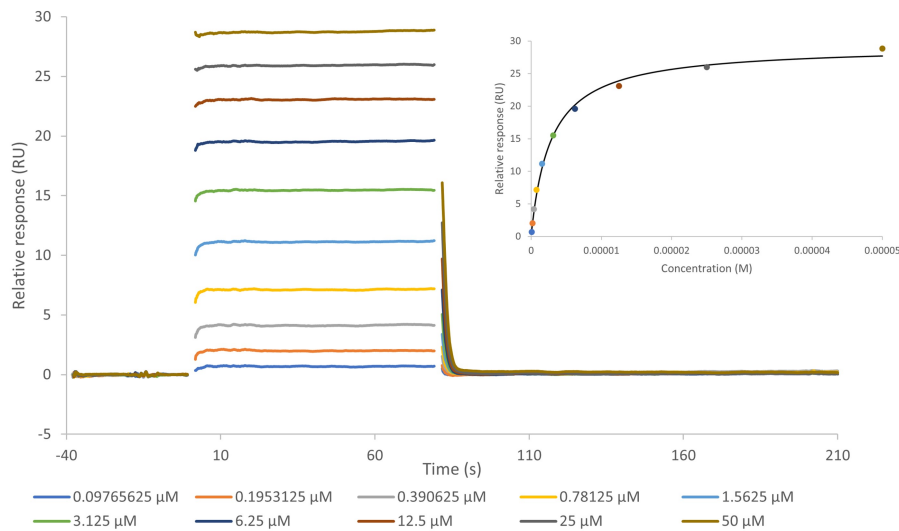


# HNMT SPR Assay Service

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

Target Overview	
Target	HNMT
HGNC Symbol	HNMT
Synonyms	Histamine N-methyltransferase
Scientific Information	<a href="https://www.uniprot.org/uniprotkb/P50135">P50135</a> ( <a href="https://www.uniprot.org/uniprotkb/P50135">https://www.uniprot.org/uniprotkb/P50135</a> )
Protein Source	Produced at Reaction Biology (aa 2-292)
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 HNMT



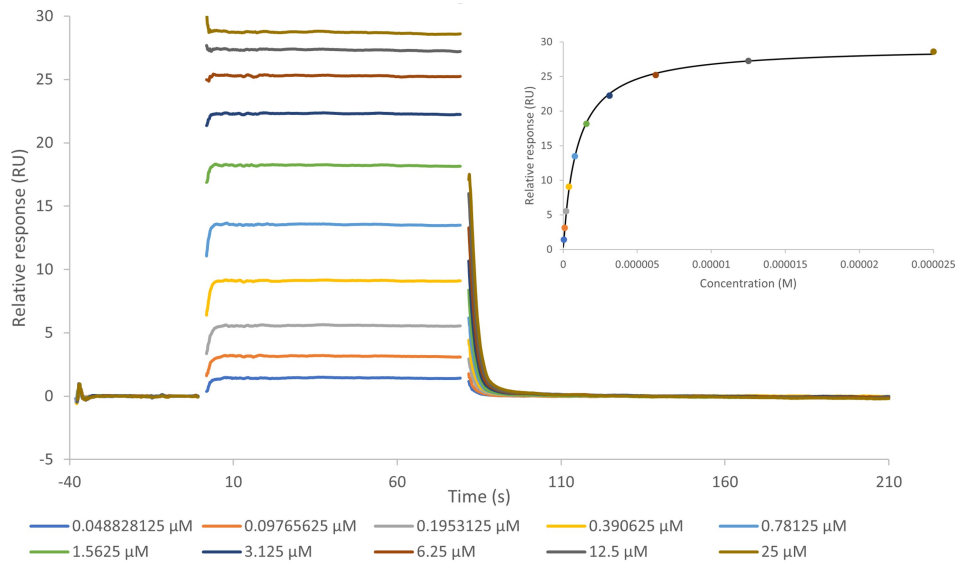
Binding Model	Target	Analyte	Chi <sup>2</sup> (RU <sup>2</sup> )	K <sub>D</sub> (M)	R <sub>max</sub> (RU)	Offset (RU)
Steady state affinity	HNMT	SAM	5.91e-1	2.85e-6	28.9	0.4

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 HNMT



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	HNMT	SAH	8.70e-2	9.53e-7	28.9	0.4

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