

# MTH1 Replication-related Assay Service

MTH1 [Replication-related Assay](#) for compound screening and profiling via quantification of enzyme activity using the PPiLight™ Inorganic Pyrophosphate Assay.

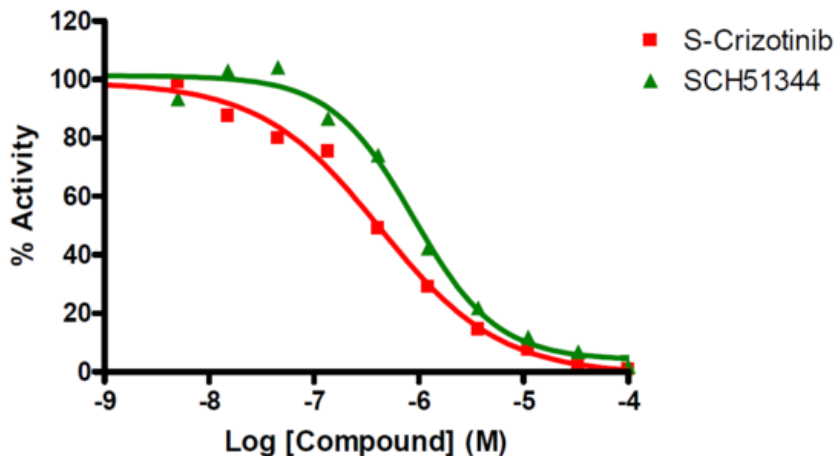
## Target Overview

CAT#	MTH1
Enzyme	Human MTH1
Synonyms	NUDT1, 7,8-dihydro-8-oxoguanine triphosphatase, Nucleoside diphosphate-linked moiety X motif 1
Substrate	15 μM 8-Oxo-2'-dGTP

## Assay Properties

Reaction	8-oxo-dGTP + H <sub>2</sub> O = 8-oxo-dGMP + diphosphate
Readout	Enzymatic activity via PPiLight™ Inorganic Pyrophosphate Assay (Lonza)
Reference Compound IC50s	<b>Compounds IC50 (nM)</b> S-Crizotinib 428 SCH51344 894
Scientific Information	P36639 ( <a href="http://www.uniprot.org/uniprot/P36639">http://www.uniprot.org/uniprot/P36639</a> )

## Reference compound IC50 for MTH1



Let's Discover Together.

© Jul 2022 Reaction Biology. All rights reserved.

[https://www.reactionbiology.com/datasheet/MTH1\\_rep\\_Malvern](https://www.reactionbiology.com/datasheet/MTH1_rep_Malvern)

