



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 Assay Buffer***

**Catalog #:** 73002

**Lot #:** 150708

**DESCRIPTION:** IDO1 Assay Buffer is an optimized, proprietary formulation designed for diluting IDO1 (BPS Cat. #71182) for *in vitro* enzyme activity assays. This product is intended to be used with the BPS IDO1 Inhibitor Screening Assay Kit (BPS Cat. #72021) and/or with purified, active IDO1, His-tag (BPS Cat. #71182) in combination with IDO1 Reaction Solution (BPS Cat. #73001).

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

**QUANTITY:** 1 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. #72021.

#### **RELATED PRODUCTS:**

| <b><u>Product Name</u></b>         | <b><u>Catalog #</u></b> | <b><u>Size</u></b> |
|------------------------------------|-------------------------|--------------------|
| IDO1 Inhibitor Screening Assay Kit | 72021                   | 100 reactions      |
| IDO1 Reaction Solution             | 73001                   | 10 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)