

IDH2 (R172Q)

Isocitrate dehydrogenase assay, Diaphorase/Resazurin-coupled assay

CAT#	IDH2 (R172Q)								
Enzyme	Human IDH2 (R172Q) mutant								
Synonym	Isocitrate dehydrogenase [NADP], mitochondrial (EC:1.1.1.42); ICD-M; Oxalosuccinate decarboxylase								
Substrate	50 μ M α -ketoglutarate, 15 μ M NADPH								
Reaction	α KG + NADPH = 2HG + NADP								
Readout	Enzymatic activity								
Reference Compound IC50s	<table border="0"> <tr> <td>Compound</td> <td>IC50 (nM)</td> </tr> <tr> <td>IDH2-C100</td> <td>83</td> </tr> <tr> <td>AG-221</td> <td>19</td> </tr> <tr> <td>AGI-6780</td> <td>10</td> </tr> </table>	Compound	IC50 (nM)	IDH2-C100	83	AG-221	19	AGI-6780	10
Compound	IC50 (nM)								
IDH2-C100	83								
AG-221	19								
AGI-6780	10								
Scientific Information	P48735 (http://www.uniprot.org/uniprot/P48735)								
Screening Location	Malvern, USA								

Reference compound IC50 for IDH2 (R172Q)

