

## Cyp3A4

(Cytochrome P450, subfamily IIIA, polypeptide 4)

- **CAT #:** Cyp3A4
- **Enzyme:** Cyp3A4 + NADPH-P450 Reductase (Microsomes from insect cells that were infected with baculovirus containing the human Cyp3A4 cDNA and rabbit Cytochrome P450 cDNA)
- **Substrate:** Florescent Substrate
- **Reaction:**  $RH + O_2 + NADPH + H^+ \rightarrow ROH + H_2O + NADP^+$

- **Reference Compound IC50s:**

Compound	IC50(microM)
Ketoconazole	0.04
Proadifen(SKf525A)	0.28
Aminobenzotriazole	6.38

- **Scientific information:** [NP\\_059488](http://www.ncbi.nlm.nih.gov/entrez/viewer.fcgi?db=protein&val=13435386)  
(<http://www.ncbi.nlm.nih.gov/entrez/viewer.fcgi?db=protein&val=13435386>)

### IC50s of Control inhibitors for Cyp3A4

