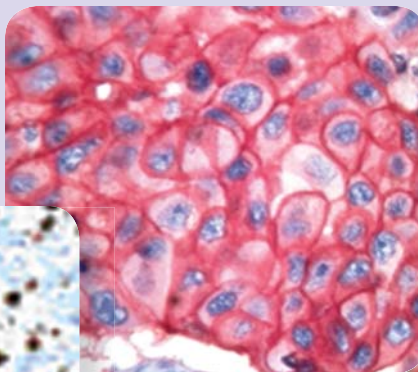
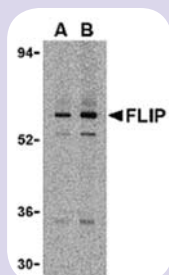


Monoclonal & Polyclonal Antibodies

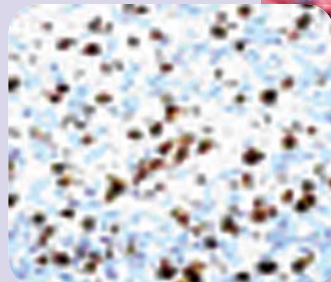
Economy research-grade size antibodies with the affordable prices
 All antibodies are tested on ELISA and Western Blot
 Applications: Flow Cytometry, IHC, Western Blot and ELISA
 Innovative products of exceptional quality



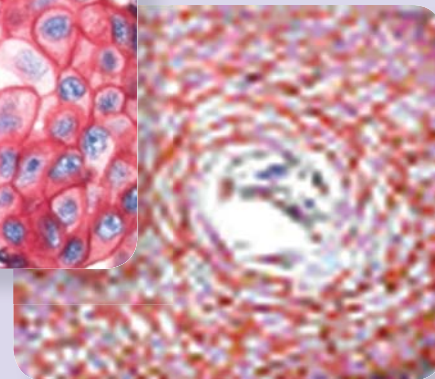
Left: Human breast cancer stained with anti-ErbB2 (Her-2) polyclonal antibody. Below: Human breast cancer stained with anti-Cadherin.



Western blot analysis of FLIP in K562 cell lysate with FLIP antibody at (A) 1 and (B) 2 µg/ml.



Left: Human tonsil stained with anti-Ki-67.



Alzheimers' Disease Related Antibodies

Antibody	Poly/Monoclonal	Size	Catalog #
Anti-β-Amyloid (1-40)	P	50 µg	54295
Anti-APP (NT), Anti-APP (CT)	P	50 µg	54095, 54096
Anti-APP [pT668]	P	50 µg	54296
Anti-BACE-1, Anti-BACE-2	P	50 µg	54101, 54100
Anti-caspase-4; Anti-caspase-7	P	50 µg	54206, 54211
Anti-Presenilin-1; Anti-Presenilin-2	P	50 µg	54297, 54298

Alzheimers' Disease β-Amyloid Protein Detection Set	Poly/Monoclonal	Size	Catalog #
Anti-APP (NT), Anti-APP(CT); Anti-BACE, Anti-BACE-2	P	N/A	54294

Neuroscience Antibodies

Product	Poly/Monoclonal	Size	Catalog #
Anti-Myelin Basic Protein (1-11)	P	50 µg	54006
Anti-Neurofilament	M	50 µg	53705
Anti-Neuron Specific Endolase (NSE)	P	50 µg	53724
Anti-AKT (pan)	P	50 µg	54299
Anti-Akt [pS473]	P	50 µg	28101
Anti-Akt/PKC[pT308]	P	50 µg	28103
Anti-PLP (139-151) and Anti-PLP (178-191)	P	50 µg	54004, 54005
Anti-MOG (35-55)	P	50 µg	54003

Adipokines, Adiponectin, and CTRPs 1-7 Detection Set	Poly/Monoclonal	Size	Catalog #
Adiponectin, CTRP1, CTRP2, CTRP3, CTRP4, CTRP5, CTRP6, and CTRP7	P	N/A	54293

Cancer Research Antibodies

Product	Poly/Monoclonal	Size	Catalog #
Anti ErbB-2(Her-2) (pY1222) NEW t Anti-ErbB - 2NEW	P	50 µg	29613, 29614
Anti-Actin, smooth muscle	P	50 µg	29553, 29553
Anti-MMP-1, Anti-MMP-2	P	50 µg	53667, 53674
Anti-MMP-3, Anti-MMP-7. Ask for more	P	50 µg	53676, 53677
Anti-Melanoma (gp100)	M	50 µg	53639
Anti-p27kip1	P	50 µg	53744
Anti-EGFR	M	50 µg	53449
Anti-Estrogen Receptor	P	50 µg	53467
Anti-P53 (pSer15)	P	50 µg	28010, 28010
Anti-Cadherin (Pan)	P	50 µg	53279
Anti-MAGE (IN)	P	50 µg	28004, 28004
Anti-Ki-67	P	50 µg	53611
Anti-MDM-2	M	50 µg	53631

