

CYP2D6 Inhibition Assay Service

CYP2D6 for compound screening and profiling via fluorescence-based quantification of enzyme activity.

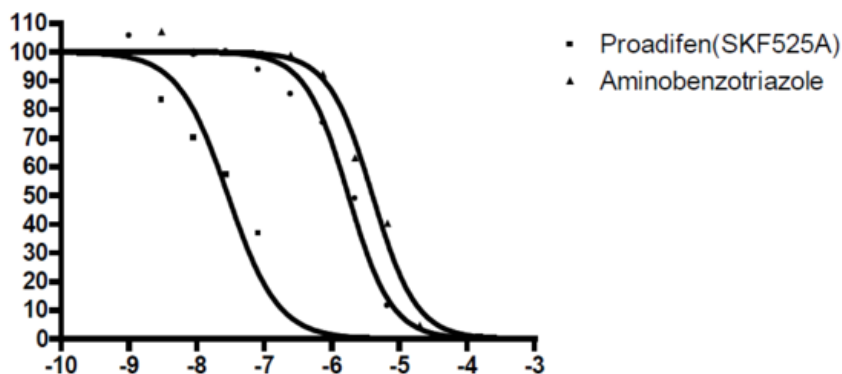
Target Overview

CAT#	Cyp2D6
Enzyme	Cyp2D6 + NADPH-P450 Reductase (microsomes from insect cells that were infected with baculovirus containing the human Cyp2D6 cDNA and rabbit Cytochrome P450 cDNA)
Synonym	Cytochrome P450, family 2, subfamily D, polypeptide 6
Substrate	Fluorescent substrate

Assay Properties

Reaction	$RH + O + NADPH + H = ROH + H_2O + NADP$	
Readout	Enzymatic activity	
Reference Compound IC50s	Compound	IC50 (μM)
	Ketoconazole	1.76
	Proadifen (SKF-525A)	0.03
	Aminobenzotriazole	4.00
Scientific Information	NP_000097 (http://www.ncbi.nlm.nih.gov/entrez/viewer.fcgi?db=protein&val=40805836)	
Screening Location	Malvern, USA	
Further Information	More information can be found on our website Cytochrome P450 Assay Services .	

Reference compound IC50 for Cyp2D6



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/cyp2d6_malvern

