

Antibody Specificity Tables Human P450 Inhibitory Antibodies

The following is a summary of the crossreactivities of antibodies which have applications to the immunoinhibition of human cytochrome P450 enzymes. Antibodies are polyclonal unless otherwise noted.

Human Cytochrome P450														
Anti-serum	1A1	1A2	2A6	2B6	2C8	2C9	2C18	2C19	2D6	2E1	3A4	3A5	OR	Cat. No.
Anti-CYP1A2	-	+++	-	-	-	-	-	-	-	-	-	-	-	458312
Anti-CYP2A6*	-	-	+++	-	-	-	-	-	-	+	-	-	-	458106
Anti-CYP2A6*			+++	-	-	-	-	-	-	-	-	-	-	458306
Anti-CYP2B6*	-	-	-	+++	-	-	-	-	-	-	-	-	-	458326
Anti-CYP2C8*	-	-	-	-	+++	-	-	-	-	-	-	-	-	458328
Anti-CYP2D6*	-	-	-	-	-	-	-	-	+++	-	-	-	-	458366
Anti-CYP2E1*	-	-	-	-	-	-	-	-	-	+++	-	-	-	458321
Anti-CYP3A4*	-	-	-	-	-	-	-	-	-	-	+++	+++	-	458334

*: Monoclonal Antibody
 +++: Very Strongly Crossreactive
 ++: Strongly Crossreactive
 +: Crossreactive
 (+): Weakly Crossreactive
 -: Crossreactivity Not Detected
 lm: Lower Mobility
 hm: Higher Mobility