

Cyp2C19

(Cytochrome P450, family 2, subfamily C, polypeptide 19)

- **CAT #:** Cyp2C19
- **Enzyme:** Cyp2C19 + NADPH-P450 Reductase (Microsomes from insect cells that were infected with baculovirus containing the human Cyp2C19 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	2.05
Proadifen(SKF525A)	3.15

- **Scientific information:** [NP_000760](http://www.ncbi.nlm.nih.gov/entrez/viewer.fcgi?db=protein&val=4503219)
(<http://www.ncbi.nlm.nih.gov/entrez/viewer.fcgi?db=protein&val=4503219>)

IC50s for Control inhibitors for Cyp2C19

