDO1 – HEK293 Cells with induction plus all assay components, hIDO1 – HEK293 Cells with induction plus all assay components and INCB024360.

**Figure 2. Dose response of hIDO1 activity in hIDO1-HEK293 cells to reference inhibitor INCB024360.**



The results are shown as percentage of absorbance. The normalized absorbance for hIDO1 expressed cells without inhibitor treatment was set at 100%. The  $IC_{50}$  of INCB024360 is ~ 8.7 nM.

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

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



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

### Vector and sequence

N-terminal FLAG-tagged human IDO1 (accession number NM\_002164) was cloned into a tetracycline-regulated expression vector.

Polylinker: CMV-tetracycline operator (x 2)-EcoRI-FLAG-IDO1-XhoI---SV40-neomycin<sup>R</sup>

hIDO1 sequence (accession number NM\_002164)

MDYKDDDDKAHAMENSWTISKEYHIDEVGFALPNPQENLPDFYNDWMFIAKHLPLDIESGQL  
RERVEKLNMLSIDHLTDHKSQRLARLVLCITMAYVWGKGHGDVRKVLPRNIAVPYCQLSKKL  
ELPPILVYADCVLNWKKKDPNKLPTYENMDVLFSDRGDCSKGFFLVSLLEIAAASAIKVIPTV  
FKAMQMQRDRTLKALLEIASCLEKALQVFHQIHDVNPKAFFSVLRILSGWKGPNQLSDGLV  
YEGFWEDPKFAGGSAGQSSVFQCFDVLLGIQQTAGGGHAAQFLQDMRRYMPPAHRNFLCS  
LESNPSVREFVLSKGDAGLREAYDACVKALVSLRSYHLQIVTKYLIPASQQPKENKTSSEDPKSL  
EAKGTGGTDLMNFKTVRSTTEKSLLKEG

### References

1. Yue, E., *et al.*, *J. Med. Chem.* 2009; **52**: 7364–7367.
2. Liu, X., *et al.*, *Blood*. 2010; **115(17)**: 3520-3530.

### License Disclosure:

Purchase of this cell line grants you with a 10-year license to use this cell line in your immediate laboratory, for research use only. This license does not permit you to share, distribute, sell, sublicense, or otherwise make the cell line available for use to other laboratories, departments, research institutions, hospitals, universities, or biotech companies. The license does not permit the use of this cell line in humans or for therapeutic or drug use. Inappropriate use or distribution of this cell line will result in revocation of the license and result in an immediate cease of sales and distribution of BPS products to your laboratory. Publications using this cell line should reference BPS Bioscience, Inc., San Diego.

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](mailto:info@bpsbioscience.com)

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