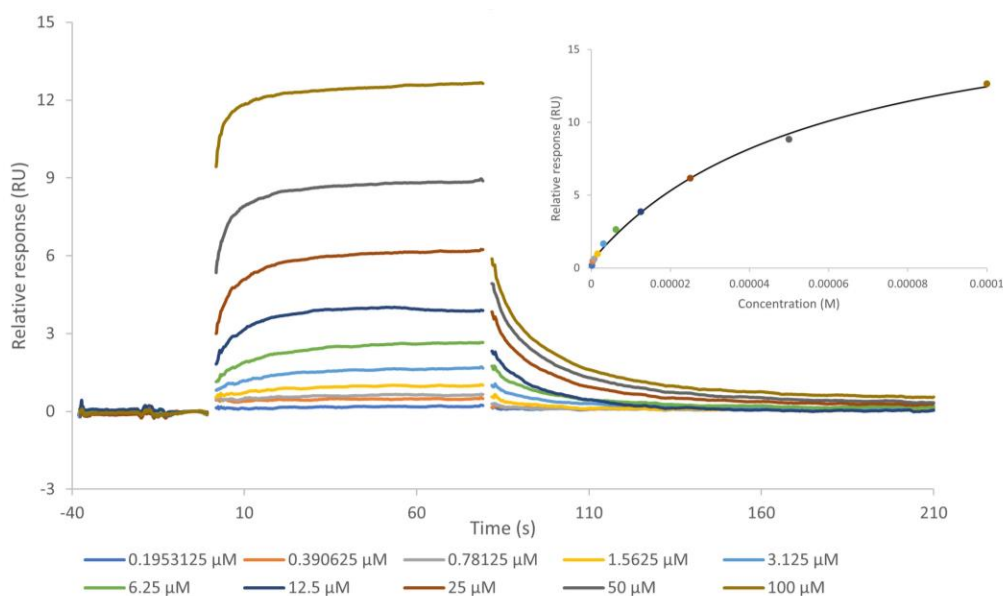


INMT SPR Assay Service

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

Target Overview	
Target	INMT
HGNC Symbol	INMT
Synonyms	Indolethylamine N-methyltransferase
Scientific Information	O95050 (https://www.uniprot.org/uniprotkb/O95050)
Protein Source	Produced at Reaction Biology (aa 2-263)
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website SPR Assay Services for Drug Discovery .
SAM binding to INMT	



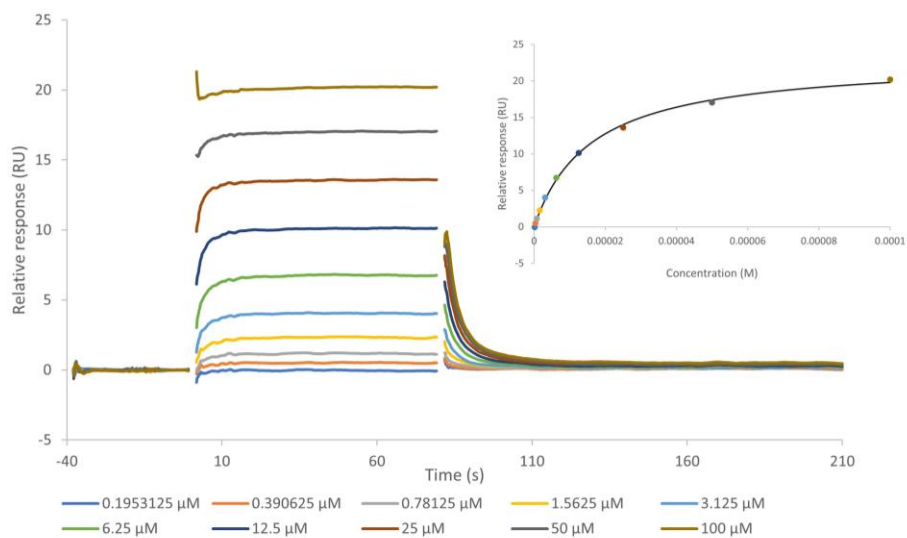
Binding Model	Target	Analyte	Chi ² (RU ²)	K _D (M)	R _{max} (RU)	Offset (RU)
Steady state affinity	INMT	SAM	7.07e-2	5.80e-5	19.0	0.5

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 INMT



Binding Model	Target	Analyte	Chi ² (RU ²)	K _D (M)	R _{max} (RU)	Offset (RU)
Steady state affinity	INMT	SAH	1.07e-1	1.60e-5	22.9	0.0

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