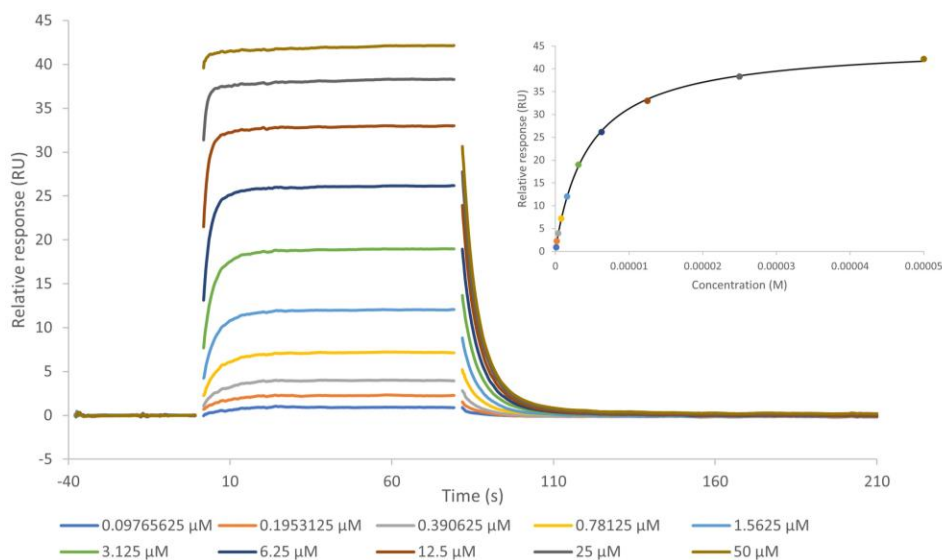


# PNMT SPR Assay Service

PNMT for direct binding screening and profiling of test substance with [surface plasmon resonance \(SPR\)](#) technique.

Target Overview	
Target	PNMT
HGNC Symbol	PNMT
Synonyms	Phenylethanolamine N-methyltransferase
Scientific Information	<a href="https://www.uniprot.org/uniprotkb/P11086">P11086</a> ( <a href="https://www.uniprot.org/uniprotkb/P11086">https://www.uniprot.org/uniprotkb/P11086</a> )
Protein Source	Produced at Reaction Biology (aa 2-282)
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website <a href="#">SPR Assay Services for Drug Discovery</a> .
SAM binding to PNMT	



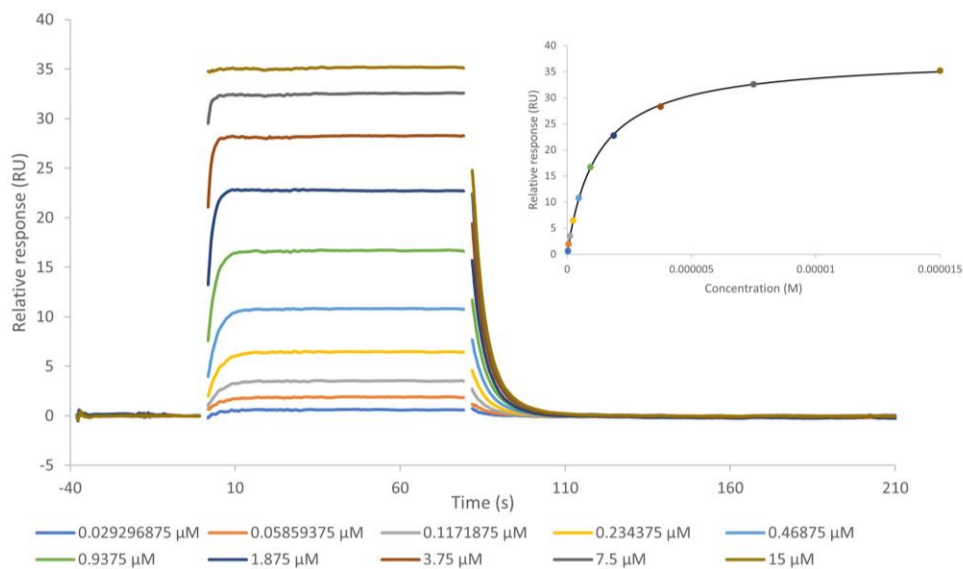
Binding Model	Target	Analyte	Chi <sup>2</sup> (RU <sup>2</sup> )	K <sub>D</sub> (M)	R <sub>max</sub> (RU)	Offset (RU)
Steady state affinity	PNMT	SAH	7.17e-2	1.21e-6	37.7	0.1

The data were collected on a Biacore 8K using multi-cycle kinetics. The data were analyzed using a steady state affinity model.

Let's Discover Together.

© Jan 2024. Reaction Biology. All rights reserved.

## SAH binding to PNMT



Binding Model	Target	Analyte	Chi <sup>2</sup> (RU <sup>2</sup> )	K <sub>D</sub> (M)	R <sub>max</sub> (RU)	Offset (RU)
Steady state affinity	PNMT	SAM	1.17e-1	4.58e-6	45.1	0.4

The data were collected on a Biacore 8K using multi-cycle kinetics. The data were analyzed using a steady state affinity model.