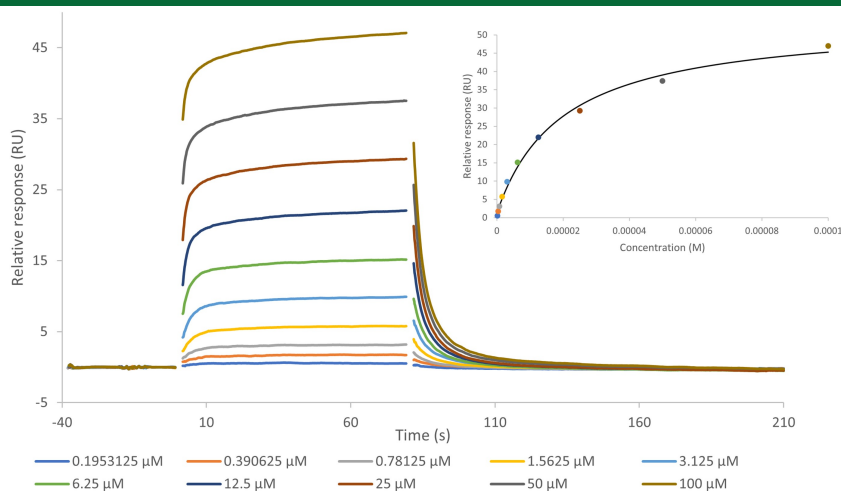


GNMT SPR Assay Service

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

Target Overview	
Target	GNMT
HGNC Symbol	GNMT
Synonyms	Glycine N-methyltransferase
Scientific Information	Q14749 (https://www.uniprot.org/uniprotkb/Q14749)
Protein Source	Produced at Reaction Biology (aa 2-295)
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website SPR Assay Services for Drug Discovery .

SAM binding to GNMT



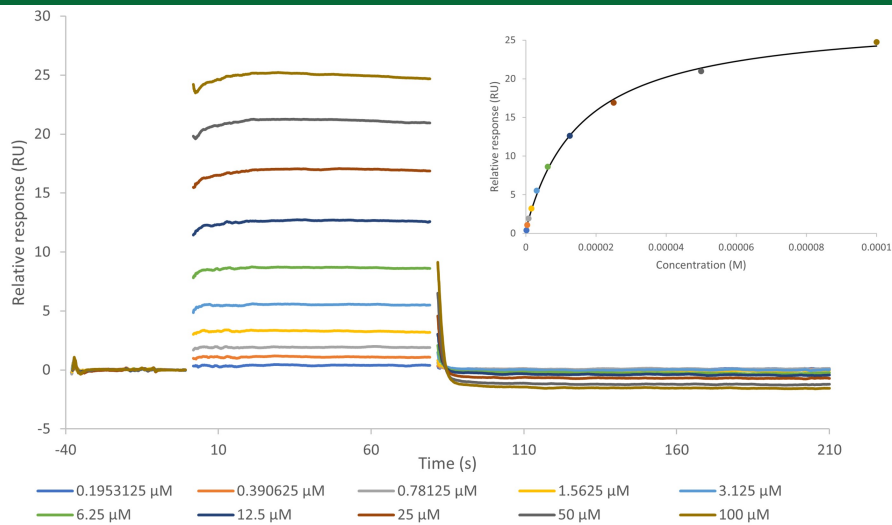
Binding Model	Target	Analyte	Chi ² (RU ²)	K _D (M)	R _{max} (RU)	Offset (RU)
Steady state affinity	GNMT	SAM	1.88e+0	1.99e-5	52.7	1.3

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 GNMT



Binding Model	Target	Analyte	Chi ² (RU ²)	K _D (M)	R _{max} (RU)	Offset (RU)
Steady state affinity	GNMT	SAH	1.72e-1	1.58e-5	27.4	0.5

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