

CYP1A2 Inhibition Assay Service

CYP1A2 for compound screening and profiling via fluorescence-based quantification of enzyme activity.

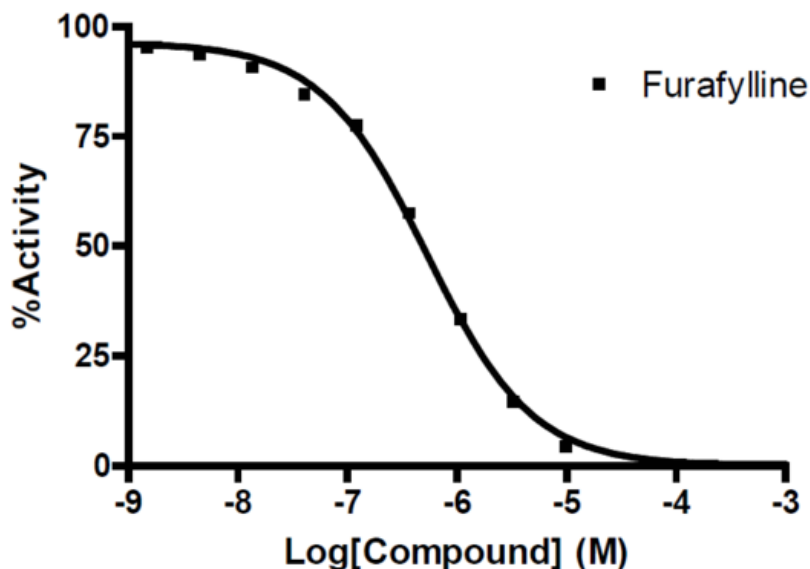
Target Overview

CAT#	Cyp1A2
Enzyme	Cyp1A2 + NADPH-P450 Reductase (microsomes from insect cells that were infected with baculovirus containing the human Cyp1A2 cDNA)
Synonym	cytochrome P450, family 1, subfamily A, polypeptide 2, CYP 1A2
Substrate	Fluorescent substrate

Assay Properties

Reaction	$RH + O_2 + NADPH + H^+ = ROH + H_2O + NADP$	
Readout	Enzymatic activity	
Reference Compound IC50s	Compound	IC50 (μM)
	Ketoconazole	2.05
	Proadifen (SKF525A)	3.15
Scientific Information	NP_000752 (http://www.ncbi.nlm.nih.gov/entrez/viewer.fcgi?db=protein&id=73915100)	
Screening Location	Malvern, USA	
Further Information	More information can be found on our website Cytochrome P450 Assay Services .	

Reference compound IC50 for Cyp1A2



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/cyp1a2_malvern

