

# IDH2 wt Isocitrate Dehydrogenase Assay Service

IDH2 wt Isocitrate Dehydrogenase Assay for compound screening and profiling via [Diaphorase/Resazurin-coupled assay](#) for quantification of enzyme activity.

<b>CAT#</b>	IDH2 wt
<b>Enzyme</b>	Human IDH2 wild type
<b>Synonym</b>	Isocitrate dehydrogenase [NADP], mitochondrial (EC:1.1.1.42); ICD-M; Oxalosuccinate decarboxylase)
<b>Substrate</b>	150 $\mu$ M isocitrate , 25 $\mu$ M NADP
<b>Reaction</b>	isocitrate + NADP = a-KG + NADPH
<b>Readout</b>	Enzymatic activity
<b>Reference Compound IC50s</b>	<b>Compounds</b> <b>IC50 (nM)</b> IDH-C100 227
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/P48735">P48735</a> (http://www.uniprot.org/uniprot/P48735 )
<b>Screening Location</b>	Malvern, USA

Reference compound IC50 for IDH2 WT

