

# COMT Methyltransferase Assay Service

Activity-based COMT Methyltransferase Assay for compound screening and profiling via [FlashPlate methyltransferase assay](#).

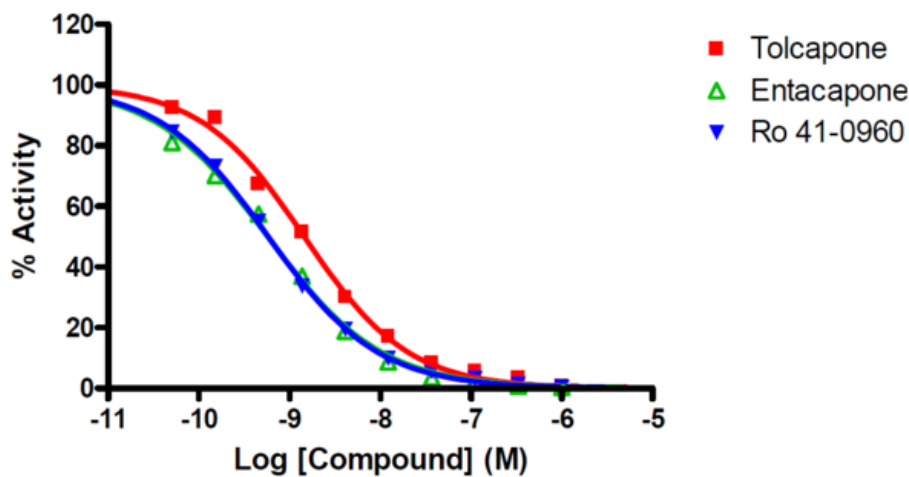
## Target Overview

CAT#	COMT
Enzyme	Human COMT
Synonym	Catechol O-methyltransferase
Substrates	COMT-S01, 0.5 $\mu$ M SAM, 1 $\mu$ M

## Assay Properties

Reaction	S-adenosyl-L-[methyl- $^3$ H]methionine + a catechol = S-adenosyl-L-homocysteine + a guaiacol [methyl- $^3$ H]
Readout	Enzymatic activity
Reference Compound IC50s	<b>Compounds IC50 (nM)</b> Tolcapone 1.5 Entacapone 0.6 Ro 41-0960 0.6
Scientific Information	<a href="http://www.uniprot.org/uniprot/P21964">P21964</a> ( <a href="http://www.uniprot.org/uniprot/P21964">http://www.uniprot.org/uniprot/P21964</a> )
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website <a href="#">Methyltransferase Assays</a> .

## Reference compound IC50 for COMT



Let's Discover Together.

© Oct 2022 **Reaction Biology**. All rights reserved.

[https://www.reactionbiology.com/datasheet/comt\\_methyl\\_malvern](https://www.reactionbiology.com/datasheet/comt_methyl_malvern)

