

Antibody Cross Reactivity Table

B= bovine, C= chicken, D= dog, H= human,
M= mouse, P= porcine, R= rat, Rb= rabbit NA= not applicable

| Antibody, Catalog #, Clone, | Quantity | Immunogen | Immunized host animal | Cross reactivity | No cross reactivity |
|--------------------------------|----------|-----------------------|-----------------------|------------------|---------------------|
| Bromodeoxyuridine | | | | | |
| TAK M050 BU6-4 | 100 µg | BSA-Bromodeoxyuridine | M | | |
| E-cadherin | | | | | |
| TAK M106 HECD-1 | 100 µg | H | M | | B,C,D,R |
| TAK M126 SHE78-7 | 100 µg | H | M | | R |
| TAK M107 ECCD-1 | 100 µg | M | R | | B,C,D,H |
| TAK M108 ECCD-2 | 100 µg | M | R | H, D | B,C |
| P-cadherin | | | | | |
| TAK M109 PCD-1 | 100 µg | M | R | | C,H |
| TAK M127 NCC-CAD-299 | 100 µg | H | M | | |
| N-cadherin | | | | | |
| TAK M110 NCD-2 | 100 µg | C | R | | H,M |
| Polyclonal TAK M142 | 400 µg | H (peptide) | Rb | B,C,M | |
| P, E, N-cadherin | | | | | |
| Polyclonal TAK M143 | 400 µg | H (peptide) | Rb | C,M | |
| Candida albicans | | | | | |
| Polyclonal TAK M170 | 10 µg | C. albicans | Rb | | |
| α-Catenin | | | | | |
| Polyclonal TAK M133 | 400 µg | H (peptide) | Rb | P, M | |
| Dentin Matrix Protein 1 | | | | | |
| Polyclonal TAK M176 | 100 µg | R (peptide) | Rb | H, M | |
| Enolase, C. albicans | | | | | |
| Polyclonal TAK M169 | 200 µg | C. albicans | Rb | | |

B= bovine, C= chicken, D= dog, H= human,
M= mouse, P= porcine, R= rat, Rb= rabbit NA= not applicable

| Antibody, Catalog #, Clone, | Quantity | Immunogen | Immunized host animal | Cross reactivity | No cross reactivity |
|--------------------------------|----------|--|-----------------------|---|-----------------------------------|
| Fibronectin | | | | | |
| Tak M001 FN8-12 | 400 µg | H | M | | B, P, R, Rb |
| TAK M002 FN12-8 | 400 µg | H | M | B | P, R, Rb |
| TAK M009 FN9-1 | 400 µg | H | M | B, P, Rb | R |
| TAK M010 FN30-8 | 400 µg | H | M | | B, P, R, Rb |
| TAK M013 FN1-1 | 400 µg | H | M | B, P, Rb | R |
| TAK M014 FN21-1 | 400 µg | H | M | P, Rb | B, R |
| TAK M114 FNC4-4 | 400 µg | H | M | B, P, Rb | R |
| TAK M115 FNH3-8 | 400 µg | H | M | Rb | B, P, R |
| GMP-140 | | | | | |
| TAK M062 WGA-1 | 100 µg | H (platelets) | M | | D, P, R, Rb |
| TAK M063 PL7-6 | 100 µg | H (platelets) | M | | D, P, R, Rb |
| GP Ib | | | | | |
| TAK M064 WGA-3 | 100 µg | H (platelets) | M | | |
| TAK M123 GUR20-5 | 100 µg | H (platelets) | M | | |
| GP IIb/IIIa Complex | | | | | |
| TAK M122 PT25-2 | 100 µg | H (platelets) | M | | |
| Heme Oxygenase-1 | | | | | |
| TAK M174 GST-1 | 100 µg | R | M | H, M | Rb |
| TAK M175 GST-3 | 100 µg | R | M | H | Rb, M |
| Influenza | | | | | |
| TAK M145 C179 | 100 µg | Influenza A virus A/Osaka/57 strain (H2N2) | M | H2N3 of chicken influenza | HA in H2N2 subtype |
| TAK M146 F49 | 100 µg | Influenza A virus A/India/055/1957 strain (H3N2) | M | A/India/055/1957 (H3N2) and A/Moscow/10/59 (H3N2) | HA in H1N1 and H2N2 subtypes |
| TAK M147 C111 | 100 µg | Influenza A virus A/Osaka/57 strain (H2N2) | M | | B virus |
| TAK M148 9D6 | 100 µg | Influenza B virus B/Hagaoka/1/57 strain | M | | all subtypes of Influenza A virus |
| Polyclonal TAK M149 | 400 µg | Vaccine strains of Influenza A/Hongkong/229E(H1N1), A/Wuhan/359/05(H3N2), B/Moscow/10/59 and B/Guangdong/05/99 | Rb | Influenza A & B virus | |
| Insulin | | | | | |
| TAK M056 IS11-1 | 200 µg | H | M | B, P | |
| Laminin | | | | | |
| TAK M020 LN82-13 | 100 µg | H | M | Rb | M |
| Mal f2 | | | | | |
| TAK M172 5B4 | 200 µg | M. furfur | BALB/c mouse | | |
| MnSOD, Candida albicans | | | | | |
| Polyclonal TAK M168 | 200 µg | C. albicans | Rb | | |

