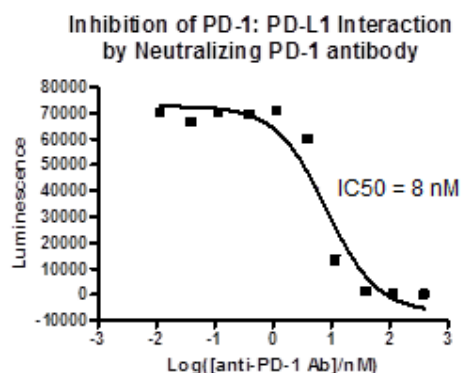


# Immunotherapy Screening Services

Immunotherapy is defined as the treatment of diseases by inducing, enhancing, or suppressing the immune system. Since the immune system plays a role in virtually all human diseases and immune modulators have shown clinical efficacy, immunotherapy is becoming an increasingly prominent area of pharmaceutical focus. Most of these therapeutics target a variety of immune checkpoint pathways, such as the PD-1:PD-L1/L2 signaling axis, to either enhance or inhibit immune function.

BPS Bioscience's Immunotherapy Screening Services allow you to screen for inhibitors of any of the following targets in our biochemical assays:

Biochemical Assays	
○ BTLA:HVEM	○ IL-17A: IL-17RA
○ CD28:B7-1	○ OX40: OX40L
○ CD47:Sirpa	○ PD-1:PD-L1
○ CTLA4:B7-1	○ PD-1:PD-L2
○ COX1	○ PD-1: B7-1
○ COX2	○ TIGIT: CD155
○ IDO-1	○ TPH1
○ IDO-2	○ TPH2



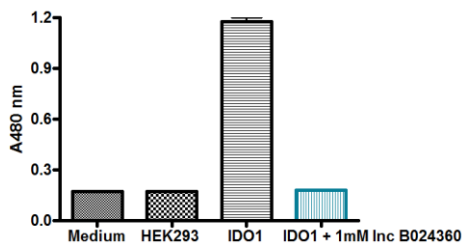
**Figure 1.** Inhibition of PD-1:PD-L1 binding by PD-1 Neutralizing Antibody (BPS Catalog #71120), measured with BPS's PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit, Catalog #72003.

- Fast turn-around time – typically under 2 weeks
- Determine IC50 values (10 point, duplicate) or screen a single concentration over our full panel of Immunotherapy targets
- Fully customizable to meet your needs
- Choice of dose range and concentrations

## Cell-Based Assays

BPS also offers cell-based assays for the popular IDO1 and TDO drug targets. These enzymes catalyze the first and rate-limiting step in tryptophan catabolism down the kynurenine pathway. Their activity and expression has been found to be up-regulated in numerous cancers.

Let BPS' experts help you screen for IDO1 and/or TDO inhibitors in our cell-based assays.



**Figure 2:** Demonstration of IDO1 activity and IDO1 inhibition in BPS' IDO1 Cell-Based Screening Assay. Conditions from left to right: medium only (no cells), HEK293 cells plus all assay components, IDO1-expressing HEK293 plus all assay components, IDO1-expressing HEK293 plus all assay components and 1  $\mu$ M INCB024360.

*We also offer custom assays for other immunotherapy targets!*

Contact us for a free quotation or information on our Immunotherapy Screening Services by visiting:

[www.bpsbioscience.com/contact-us](http://www.bpsbioscience.com/contact-us)

\*To review sample data for the Immunotherapy assays listed above, click on the following link or copy and paste it into your web browser: [http://www.bpsbioscience.com/images/pdf/Immunotherapy\\_sample\\_data.pdf](http://www.bpsbioscience.com/images/pdf/Immunotherapy_sample_data.pdf)

BPS Bioscience, Inc. • 6044 Cornerstone Court West, Suite E, San Diego, CA 92121

Tel: 858.829.3082 • Fax: 858.481.8694 • Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com) • Website: [www.bpsbioscience.com](http://www.bpsbioscience.com)