



1	TAATACGACT	CACTATAGGG	GAATTGTGAG	CGGATAACAA	TTCCCTCTA	GAATAATTT	TGTTAACTT	TAAGAAGGAG	ATATACATAT	GGGGCGAAA	TCTAACGGTG
111	AGAAGAAGTA	CATGTGTGGT	TTCAAGCAAG	GCTTTAAGAG	CTGTGCGAAG	AAAGAGGATG	TGATCTCCGA	GAAAGGTGGT	AAGCTGCAGA	AATGCTTCAA	ATACGTCGAC
221	GCGGCGAGCG	CAACCCTGAA	TGAGAAAGCA	GTGGAAGAAC	TGAAGAAAGA	TCCGAGCGTT	GCCTATGTCC	AAGAGGACAA	GCTCTTCAA	GCTTTGACTA	GTACCATGGC
331	GGGATCCGG	TGCGAATTC	TCGAGGCGGC	CGCATAAGCC	CGAAAGGAAG	CTGAGTTGGC	TGCTGCCACC	GCTGACCAAT	AACTAGCATA	ACCCCTTGGG	GCCTCTAAAC
441	GGGTCTTGG	GGTATTTTC	CTGAAAGGAG	GAACATATAC	CGGATATCCC	GCAAGAGGCC	CGCGAGTACC	GGCATAACCA	AGCCTATGCC	TACAGCATCC	AGGTGTACGG
551	TGCCGAGGAT	GACGATGAGC	GCATTGTTAG	ATTTTCATACA	CGGTGCCTGA	CTGCGTTAGC	AATTTAACTG	TGATAAACTA	CCGCATTAAT	GCCTATCGAT	GATAAGCTGT
661	CAACATGAG	ATTTCTTGAA	GACGAAAGGG	CCTCGTGATA	CGCCTATTTT	TATAGGTTAA	TGTCATGATA	ATAATGGTFT	CTTAGACGTC	AGGTGGCACT	TTTCGGGGAA
771	ATGTGCGCG	AACCCCTAT	TGTTTATTTT	TCTAAATACA	TTCAAATATG	TATCCGCTCA	TGAGACAATA	ACCCGTATAA	ATGCTTCAAT	AATATTGAAA	AAGGAAGAGT
881	ATGAGTATTC	AACATTTCCG	TGTCGCCCTT	ATTCCTTTT	TTGCGGCATT	TTGCCTTCT	GTTTTTGCTC	ACCCGAGAAC	GCTGGTGAAA	GTAAAAGATG	CTGAAGATCA
991	GTTGGGTGCA	CGATGGGTT	ACATCGAAT	GGATCTCAAC	AGCGGTGAAG	TCCTTGAGAG	TTTTTCGCC	GAAGAACGTT	TTCCAATGAT	GAGCATTTTT	AAAGTTCTGC
1101	TATGTGGCG	GGTATTATC	CGTGTGACG	CCGGGCAAGA	GCAACTCGGT	CGCCGCATAC	ACTATTCTCA	GAATGACTTG	GTGTGACT	CACCACTGAT	AGAAAAGCAT
1211	CTTACGGATG	GCATGACAGT	AAGAGAATTA	TGCAGTGCTG	CCATAACCAT	GAGTGATAAC	ACTGCGGCCA	ACTTACTTCT	GACAACGATC	GGAGGACCGA	AGGAGCTAAC
1321	CGCTTTTTTG	CACAACATGG	GGGATCATGT	AACTCGCCCT	GATCGTTGGG	AACCGGAGCT	GAATGAAGCC	ATACCAAACG	ACGAGCGTGA	CACCACGATG	CCTCGACCAA
1431	TGGCAACAAC	GTTCGCGAAA	CTATTAACAG	GCGAAGTACT	TACTCTAGCT	TCCGCGCAAC	AATTAATAGA	CTGGATGGAG	GCGGATAAAG	TTGAGGACC	ACTTCTCGGC
1541	TCGGCCCTTC	CGGTGGGCTG	GTTTATTGCT	GATAAATCTG	GAGCCGGTGA	GCGTGGGTCT	CGCGGTATCA	TTGCGAGCACT	GGGGCCAGAT	GGTAAGCCCT	CCGCTATCGT
