

## Data Sheet

### PPID (CYP-40)

Human, recombinant, N-terminal GST-tag

**Catalog #:** 71095

Lot #: 130703      Conc.: 1.87 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8, 110 mM NaCl, 2.2 mM KCl, 16 mM glutathione, and 20% glycerol.

**Stability:** >6 months at -80°C

#### References:

1. Tavecchio, M., *et al.*, *J. Biol. Chem.* 2013;**288(8)**:5553-5561.
2. Ward, B.K., *et al.*, *J. Biol. Chem.* 2002;**277(43)**:40799-40809.

#### Description:

Human Peptidyl-prolyl cis-trans isomerase D (PPID), also known as peptidyl-prolyl cis-trans isomerase Cyclophilin-40 (CYP-40, CyP-D), GenBank Accession No. NM\_005038, a.a. 2-370(end), with N-terminal GST-tag, MW 67.5 kDa, expressed in an *E. coli* cell expression system.

#### Application:

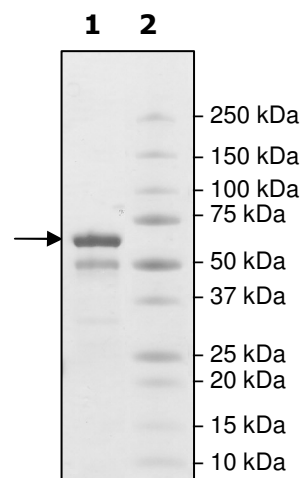
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance

#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
3.7 µg PPID  
**Lane 2:**  
Protein Marker

**MW:** 67.5 kDa  
**Purity:** ≥73%



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)



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)

**Sequence:**

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI  
DGDVKLTQSMARIYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSKDFETLK  
VDFLSKLPPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV  
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDLEVLFQGPLGSSHPSP  
QAKPSNPSNPRVFFDVDIGGERVGRIVLELFADIVPKTAENFRALCTGEKGIGHTTGKP  
LHFKGCPFHRIKKFMIQGGDFSNQNGTGGESIYGEKFEDENFHYKHDREGLLSMANA  
GRNTNGSQFFITTVPTPHLDGKHVVFVGQVIKIGIVARILENVEVKGEKPAKLCVIAECG  
ELKEGDDGGIFPKDGS GDSHPDFPEDADIDLKDVDKILLITEDLKNIGNTFFKSQNWEM  
AIKKYAEVLRVYDSSKAVIETADRAKLQPIALSCVLNIGACKLKMSNWQGAIDSCLEAL  
ELDPSNTKALYRRAQGWQGLKEYDQALADLKKAQGIAPEDKAIQAELLKVKQKIKAQ  
KDKEKAVYAKMFA

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)