



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

Data Sheet

TDO Fluorogenic Reaction Solution

Catalog #: 73010

Lot #: 150924

DESCRIPTION: TDO Fluorogenic 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 Human TDO Fluorogenic Inhibitor Screening Assay Kits (BPS Cat. #72039 & Cat. #72049) 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. #72039 & Cat. #72049.

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
Human TDO Fluorogenic Inhibitor Screening Assay Kit	72039	96 reactions
Human TDO Fluorogenic Inhibitor Screening Assay Kit	72049	384 wells
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

Please visit our website at: www.bpsbioscience.com