1651	AGTTATCTAC	ACGACGGGGA	GTCAGGCAAC	TATGGATGAA	CGAAATAGAC	AGATCGCTGA	TACAGGTGCC	TCAGGTGATTA	AGCATTTGTA	ACTGTCAGAC	CAAGTTTACT
1761	CATATATACT	TTAGATTGAT	TTAAAACCTC	ATTTTTAAT	TAAAAGGATC	TAGGTGAAGA	GCATTTTTGA	TAATCTCATG	ACCAAATCC	CTTAACGCTGA	GTTTTGCTTC
1871	CACTGAGCGT	CAGACCCCGT	AGAAAAGATC	AAAGGATCTT	CTTGAGATCC	TTTTTTTCTG	CGCGTAATCT	GCTGCTTGCA	AACAAAAAAA	CCACCGCTAC	CAGCGTGTGT
1981	TTGTTTGGCG	GATCAAGAGC	TACCAACTCT	TTTTCCGAG	GTAAGTGGCT	TCAGCAGAGC	GCAGATACCA	AATACTGTCC	TTCTAGTGA	GCCGTAGTTA	GGCCACCACT
2091	TCAAGAACTC	TGTAGCACCG	CTACATATAC	TCGCTCTGCT	AATCCTGTTA	CCAGTGCGTG	CTGCCAGTGG	CGATAAGTCG	TGCTTACCG	GGTTGGACTC	AAGACGATAG
2201	TTACCCGGATA	AGGCGCAGCG	GTCGGGCTGA	ACGGGGGGTT	CGTGACACCA	GCCCAGCTTG	GAGCGAACGA	CCTACACCGA	ACTGAGATA	CTACAGCGTG	AGCTATGAGA
2301	AAGCGCCACG	CTTCCCGAAG	GGAGAAAGGC	GGACAGGTAT	CCGTAAGCC	GCAGCGAGG	AACAGGAGAG	CGCAGGAGG	AGCTTCCAGC	AGGAAACGCC	TGGTATCTTT
2421	ATAGTCTGTG	GGGTTTFCG	CACCTCTGAC	TTGAGCGCTG	ATTTTTGTGA	TACCTCTGCG	GGGGCGGAG	CCATATGAAA	AGCCGACGCA	ACCGCCCTT	TTTACGGTTT
2531	CTGGCCTTTT	GCTGGCCTTT	TGCTCACATG	TTCTTCTCTG	CGTTATCCCC	TGATCTGTG	GATAACCGTA	TTACCCGCTT	TGAGTGAGCT	GATACCGCTC	CGCCGAGCCG
2641	AACGACCGAG	CGCAGCGAGT	CAGTGAGCGA	GGAAGCGGAA	GAGGGCCTGA	TGCGGTATTT	TCTCCTTACG	CATCTGTGCG	GTATTTTACA	CCGCATATAT	GGTGCACTCT
2751	CAGTACAATC	TGCTCTGATG	CGCATAGT	AAGCCAGTAT	ACACTCCGCT	ATCGCTACGT	GACTGGGTCA	TGGCTGCGCC	CCGACACCCG	CCAACACCCG	CTGACGCGCC
2861	CTGACGGGCT	TGCTGCTCC	CGGCATCCGC	TTACAGACAA	GCTGTGACCG	TCTCCGGGAG	CTGCATGTGT	CAGAGGTTTT	CACCGTCACT	ACCGAAACGC	GCGAGGCGAG
2970	TGCGGTAAAG	CTCATCAGCG	TGGTCTGAA	CGGATTCACA	GATGCTGCGC	TGTTTATCCG	CGTCCAGCTC	GTTGAGTTTC	TCCAGAAGCG	TTAATGTCTG	GCTTCTGATA
3081	AAGCGGGCCA	GGTTAAGGGC	GGTTTTTTCG	TGTTTGGTGA	CTGATGCGTC	CTGTGAAGGG	GCTATTTTGT	TCATGGGGGT	AATGATCCG	AGTAAACGAG	GAGAGGATGCT
3191	CACGATACGG	GTTACTGATG	ATGAACATGC	CCGGTACTG	GAACGTTGTG	AGGGTAAACA	ACTGGCGGTA	TGGATGCGCG	GGGACCAGAG	AAAAATCACT	CAGGGTCAAT
3301	GCCAGCGCTT	CGTTAATACA	GATGTAGGTG	TTCCACAGGG	TAGCCAGCAG	CATCTGCGA	TGCAGATCCG	GAACATAATG	GTGACGGGCG	CTGACTCCG	CGTTTCCAGA
