

Antibody Specificity Tables

Human P450 Immunoblotting Antibodies

The following is a summary of the immunoblotting crossreactivities of antibodies which have applications in the study of cytochrome P450 enzymes in human tissues. Antibodies are polyclonal unless otherwise noted.

Human Cytochrome P450																
Anti-serum	1A1	1A2	1B1	2A6	2B6	2C8	2C9	2C18	2C19	2D6	2E1	3A4	3A5	3A7	4A11	Cat. No.
Anti-CYP1B1	-	-	+++	-	-	-	-	-	-	-	-	-	-	-	-	458211
Anti-CYP2A6*	-	-	-	+++	-	-	-	-	-	-	+lm	-	-	-	-	458106
Anti-CYP2B6	-	-	-	-	+++	-	-	-	-	-	-	-	-	-	-	458226
Anti-CYP2C9	-	-	-	-	-	-	+++	-	-	-	-	-	-	-	-	458209
Anti-CYP2C19	-	(+)	-	(+)	(+)	-	-	-	+++	-	-	-	(+)	-	-	458219
Anti-CYP2D6*	-	-	-	-	-	-	-	-	-	+++	-	-	-	-	-	458246
Anti-CYP3A4*	-	-	-	-	-	-	-	-	-	-	-	+++	+++	+++	-	458254
Anti-CYP3A4	-	-	-	-	-	-	-	-	-	-	-	+++	-	+++	-	458234
Anti-CYP3A5	-	-	-	-	-	-	-	-	-	-	-	-	+++	-	-	458235

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