



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Data Sheet**

### **hTDO-HEK293 Recombinant Cell line**

### **Catalog: #60534**

**Description:** Stable recombinant HEK293 cell line expressing human tryptophan 2,3 dioxygenase (TDO2, TDO, TPH2, TRPO), Genbank accession number NM\_005651.1.

**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.

#### **Applications**

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

#### **Format**

Each vial contains  $1.75 \times 10^6$  cells in 1 ml of 10% DMSO

#### **Storage**

Immediately upon receipt, store in liquid nitrogen.

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

- Culture Medium: Thaw Medium 1 (BPS Cat. #60187)
- Growth Medium 1B (BPS Cat. #79531)
- Geneticin
- PBS
- 0.05% Trypsin EDTA
- 6.1 N Trichloroacetic acid (Sigma #T0699)
- Acetic acid (Sigma #320099)
- TDO Cellular Activity QuickDetect™ Supplements (BPS Cat. #62002). Note: other formulations can be used, but significant optimization may be required.

#### **Mycoplasma testing**

The cell line has been screened using the PCR-based VenorGeM Mycoplasma Detection kit (Sigma-Aldrich) to confirm the absence of Mycoplasma species.

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)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**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 1B (BPS Cat. #79531):** Thaw Medium 1 (BPS Cat. #60187) plus 400 µg/ml of Geneticin (Life Technologies #11811031) to ensure that recombinant expression is maintained. hTDO-HEK293 cells should exhibit a typical cell division time of ~33 hours.

Cells should be grown at 37°C with 5% CO<sub>2</sub> using Growth Medium 1B.

**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**). Spin down cells, resuspend cells in pre-warmed Thaw Medium 1 (**no Geneticin**), and 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**), 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 1B (**contains Geneticin**)

**To passage the cells,** rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with 0.05% Trypsin/EDTA, add Growth Medium 1B and transfer to a tube. Spin down cells, resuspend cells and seed appropriate aliquots of cell suspension into new culture vessels. Subcultivation ratio: 1:10 to 1:20 weekly or twice a week.

**To freeze down the cells,** rinse cells with phosphate buffered saline (PBS), and detach cells from culture vessel with 0.05% Trypsin/EDTA. Add Thaw Medium 1 and transfer to a tube, spin down cells, and resuspend in freezing medium (10% DMSO + 90% FBS). Place at -80°C overnight and place in liquid nitrogen the next day.

### Functional validation

Human TDO has been stably expressed in HEK293 cells and its activity was characterized by an absorbance-based assay to measure the conversion of L-Trp to Kyn in cell culture medium.

When hTDO is constitutively expressed in hTDO-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 which results in production of a yellow color. The hTDO enzymatic activity in hTDO-HEK293 cells is blocked by a known hTDO inhibitor, 680C91, as shown by the drop in the absorbance signal relative to the basal level in the parental HEK293 cells.

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)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Sample protocol to determine the effect of reference inhibitor 680C91 on constitutively expressed hTDO in hTDO-HEK293 cells:**

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

- 1) On day 1, seed hTDO-HEK293 cells at a density of 30,000 cells in 100 µl of Thaw Medium 1 into each well of a tissue culture-treated 96-well plate. Incubate cells at 37°C in a CO<sub>2</sub> incubator overnight.
- 2) On the next day, prepare *Assay Medium* according to the protocol provided with BPS Cat. #62002. Briefly, after thawing, dilute Assay Supplement 1 1:50 and Assay Supplement 2 1:100 into cell culture medium.
- 3) Remove culture medium and treat cells with the test inhibitor in *Assay Medium*, in this case, we used 10 µM 680C91 in 200 µl of freshly prepared *Assay Medium*. Add 200 µl of Culture Medium containing DMSO to cell-free control wells (for determining background absorbance) and parental HEK293 cell control wells, if desired (as a negative control for any basal level of TDO from 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%.*
- 4) On day 3, remove 140 µl of medium from each well of the cell culture and transfer into a new 96-well plate. Add 10 µl of 6.1 N trichloroacetic acid to each well. Incubate the plate at 50°C for 30 minutes. 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.
- 5) During the incubation, prepare *Detection Reagent Solution* by dissolving Detection Reagent (provided with BPS Cat. #62002) 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.
- 6) Transfer 100 µl of supernatant to a transparent 96-well plate and mix with 100 µl 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.
- 7) Data analysis: in the absence of the reference inhibitor the absorbance (A<sub>t</sub>) in each should be set to 100%. The absorbance of cell-free control wells (A<sub>b</sub>) 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-A<sub>b</sub>)/(A<sub>t</sub>-A<sub>b</sub>), where A= the absorbance in the presence of the compound.