3411	CTTTACGAAA	CACGAAAC	GAAGACCATT	CATGTTGTTG	CTCAGGTCGC	AGACGTTTTG	CAGCAGCAGT	CGCTTCAAGT	TCGCTCGCGT	ATCGGTGATT	CATTTCTGCTA
3521	ACCAGTAAGG	CAACCCCGCC	AGCCTAGCCG	GGTCTCAAC	GACAGGAGCA	CGATCATGCG	CACCCGTGGC	CAGGACCCAA	CGTGCCCGCA	GATGCGCCCG	GTGCGGCTGC
3631	TGGAGATGGC	GGAGCGGCG	GATATGTTCT	GCCAAGGGTT	GGTTTGGCGA	TTACAGTTTC	TCCGCAAGAA	TGATTTGGCT	CCAATCTTGG	GAGTGGTGA	TCCCTTAGCG
3741	AGGTGCGCC	GGCTTCCATT	CAGTTCGAGG	TGCGCCGCGT	CCATGACCCG	CGCAGCAACG	GACGAGGACA	GACAAAGGTAT	AGCGCCGCGC	AGGACATCCA	TGCCAACCCG
3851	TTCCATGTGC	TCGCGGAGCG	GGCATAAATC	GCCGTGACGA	TCAGCGGTCC	AGTGATCGAA	GTTAGGCTGG	TAAGAGCCCG	GAGCGATCCT	TGAAGCTGTC	CCTGATGGTC
2961	GTCATCTACC	TGCTGGGACA	GCATGGCCTG	CAACGCGGGC	ATCCGAGTGC	CGCCGGAAGC	GAGAAGAATC	ATAATGGGGA	AGGCCATCCA	GCCTCGGTC	GCGAACGCCA
4071	GCAAGACGTA	GCCAGCGCG	TCGGCCGCCA	TGCGCGGAT	AATGGCTCGC	TCTCGCCGA	AACGTTTGGT	GGCGGGACCA	GTGACGAAGG	CTTGAGCGAG	GGCGTCAAG
4181	ATTCGGAATA	CCGCAAGCGA	CAGGCCGATC	ATCGTCCGCG	TCCAGCGAAA	GCGGTCTCG	CCGAAAATGA	CCGAGAGCCG	TGCCGGCACC	TGCTCTACGA	GTTGCATGAT
4291	AAAGAAGACA	GTCATAAGTG	CGCGCAGAT	AGTCATGCC	CGCGCCCACT	GGAAGGAGCT	GACTGGGTG	AAGGCTCTCA	AGGGCATCGG	TCGAGATCCC	GGTGCCTAAT
4401	GAGTGAGCTA	ACTTACATTA	ATTGCGTTGC	GCTCACTGCC	CGTTTTCCAG	TCGGGAAACC	TGTCGTGCCA	GCTGCATTA	TGAATCGGCC	AACGCGCGGG	GAGAGGCGGT
4511	TTGCGTATTG	GGCGCCAGGG	TGTTTTTTCT	TTTTACCAGT	GAGACGGGCA	ACAGCTGATT	GCCCTTACC	GCCTGGCCCT	GAGAGAGTTG	CAGCAAGCGG	TCCACGCTGG
4621	TTTGCCCCAG	CAGGCGAAAA	TCCTGTTTGA	TGGTGGTTAA	CGCGGGGATA	TAACATGAGC	TGCTTCCGGT	ATCGTCTGAT	CCCACTACC	AGATATCCGC	ACCAACGCGC
4731	AGCCCGGACT	CGGTAATGGC	GCGCATGGCG	CCCAGCGCCA	TCTGATCGTT	GGCAACCAGC	ATCGCAGTGG	GAACGATGCC	CTCATTCAGC	ATTTGCATGG	TTTGTGAAAA
4841	ACCGGACATG	GCATCCAGT	CGCCTTCCCG	TTCCGCTATC	GGCTGAATTT	GATTCGGAGT	GATCATTTTA	TGCCAGCCAG	CCGACGCGAG	ACGCGCCGAG	ACGAACTTA
4951	ATGGGCCCCG	TAACAGCGCG	ATTTGCTGGT	GACCAATGTC	GACCAGATGC	TCCACGCCCA	GTCCGTACC	GTCTTCATGG	GAGAAAATAA	TACTGTGATG	GGGTGCTTGG
