



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: info@bpsbioscience.com

## Data Sheet

### TDO Cellular Activity QuickDetect™ Supplements Catalog: #62002-1

**Description:** The TDO Cellular Activity QuickDetect™ Supplements are designed to complement the hTDO-HEK293 Recombinant Cell Line (BPS Cat. #60534) or other TDO-expressing cell lines. This kit contains the cell culture medium supplements necessary for activation of TDO and for the analysis and detection of tryptophan 2,3 dioxygenase (TDO)-catalyzed conversion of L-tryptophan (L-Trp) to Kynurenine (Kyn). The supplements and the detection reagents, when used as described, allow for indirect measurement of Kyn levels by analyzing absorption at 480 nm.

**Background:** L-Trp is an essential amino acid necessary for protein synthesis in mammalian cells, and the L-Trp to 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. 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

Component	Amount	Storage
TDO Assay Medium Supplement 1	400 µl	4°C
TDO Assay Medium Supplement 2	200 µl	-20°C
Detection Reagent	200 mg	Room Temp.

#### Materials Required but Not Supplied

hTDO-HEK293 Recombinant Cell Line (BPS Cat. #60534) or other TDO-expressing cell line and appropriate cell culture medium  
680C91 (Sigma Aldrich #SML0287) or other TDO inhibitor  
6.1 N (concentrated) trichloroacetic acid\*  
17.4 N (concentrated) glacial acetic acid\*

*\*Note: both trichloroacetic acid and acetic acid are strongly corrosive acids; please use gloves and appropriate protective clothing.*

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 Assay Procedure

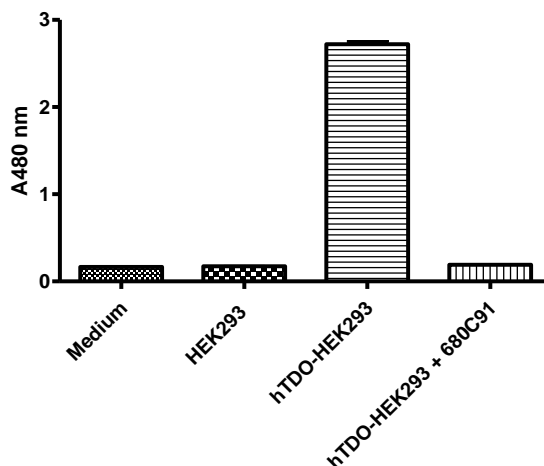
- 1) On day 1, seed cells at a density of 30,000 cells in 100  $\mu$ l of growth medium *without antibiotic* into each well of a tissue culture-treated 96-well plate. Incubate cells at 37°C in a CO<sub>2</sub> incubator overnight.
- 2) Prepare fresh *Assay Medium* by diluting **TDO Assay Medium Supplement 1** 1:50 and **TDO Assay Medium Supplement 2** 1:100 into cell culture medium.
- 3) Dilute the 680C91 (100  $\mu$ M) or test compound to the desired concentration in fresh *Assay Medium*. Remove the cell culture medium from transfected cells and replace with 200  $\mu$ l of *Assay Medium* containing the inhibitor compound. Incubate cells overnight at 37°C in a CO<sub>2</sub> incubator. *Note: The final concentration of DMSO in the cell culture should not exceed 0.3%.*
- 4) On the next day, 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.
- 5) Prepare *Detection Reagent Solution* by dissolving **Detection Reagent** 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  $\mu$ l of supernatant to a transparent 96-well plate and mix with 100  $\mu$ l of fresh *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: The total absorbance (At), without inhibitor treatment, 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. 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 all assay components, hTDO-HEK293 plus all assay components and 100  $\mu$ M 680C91.

**Reference**

1. Liu, X., *et al.*, *Blood*. 2010; **115(17)**: 3520-3530.

**Related Products**

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
hTDO-HEK293 Recombinant Cell Line	60534	2 vials
TDO Cell-Based Assay Kit	72033	100 rxns
IDO1 Cell-Based Assay Kit	72031	100 rxns
TDO Inhibitor Screening Assay Kit	72023	96 rxns
IDO1 Inhibitor Screening Assay Kit	72021	96 rxns
IDO2 Inhibitor Screening Assay Kit	72022	96 rxns
TDO, His-tag	71195	50 $\mu$ g
IDO1, His-tag	71182	50 $\mu$ g
IDO2, His-tag	71194	200 $\mu$ g

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)