

## Antibody Specificity Tables

### Human UGT Immunoblotting Antibodies

The following is a summary of the immunoblotting crossreactivities of antibodies which have applications to the study of human UDP-glucuronosyl transferase (UGT) enzymes in human tissues. All of these antibodies are polyclonal.

Specificity of Immunoblotting								
Anti-sera	Human UDP-Glucuronosyl Transferase							Cat. No.
	1A1	1A4	1A6	1A9	1A10	2B7	2B15	
Anti-UGT1A Subfamily	+++	+++	+++	+++	+++	-	-	458410
Anti-UGT1A1	+++	-	-	-	-	-	-	458411
Anti-UGT1A6	-	-	+++	-	-	-	-	458416
Anti-UGT2B7	-	-	-	-	-	+++	-	458427

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