

## Cyp2C9

(Cytochrome P450, family 2, subfamily C, polypeptide 9)

- **CAT #:** Cyp2C9
- **Enzyme:** Cyp2C9 + NADPH-P450 Reductase (Microsomes from insect cells that were infected with baculovirus containing the human Cyp2C9 cDNA and rabbit Cytochrome P450 cDNA)
- **Substrate:** Vivid OOMR Substrate
- **Reaction:**  $\text{RH} + \text{O}_2 + \text{NADPH} + \text{H}^+ \rightarrow \text{ROH} + \text{H}_2\text{O} + \text{NADP}^+$

- **Reference Compound IC50s:**

Compound	IC50(microM)
Ketoconazole	0.71

- **Scientific information:** [AAH70317](#)  
(<http://www.ncbi.nlm.nih.gov/entrez/viewer.fcgi?db=protein&val=47682790>)

### IC50 of Ketoconazole for Cyp2C9