B= bovine, C= chicken, D= dog, H= human,
M= mouse, P= porcine, R= rat, Rb= rabbit NA= not applicable

| Antibody, Catalog #, Clone, | Quantity | Immunogen | Immunized host animal | Cross reactivity | No cross reactivity |
|-------------------------------|----------|-----------------|-----------------------|-------------------|---------------------|
| Osteocalcin | | | | | |
| TAK M041 OC4-30 | 100 µg | B | M | C, D, H, P, Rb, R | M |
| TAK M042 OCG2 | 100 µg | B | M | H | M, R |
| TAK M043 OCG3 | 100 µg | B | M | H, P, R, Rb | M |
| TAK M044 OCG4 | 100 µg | B | M | C, D, H, P, Rb | M, R |
| TAK M171 GluOC4-5 | 100 µg | H | M | B, D, P, R, Rb | |
| Polyclonal TAK M173 | 100 µg | M (peptide) | Rb | | R |
| Osteonectin | | | | | |
| TAK M124 OSN4-2 | 100 µg | H | M | B, P, Rb | |
| TAK M125 ON1-1 | 100 µg | B | M | H, P, Rb | |
| Phosphotyrosine | | | | | |
| TAK M138 PY20 | 100 µg | Phosphotyrosine | M | H, C, D, M, R | |
| TAK M141 | 100 µg | | | | |
| PIP | | | | | |
| TAK M011 PC5-5 | 100 µg | H | M | B, D | R, Rb |
| TAK M012 PC8-7 | 100 µg | H | M | B, D | R, Rb |
| SAP2, Candida albicans | | | | | |
| TAK M166 9H3C7 | 200 µg | C. albicans | M | | |
| Polyclonal TAK M167 | 200 µg | C. albicans | Rb | | |
| Tag DNA polymerase | | | | | |
| TAK 9002A | 250 U | NA | NA | NA | NA |
| TAK 9002B | 1000 U | | | | |
| Vitronectin | | | | | |
| TAK M017 VN58-1 | 200 µg | H | M | | B |
| von Willebrand Factor | | | | | |
| TAK M025 VW28-1 | 200 µg | H | M | | B |
| TAK M027 VW40-1 | 200 µg | H | M | | B |
| TAK M028 VW53-2 | 200 µg | H | M | | B |
| TAK M029 VW92-3 | 200 µg | H | M | | B |
| TAK M116 VW1-2 | 200 µg | H | M | | |

Note: For research use only. Not for use in diagnostic or therapeutic procedures.