

PTPN6/SHP1 Phosphatase Assay Service

PTPN6/SHP1 [Phosphatase Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

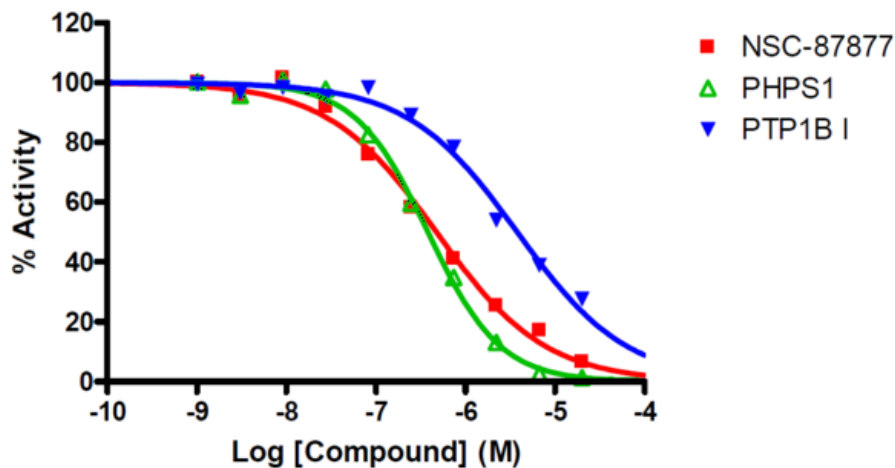
Target Overview

CAT#	PTPN6/SHP1
Enzyme	Human PTPN6/SHP1
Synonyms	Tyrosine-protein phosphatase non-receptor type 6, Hematopoietic cell protein-tyrosine phosphatase Protein-tyrosine phosphatase 1C, Protein-tyrosine phosphatase SHP-1, HCP, PTP1C
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>470</td> </tr> <tr> <td>PHPS1</td> <td>380</td> </tr> <tr> <td>PTP1B Inhibitor</td> <td>3,800</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	NSC-87877	470	PHPS1	380	PTP1B Inhibitor	3,800
Compounds	IC50 (nM)								
NSC-87877	470								
PHPS1	380								
PTP1B Inhibitor	3,800								
Scientific Information	P29350 (http://www.uniprot.org/uniprot/P29350)								
Screening Location	Malvern, USA								

Reference Compound IC50 for PTPN6/SHP1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/ptpn6_shp1_malvern_phos

