



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

### ***IDO1 Fluorogenic Reaction Solution***

**Catalog #:** 73009

**Lot #:** 150924

**DESCRIPTION:** IDO1 Fluorogenic Reaction Solution is an optimized, proprietary formulation containing all the necessary reaction components for IDO1-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 Human IDO1 Fluorogenic Inhibitor Screening Assay Kit (BPS Cat. #72037 & Cat. #72047) and/or with purified, active IDO1, His-tag (BPS Cat. #71182) in combination with 1x IDO1 Assay Buffer (BPS Cat. #73002).

**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. #72037 & Cat. #72047.

#### **RELATED PRODUCTS:**

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
Human IDO1 Fluorogenic Inhibitor Screening Assay Kit	72037	96 reactions
Human IDO1 Fluorogenic Inhibitor Screening Assay Kit	72047	384 wells
1x IDO1 Assay Buffer	73002	1 ml
IDO1, His-tag	71182	50 µg
IDO1 Cell-based Assay Kit	72031	100 reactions
INCB024360	27338-1	10 mg
INCB024360	27338-2	100 mg
NLG919	27337	10 mg
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)