

## Cyp1A2

(cytochrome P450, family 1, subfamily A, polypeptide 2)

- **CAT #:** Cyp1A2
- **Enzyme:** Cyp1A2 + NADPH-P450 Reductase (Microsomes from insect cells that were infected with baculovirus containing the human Cyp1A2 cDNA)
- **Substrate:** Florescent Substrate
- **Reaction:**  $RH + O_2 + NADPH + H^+ \rightarrow ROH + H_2O + NADP^+$
- **Reference Compound IC50s:**

Compound	IC50(microM)
Ketoconazole	2.05
Proadifen(SKF525A)	3.15

- **Scientific information:** [NP\\_000752](http://www.ncbi.nlm.nih.gov/entrez/viewer.fcgi?db=protein&id=73915100)  
(<http://www.ncbi.nlm.nih.gov/entrez/viewer.fcgi?db=protein&id=73915100>)

### Furafylline IC50 for Cyp1A2

