

Antibody Specificity Tables

Rat P450 Immunoblotting Antibodies

The following is a summary of the immunoblotting crossreactivities of antibodies which have applications to the study of rat cytochrome P450 enzymes in rat tissues. The Antibody below, is an antipeptide antisera developed in rabbits and is a polyclonal antibody.

Specificity of Immunoblotting															
Anti-serum	Purified Rat Cytochrome P450 Form (50 ng)														Cat. No.
	1A1	1A2	1B1	2A1	2A2	2B1	2B2	2C6	2C11	2C13	2D1	2E1	3A2	4A1	
Anti-CYP1B1*	-	-	+++	-	-	-	-	-	-	-	-	-	-	-	458511

*Specificity was determined by Western blotting a cDNA-expressed rat cytochrome P450 panel.

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