

# SOS2 Binding Assay Service (Thermal shift)

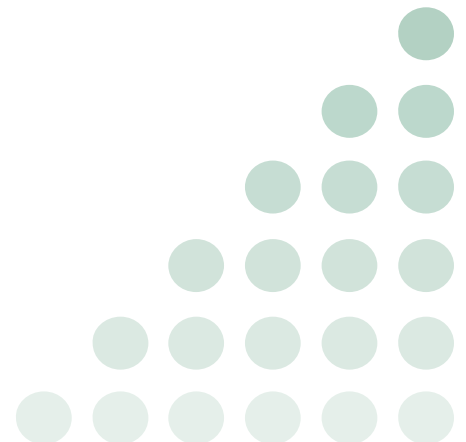
**KRAS-related Assays:** SOS2 binding assay service for compound screening and profiling using the thermal shift assay format to measure compound binding to the target

## Target Overview

<b>Target</b>	Human SOS2
<b>HGNC</b>	SOS2
<b>Protein</b>	SOS2 (aa 562-1047), strep tagged. Expressed in Ecoli.
<b>Synonyms</b>	Son of sevenless homolog 2

## Assay Properties

<b>Scientific information</b>	SOS2: <a href="https://www.uniprot.org/uniprot/Q07890">Q07890</a> ( <a href="https://www.uniprot.org/uniprot/Q07890">https://www.uniprot.org/uniprot/Q07890</a> )
<b>Screening location</b>	Malvern, PA, US
<b>Reaction</b>	Compound binding to SOS2
<b>Readout</b>	Protein stability change due to ligand binding via <a href="#">thermal shift assay</a>
<b>Concentration of protein in assay</b>	Typically tested at 0.1 mg/ml
<b>Concentration of compound in assay</b>	n/a
<b>Apo form Tm (°C)</b>	43.5
<b>Reference Compound Tms</b>	No reference compound is currently available for SOS2. No protein stabilization was observed when tested with reference compounds BAY293, ARS-1620, AMG-510 or BI-2852.



Melt peak of SOS2 protein at increasing temperatures in presence of compounds BAY-293 (pink), ARS-1620 (green), AMG-510 (blue) or BI-2852 (brown) or none (black)