5061	TCAGAGACAT	ACAAGAAATA	CCCGGAAACA	TTAGTGCAGG	CAGCTTCCAC	AGCAATGGCA	TCCTGGTCAT	CCAGCGGATA	GTAAATGATC	AGCCCATGCA	CGCGTTGCGC
5231	GAGAAGATTG	TGCACCGCGC	CTTTACAGGC	TTGCGAGCGC	CTTCTGTTCA	CCATCGACAC	CACCACGCTG	GCACCCAGTT	GATCGGCGCG	AGATTTAATC	GCCGCGACAA
5171	TTTGGACGCG	CGGTGCGAGG	GCCAGACTGG	AGGTGGCAAC	GCCAATCAGC	AACGACTGTT	TGCCCGCCAG	TTGTTGTGCC	ACGCGGTTGG	GAATGTAATT	CAGCTCCGCC
5391	ATCGCGCTT	CCACTTTTTC	CGCGTTTTTC	GCAGAAAGCT	GGCTGGCCGT	GTTACACAGC	CGGAAACCGG	TCTGATAAGA	GACACCGGCA	TACTCTGCGA	CATCGTATAA
5501	CGTTACTGCT	TTCACTTCA	CCACCTGAA	TTGACTCTCT	TCCGGGCGCT	ATCATGCCAT	ACCCGAAAG	GTTTTGCGCC	ATTCGATGGT	GTCCGGGATC	TCGACGCTCT
5611	CCCTTATGCG	ACTCTGCGAT	TAGGAAGCAG	CCCAGTAGTA	GTTTGGAGCC	GTTGAGCACC	CGCCCGCAA	GGAAATGGTG	ATGCAAGGAG	ATGGCGCCCA	ACAGTCCCCC
5721	GGCCAGGGGG	TCTGCCACCA	TACCCACGCC	GAAACAAGCG	CTCATGAGCC	GAAAGTGGCG	AGCCCGACT	TCCCATCCG	TGATGTCCGC	GATATAGGGC	CCAGCAACCCG
5831	CACCTGTGGC	GCCGGTGATG	CCGGCCACGA	TGCGTCCGGC	GTAGAGGATC	GAGATCTCGA	TCCCGGAAA	T			

Profinity eXact vectors, tags, and resins are exclusively licensed under patent rights of Potomac Affinity Proteins. This product is intended for research purposes only. The composition and/or use of the T7 expression system is claimed in one or more patents licensed to Bio-Rad by Brookhaven Science Associates, LLC. A separate license is required for any commercial use, including use of these materials for research or production purposes by any commercial entity.



**Bio-Rad**  
Laboratories, Inc.

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600  
Canada 905 364 3435 China 86 20 8732 2339 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65  
Germany 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300 Israel 03 963 6050  
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4480 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500  
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723  
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000