

SOS1 Binding Assay Service (Thermal shift)

KRAS-related Assays: SOS1 binding assay service for compound screening and profiling using the thermal shift assay format to measure compound binding to the target

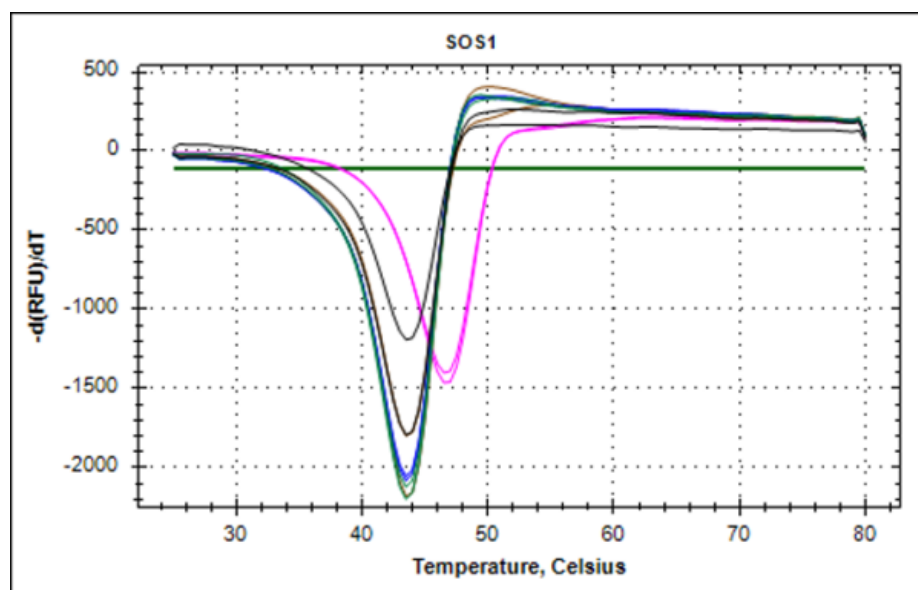
Target Overview

Target	Human SOS1
HGNC	SOS1
Protein	SOS1 (aa 564-1069), strep tagged. Expressed in Ecoli.
Synonyms	Son of sevenless homolog 1, GF1, GGF1, GINGF, HGF, NS4

Assay Properties

Scientific information	SOS1: Q07889 (https://www.uniprot.org/uniprot/Q07889)
Screening location	Malvern, PA, US
Reaction	Compound binding to SOS1
Readout	Protein stability change due to ligand binding via thermal shift assay
Concentration of protein in assay	0.15 mg/ml (~2.5 μ M)
Concentration of compound in assay	25 μ M
Apo form Tm ($^{\circ}$C)	43.5
Reference Compound Tms	BAY-293 +3 (pink) , no stabilization was observed in the presence of other control compounds (ARS-1620, AMG-510, BI-2852)

Melt peak of SOS1 protein at increasing temperatures in presence of compounds BAY-293 (pink), ARS-1620 (green), AMG-510 (blue) or BI-2852 (brown) or none (black)



Let's Discover Together.

