

PTPN11/SHP2-CD Phosphatase Assay Service

PTPN11/SHP2 Catalytic Domain [Phosphatase Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

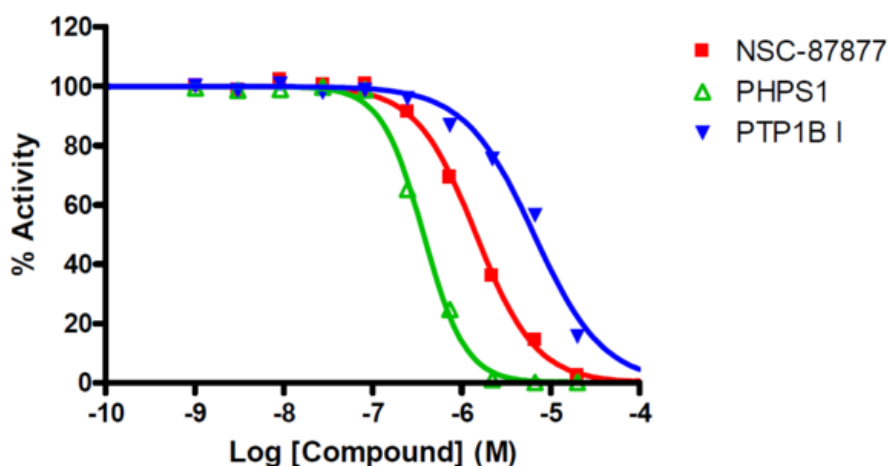
Target Overview

CAT#	PTPN11/SHP2-CD
Enzyme	Catalytic domain of human PTPN11/SHP2 (aa246-593)
Synonyms	Tyrosine-protein phosphatase non-receptor type 11, Protein-tyrosine phosphatase 1D, Protein-tyrosine phosphatase 2C, SH-PTP2, SH-PTP3, PTP2C, SHPTP2
Substrates	Fluorogenic substrate DiFMUP, 10 μ M

Assay Properties

Reaction	Protein tyrosine phosphate + H ₂ O = protein tyrosine + phosphate								
Readout	Enzymatic activity								
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>NSC-87877</td> <td>1,500</td> </tr> <tr> <td>PHPS1</td> <td>370</td> </tr> <tr> <td>PTP1B Inhibitor</td> <td>6,500</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	NSC-87877	1,500	PHPS1	370	PTP1B Inhibitor	6,500
Compounds	IC50 (nM)								
NSC-87877	1,500								
PHPS1	370								
PTP1B Inhibitor	6,500								
Scientific Information	Q06124 (http://www.uniprot.org/uniprot/Q06124)								
Screening Location	Malvern, USA								
Further information	More information can be found on our website Phosphatase Assay Services .								

Reference Compound IC50 for PTPN11/SHP2



Let's Discover Together.

© May 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/ptpn11_shp2_cd_malvern_phos

