



6044 Cornerstone Court West, Suite 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

### **TDO Reaction Solution**

**Catalog #:** 73005

**Lot #:** 150616

**DESCRIPTION:** TDO Reaction Solution is an optimized, proprietary formulation containing all the necessary reaction components for TDO-catalyzed conversion of L-tryptophan (L-Trp) to kynurenine (Kyn) in *in vitro* enzyme activity assays. This product is intended to be used with the BPS TDO Inhibitor Screening Assay Kit (BPS Cat. #72023) and/or with purified, active TDO, His-tag (BPS Cat. #71195) in combination with 1x TDO Assay Buffer (BPS Cat. #73006).

**APPLICATIONS:** Studying enzyme kinetics and HTS applications in drug discovery.

**QUANTITY:** 10 ml

**STORAGE:** -80 °C. Avoid freeze/thaw cycles.

**STABILITY:** Stable for at least one year from date of receipt when stored as directed.

**INSTRUCTIONS FOR USE:** For a detailed protocol on the use of this product, please see the protocol for BPS Cat. #72023.

#### **RELATED PRODUCTS:**

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
TDO Inhibitor Screening Assay Kit	72023	100 reactions
1x TDO Assay Buffer	73006	1 ml
TDO, His-tag	71195	50 µg
TDO Cell-based Assay Kit	72033	100 reactions
hTDO – HEK293 Cell Line	xxxxx	2 vials
N-Formylkynurenine	73000	2 mg

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

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