

# Anti-Cadherin Antibody Specification Table

Clone	Cat.#	Reactivity	No Reactivity	Application	Immunogen	Immunized Host	Subclass	Ab titer
<b>E-cadherin</b>								
HECD-1	M106	🇮🇹, GP	🐄, 🐶, 🐰, 🐷	🔍, 🌨, +, ⊖	🇮🇹 breast tumor cell line MCF-7	BALB/c 🐭	🐭 IgG <sub>1</sub>	1:10,000
SHE78-7	M126	🇮🇹	🐭	🔍, 🌨, +, ⊖	Soluble 🇮🇹 placenta E-cadherin	BALB/c 🐭	🐭 IgG <sub>2a</sub>	1:10,000
ECCD-1	M107	🐭	🇮🇹, 🐄, 🐷, 🐶	⊖	🐭 teratocarcinoma cell line F9	Wister 🐭	🐭 IgG <sub>2b</sub>	
ECCD-2	M108	🐭, 🇮🇹, 🐶	🐄, 🐷	🌨, +	🐭 teratocarcinoma and mouse liver E-cadherin fragment	Donryu 🐭	🐭 IgG <sub>2a</sub>	
<b>P-cadherin</b>								
NCC-CAD-299	M127	🇮🇹		🌨, +, ⊖	🇮🇹 epidermal carcinoma cell line A-431	BALB/c 🐭	🐭 IgG <sub>1</sub>	1:4,000
PCD-1	M109	🐭	🇮🇹, 🐷	🌨, +, ⊖	🐭 endoderm cell line PSA5-E	Donryu 🐭	🐭 IgG <sub>2a</sub>	
<b>N-cadherin</b>								
NCD-2	M110	🐷, 🐭	🇮🇹, 🐭	🌨, +, ⊖	🐷 embryonic neural retina	Wister 🐭	🐭 IgG <sub>2a</sub>	
Ncad 1-1-3	M180						IgG <sub>1</sub>	
Polyclonal	M142	🇮🇹, 🐄, 🐭, 🐷	🇮🇹 E, P-cadherin	🔍, 🌨, +	peptide of 🇮🇹 N-cadherin(808-827)-KLH conjugate	🐭		
<b>P,E,N-cadherin</b>								
Polyclonal	M143	🇮🇹 P, E, N cad. 🐷 N cad. 🐭 E cad.		🔍, 🌨, +	Inter cellular C-terminal peptide of human N-cadherin-KLH conjugate	🐭		

🇮🇹 = human	🐷 = chicken	🐴 = horse	🐭 = xenopus	🔍 = Western blotting analysis under reducing or non-reducing conditions	🐭 = Immunoprecipitation
🐄 = bovine	🐭 = mouse	🐰 = sheep	🔍 = Histology on paraffin embedded tissue	📉 = flow cytometry	🌨 = Western blotting analysis under non-reducing and non-heating condition
🐶 = dog	🐭 = rabbit	GP = guinea pig	🌨 = Histology on frozen tissue sections	⊖ = Inhibition assay dependent cell-cell contact	
🐭 = rat	🐷 = porcine	G = goat			