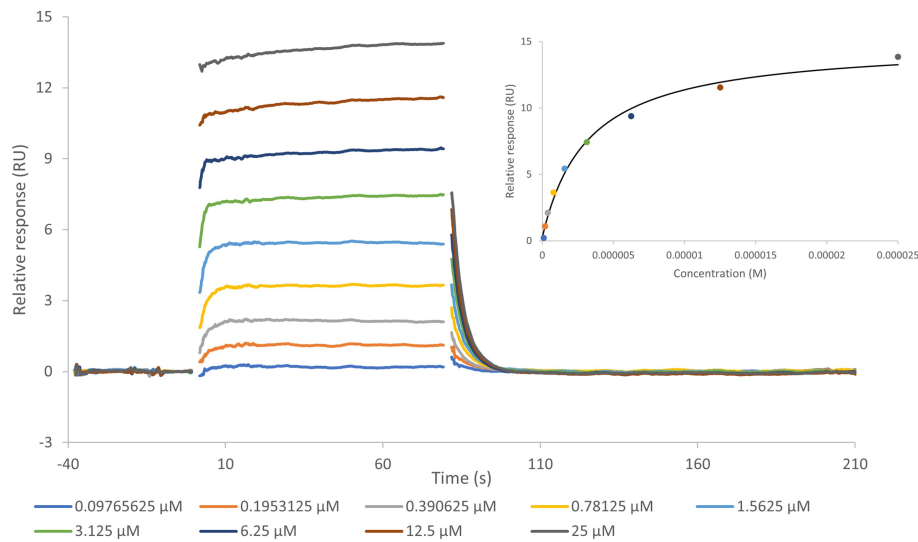


NNMT SPR Assay Service

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

Target Overview	
Target	NNMT
HGNC Symbol	NNMT
Synonyms	Nicotinamide N-methyltransferase
Scientific Information	P40261 (https://www.uniprot.org/uniprotkb/P40261)
Protein Source	Produced at Reaction Biology. Catalog # MT-11-358
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website SPR Assay Services for Drug Discovery .

SAM binding to NNMT



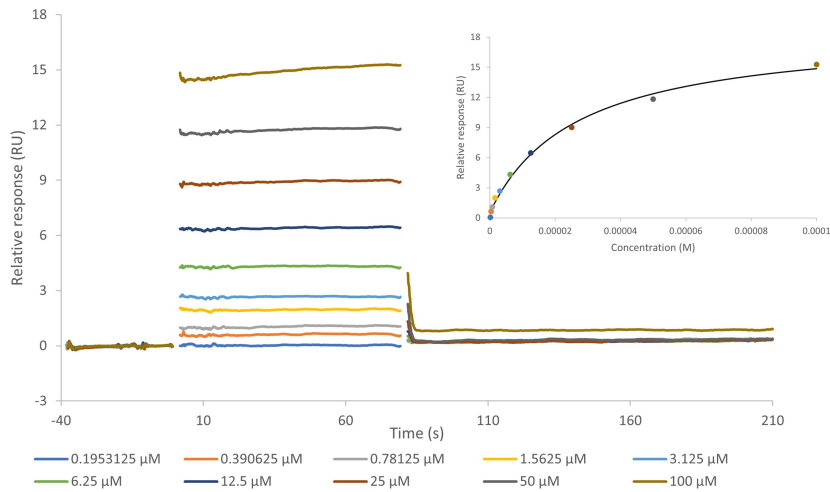
Binding Model	Target	Analyte	Chi ² (RU ²)	K _D (M)	R _{max} (RU)	Offset (RU)
Steady state affinity	NNMT	SAM	2.51e-1	3.21e-6	14.6	0.3

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.

© Mar 2024. Reaction Biology. All rights reserved.

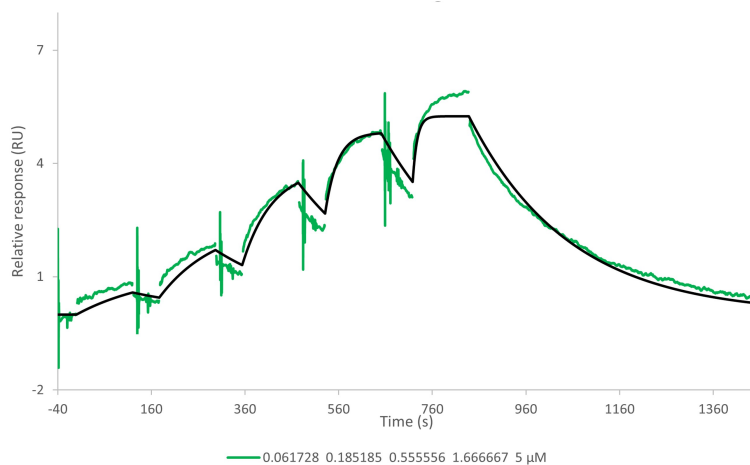
SAH binding to NNMT



Binding Model	Target	Analyte	Chi ² (RU ²)	K _D (M)	R _{max} (RU)	Offset (RU)
Steady state affinity	NNMT	SAH	1.95e-1	2.68e-5	18.1	0.6

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

JBSNF binding to NNMT



Binding Model	Target	Analyte	Chi ² (RU ²)	k _a (1/Ms)	k _d (1/s)	K _D (M)	R _{max} (RU)
1:1 Kinetic	NNMT	JBSNF-000088	7.68e-2	2.01e+4	4.66e-3	2.32e-7	5.5

The data were collected on a Biacore 8K using single-cycle kinetics. The data were analyzed using a 1:1 kinetic binding model.

Let's Discover Together.

© Mar 2024. Reaction Biology. All rights reserved.