

## Cyp2D6

(Cytochrome P450, family 2, subfamily D, polypeptide 6)

- **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)
- **Substrate:** Vivid EOMCC Substrate
- **Reaction:**  $RH + O_2 + NADPH + H^+ \rightarrow ROH + H_2O + NADP^+$
- 
- **Reference Compound IC50s:**

Compound	IC50 (microM)
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)  
(<http://www.ncbi.nlm.nih.gov/entrez/viewer.fcgi?db=protein&val=40805836>)

IC50s for Control inhibitor for Cyp2D6

