

MTHFD1 Replication-related Assay Service

MTHFD1 [Replication-related Assay](#) for compound screening and profiling via quantification of enzyme activity.

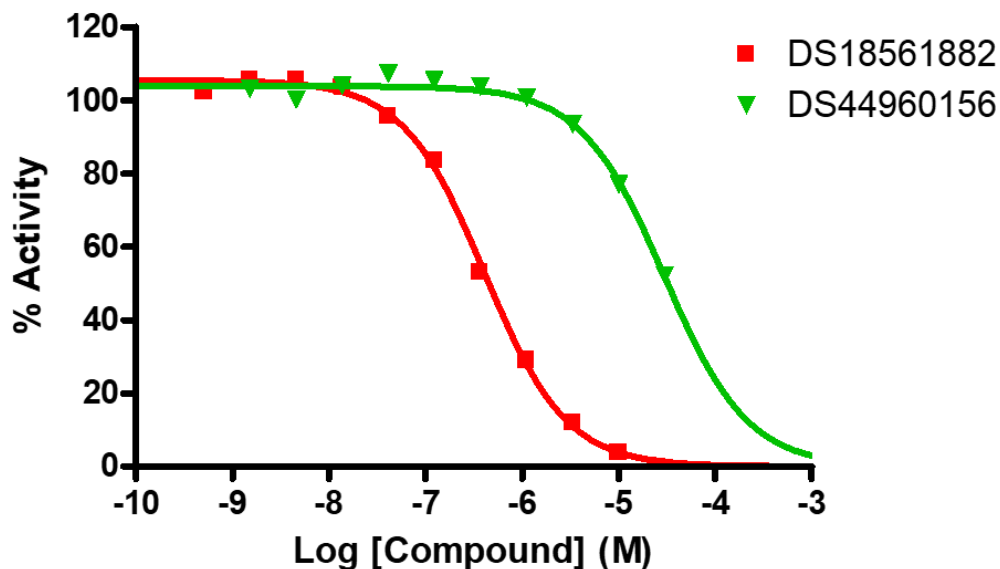
Target Overview

CAT#	MTHFD1
Enzyme	Human C-1-tetrahydrofolate synthase, cytoplasmic
Synonyms	C1-THF synthase, Methylenetetrahydrofolate dehydrogenase, Epididymis secretory sperm binding protein, MTHFC, MTHFD
Substrate	50 μ M Folitixorin (5,10-Methylenetetrahydrofolic Acid), 100 μ M NADP

Assay Properties

Reaction	(6R)-5,10-methylene-5,6,7,8-tetrahydrofolate+ NADP = 5,10-methenyltetrahydrofolate + NADPH End point measurement, Diaphorase/Resazurin coupled detection, Resorufin fluorescence measurement
Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (μM) DS18561882 0.41 DS44960156 29.8
Scientific Information	P11586 (http://www.uniprot.org/uniprot/P11586)
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website Replication-related Assays

Reference compound IC50 for MTHFD1



Let's Discover Together.

© Sep 2023 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/mthfd1_repl_malvern