Signal Transduction Antibodies

Product	Poly/Monoclonal	Size	Catalog #
Anti-SOCS-1 (CT) Anti-SOCS-3 (CT)	P	50 µg	28132; 28134
Anti-IRS-2 (pSer731)	P	50 µg	28122
Anti-Sox9 (paired), Anti-Sox9 (p Ser181)	P	50 µg	28139, 28138
Anti-CREB	P	50 µg	53410
Anti-NFκBp65 (pSer276)	P	50 µg	29775
Anti-NOS (Paired116)	P	50 µg	28130
Anti-Rho-Kinase (pThr396)	P	50 µg	29648
Anti-Glucagon	P	50 µg	29667
Anti-IRAK	P	50 µg	53592
Anti-CXCR4 (NT); Anti-CXCR4 (EL)	P	50 µg	28148, 54007
Anti-CCR5 (NT)	P	50 µg	28143

Apoptosis Related Antibodies

Product	Poly/Monoclonal	Size	Catalog #
Anti-BAD	P	50 µg	29753
Anti-CAD/DFF40 (I17)	P	50 µg	28035
Anti-BIM (IN)	P	50 µg	54013
Anti-Trail	P	50 µg	54008
Anti-FLIP (NT)	P	50 µg	54011
Anti-DR5	P	50 µg	28065
Anti-ADAM10 (CT)	P	50 µg	54012
Anti-PUMA (CT), Anti-PUMA (NT)	P	50 µg	54019, 54020
Anti-Survivin (NT)	P	50 µg	54016

Tag Antibodies

Product	Poly/Monoclonal	Size	Catalog #
Anti-GFP Tag	P	50 µg	29779
Anti-HisTag	P	50 µg	29673

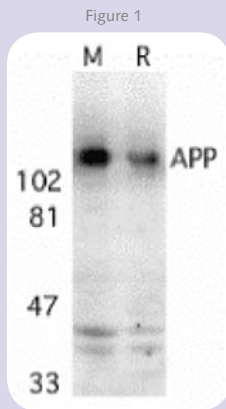


Figure 1

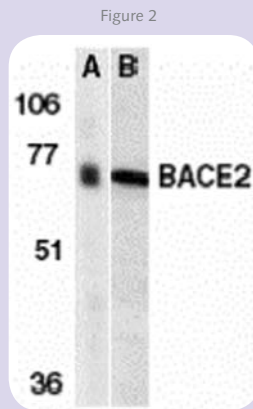


Figure 2

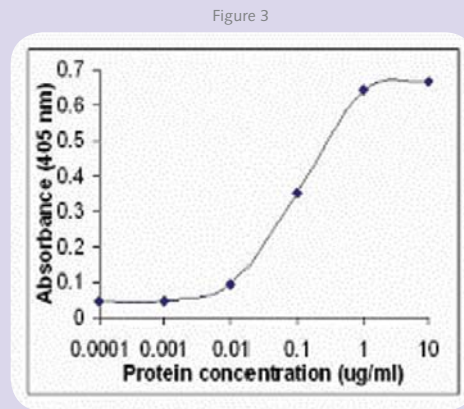


Figure 3

Figure 1. Western blot analysis of APP in mouse (M) and rat (R) brain tissue lysates with APP antibody at 2 µg/ml. Figure 2. Western blot analysis of BACE2 in human (A) and mouse (B) heart tissue lysates with BACE2 antibody at 1 µg/ml. Figure 3. AKT2 IgY Antibody

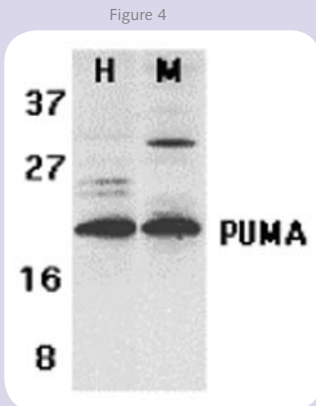


Figure 4

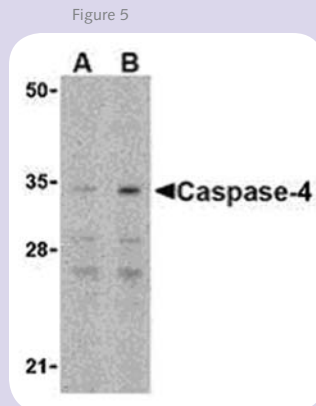


Figure 5

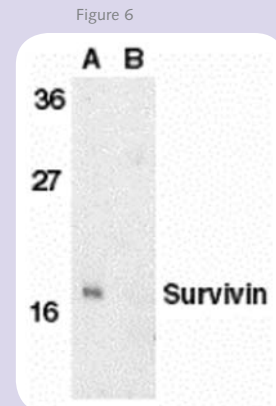


Figure 6

Figure 4. Western blot analysis of PUMA expression in human (H) K562 and mouse (M) 3T3 cell lysates with PUMA antibody at 2g/ml. Figure 5. Western blot analysis of caspase-4 in Ramos cells with caspase-4 antibody at (A) 0.5 and (B) 1 µg/ml. Figure 6. Western blot analysis of survivin in MOLT4 cell lysate in the absence (A) or presence (B) of blocking peptide with survivin antibody at 1 µg/ml.