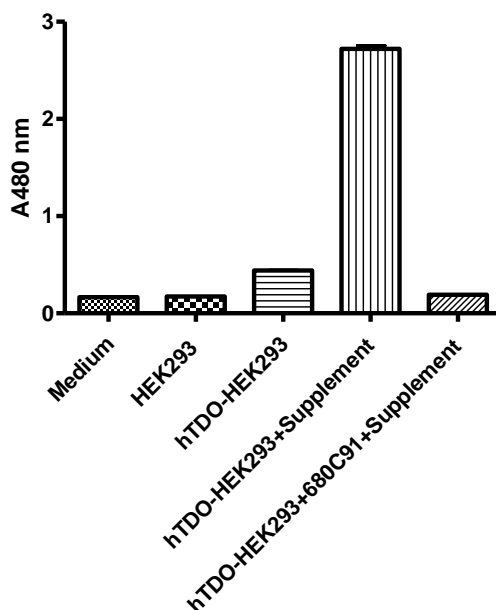
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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

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

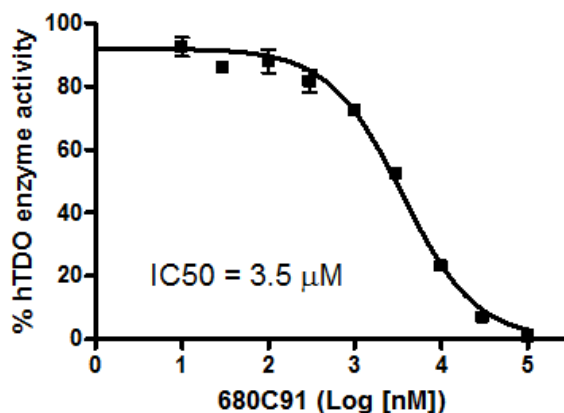
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

**Figure 1. hTDO-catalyzed Kyn production from L-Trp in hTDO-HEK293 Recombinant Cell Line (BPS Cat. #60534).**



680C91 completely blocks hTDO enzyme activity at a concentration of 100  $\mu$ M. The results are shown as raw absorbance data at 480 nm. Conditions from left to right: medium only (no cells), parental HEK293 cells plus all assay components, hTDO-HEK293 plus growth medium (no assay components), hTDO-HEK293 plus all assay components, and hTDO-HEK293 plus all assay components and 100  $\mu$ M 680C91.

**Figure 2. Dose response of hTDO activity to reference inhibitor 680C91 in hTDO-HEK293 cells.**



The results are shown as percentage of absorbance. The normalized absorbance for hTDO transfected cells without inhibitor treatment was set at 100%. The  $IC_{50}$  of 680C91 is ~ 3.5  $\mu$ M

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)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

#### **Vector and sequence**

N-terminal FLAG-tagged human TDO (accession number NM\_005651) was cloned into pcDNA3.1 vector (LifeTechnology).

Polylinker: CMV-HindIII-KpnI-BamHI-FLAG-**TDO**-XhoI-XbaI-ApaI-----SV40-neomycin<sup>R</sup>

Flag-hTDO sequence (accession number NM\_005651)

MDYKDDDDKSGCPFLGNNFGYTFFKKLPVEGSEEDKSQGTGVNRASKGGLIYGNYLHLEKVLNA  
QELQSETKGNKIHDEHLFIITHQAYELWFKQLWELDSVREIFQNGHVRDERNMLKVVSRRMHRV  
SVILKLLVQQFSILETMTALDFNDFREYLSPASGFQSLQFRLLNKIGVLQNMVRPYNRRHYRD  
NFKGEENELLLKSEQEKTLLELVEAWLERTPGLEPHGFNFWGKLEKNITRGLLEEFIRIQAKEES  
EEKEEQVAEFQKQKEVLLSLFDEKRHEHLLSKGERRLSYRALQGALMIYFYREEPRFQVPFQLL  
TSLMDIDSLMTKWRYNHVCMVHRMLGSKAGTGGSSGYHYLRSTVSDRYKVFVDLNLSTYLIP  
RHWIPKMNPTIHKFLYTAEYCDSSYFSSDESD

#### **References**

1. 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 **1.858.202.1401** 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)