



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet
TPH1 Inhibitor Screening Assay Kit
Catalog: #72056
Size: 384 reactions

BACKGROUND: Tryptophan 5-hydroxylase 1 and 2 (TPH1 and TPH2) are enzymes that catalyze the monooxygenation of tryptophan to 5-hydroxytryptophan (5-HTP) which is subsequently converted to serotonin. By catalyzing the rate-limiting step in the biosynthesis of serotonin, these enzymes also play key roles in the signaling of other neurotransmitters downstream of serotonin synthesis, like melatonin. Additionally, TPH enzymes may play a role in immune regulation through tryptophan depletion.

DESCRIPTION: The TPH1 Inhibitor Screening Assay Kit is designed to measure enzyme inhibition in a 384 reaction format. This fluorescence based assay kit is especially suitable for high throughput screening applications. The procedure is straightforward and simple, requiring only the mixing of two solutions, incubation, addition of quench solution, and measurement of fluorescence.

COMPONENTS:

| Catalog # | Component | Amount | Storage | |
|-----------|-----------------------|-----------|-----------|---|
| 71192 | TPH1 | 2 x 50 µg | -80°C | (Avoid freeze/ thaw cycles!) |
| | TPH Enzyme Solution | 10 ml | -80°C | |
| | TPH Reaction Solution | 12 ml | -80°C | |
| | TPH Quench Solution | 3 x 1 ml | -20°C | |
| | Black, 384 Well Plate | 1 | Room Temp | |

MATERIALS REQUIRED BUT NOT SUPPLIED:

Microplate reader capable of measuring Fluorescence

APPLICATIONS: Great for screening small molecular inhibitors for drug discovery and HTS applications.

STABILITY: At least 6 months from date of receipt when stored as directed.

REFERENCE:

1. Moran G. R., et al. J. Biol. Chem. 1998; 273:12259-12266.
2. Nowak E. C., et al., J. Exp. Med. 2012; 209:2127-2135.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
 San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 1) Thaw **TPH Enzyme Solution** on ice. Take only a sufficient quantity needed for the assay. Keep solution on ice. Store remaining solution in aliquots at -80°C.
- 2) Thaw **TPH Reaction Solution** on ice. Take only a sufficient quantity needed for the assay. Keep solution on ice. Store remaining solution in aliquots at -80°C.
- 3) Keep plate cold on ice. Add 5 µl of inhibitor solution to each well labeled “Test Inhibitor”. Add 5 µl of the same solution without inhibitor (Inhibitor Buffer) to each well labeled “Negative Control” and “Positive Control”.

| | Negative Control* | Positive Control | Test Inhibitor |
|---------------------------------|-------------------|------------------|----------------|
| Test Inhibitor | — | — | 5 µl |
| Inhibitor Buffer (no inhibitor) | 5 µl | 5 µl | — |
| Enzyme Solution | 20 µl | — | — |
| TPH1 Solution (11 ng/µl) | — | 20 µl | 20 µl |
| Reaction Solution | 25 µl | 25 µl | 25 µl |
| Total | 50 µl | 50 µl | 50 µl |

- 4) Add 20 µl of **TPH Enzyme Solution** to each well labeled “Negative Control”.
- 5) Thaw **TPH1** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Aliquot **TPH1** into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C. Note: **TPH1** is very sensitive to freeze/thaw cycles. Do not reuse thawed aliquots or diluted enzyme.
- 6) Dilute **TPH1** to 11 ng/µl (220 ng/reaction) in **TPH Enzyme Solution**. Keep on ice until ready to use.
- 7) Add 20 µl of diluted **TPH1** to each well labeled “Positive Control” and “Test inhibitor”.
- 8) Initiate reaction by adding 25 µl of **TPH Reaction Solution** to each well. Remove plate from ice and incubate for 4 hours at 4°C.
- 9) After incubation, add 5 µl of **TPH Quench Solution** to each well. After quenching, plate can be handled at room temperature.
- 10) Read the fluorescent intensity in a microtiter-plate reader.

Instrument Settings

| | |
|-----------------------|--------|
| Reading Mode | λ (nm) |
| Excitation Wavelength | 300 |
| Emission Wavelength | 360 |

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

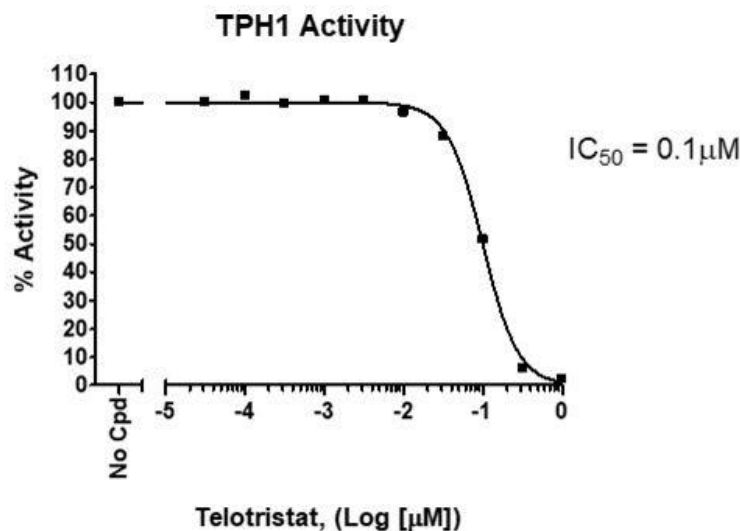


6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

CALCULATING RESULTS:

Subtract average "Negative Control" value from average "Positive Control" value to obtain total Δ Fluorescence. Subtract average "Negative Control" value from each "Test Inhibitor" value to obtain Δ Fluorescence of test compounds.

EXAMPLE OF ASSAY RESULTS:



Inhibition of TPH1 by Telotristat, measured using the *TPH1 Assay Kit*, BPS Bioscience # 72056. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

RELATED PRODUCTS:

| <u>Product</u> | <u>Catalog #</u> | <u>Size</u> |
|--|-------------------------|--------------------|
| TPH1, His-tag | 71192 | 50 µg |
| TPH2, His-tag | 71193 | 50 µg |
| TPH1 Inhibitor Screening Assay Kit | 72053 | 96 rxns. |
| TPH2 Inhibitor Screening Assay Kit | 72054 | 96 rxns. |
| TPH2 Inhibitor Screening Assay Kit | 72057 | 384 rxns. |
| Human TDO, His-tag | 71195 | 50 µg |
| Mouse TDO, His-tag | 71241 | 50 µg |
| hTDO-HEK293 Recombinant Cell line | 60534 | 2 vials |
| TDO Cellular Activity QuickDetect™ Supplements | 62002-1 | 100 rxns. |
| Human TDO Inhibitor Screening Assay Kit | 72023 | 96 rxns. |

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com