

1. Vaheri, A. and Pagano, J.S. (1965) *Virology* **27**, 434.
2. Graham, F.L. and van der Eb, A.J. (1973) *Virology* **52**, 456.
3. Melton, D. *et al.* (1984) *Nucl. Acids. Res.* **12**, 7035.
4. Wong, T.K. and Neumann, E. (1982) *Biochem. Biophys. Res. Commun.* **107**, 584.
5. Fraley, R. *et al.* (1980) *J. Biol. Chem.* **255**, 10431.
6. Gorman, C.M., Moffat, L.F. and Howard, B.H. (1982) *Mol. Cell. Biol.* **2**, 1044.
7. Groskreutz, D. and Schenborn, E.T. (1997) In: *Methods in Molecular Biology* **63**, 11 ed. R. Tuan, Humana Press, NJ.
8. Southern, P.J. and Berg, P. (1982) *J. Mol. Appl. Gen.* **1**, 327.
9. McCutchan, J.H. and Pagano, J.S. (1968) *J. Natl. Cancer Inst.* **41**, 351.
10. Gluzman, Y. (1981) *Cell* **23**, 175.
11. Kawai, S. and Nishizawa, M. (1984) *Mol. Cell. Biol.* **4**, 1172.
12. Boussif, O. *et al.* (1995) *Proc. Natl. Acad. Sci. USA* **92**, 7297.
13. Haensler, J. and Szoka, F.C. (1993) *Bioconj. Chem.* **4**, 372.
14. Kukowska-Latallo, J.F. *et al.* (1996) *Proc. Natl. Acad. Sci. USA* **93**, 4897.
15. Loytner, S., Scangos, G.A. and Ruddle, F.H. (1982) *Proc. Natl. Acad. Sci. USA* **79**, 422.
16. Felgner, P.L. *et al.* (1987) *Proc. Natl. Acad. Sci. USA* **84**, 7413.
17. Capaccioli, S. *et al.* (1993) *Bioch. Biophys. Research Communications* **197**, 818.
18. Felgner, J., Bennett, F. and Felgner, P.L. (1993) *Methods* **5**, 67.
19. Lee, J.T. and Jaenisch, R. (1996) *Nucl. Acids Res.* **24**, 5054.
20. Lamb, B.T. and Gearhart, J.D. (1995) *Current Opinion in Genetics and Development* **5**, 342.
21. Malone, R.W., Felgner, P.L. and Verma, I.M. (1989) *Proc. Natl. Acad. Sci. USA* **86**, 6077.
22. Debs, R.J. *et al.* (1990) *J. Biol. Chem.* **265**, 10189.
23. Felgner, P.L. *et al.* (1995) *Ann. NY Acad. Sci.* **772**, 126.
24. Farhood, H., Servina, N.S. and Huang, L. (1995) *Biochim. Biophys. Acta* **1235**, 289.
25. Cappechi, M.R. (1980) *Cell* **22**, 479.
26. Shigekawa, K. and Dower, W.J. (1988) *BioTechniques* **6**, 742.
27. Ye, G.N., Danielle, H. and Sanford, J.C. (1990) *Plant Molec. Biol.* **15**, 809.
28. Klein, T.M. *et al.* (1987) *Nature* **327**, 70.
29. Ausubel, F. M. *et al.* (1995) *Current Protocols in Molecular Biology*. Wiley Interscience and Greene Publishing Associates.
30. Sambrook, J., Fritsch, E.F. and Maniatis, T. (1989) *Molecular Cloning: A Laboratory Manual*, Second Edition. Cold Spring Harbor Laboratory, Cold Spring Harbor, NY.
31. Zasloff, M., Ginder, G.D. and Felsenfeld, G. (1978) *Nucl. Acids Res.* **5**, 1139.
32. Sarver, N.P. *et al.* (1981) *Mol. Cell. Biol.* **1**, 486.
33. Blochlinger, K. and Diggelmann, H. (1984) *Mol. Cell. Biol.* **4**, 2929.
34. Stark, G.R., and Wahl, G.M. (1984) *Ann. Rev. Biochem.* **53**, 337.
35. Schimke, R.T. (1988) *J. Biol. Chem.* **263**, 5989.
36. Sessa, G. and Weissmann, G. (1968) *J. Lipid Res.* **9**, 310.
37. Felgner, P.L. *et al.* (1994) *J. Biol. Chem.* **269**, 2550.
38. Wheeler, C.J. *et al.* (1996) *Biochim. Biophys. Acta* **1280**, 1.
39. Kabanov, A.V. and Kabanov, V.A. (1995) *Bioconjugate Chem.* **6**, 7.
40. Moleur, F.L. *et al.* (1996) *Gene Therapy* **3**, 1010.
41. Gao, X. and Huang, L. (1995) *Gene Therapy* **2**, 710.
42. Wilson, T., Papahadjopoulos, D. and Taber, R. (1979) *Cell* **17**, 77.
43. Behr, J.-P. *et al.* (1989) *Proc. Natl. Acad. Sci. USA* **86**, 6982.
44. Loeffler, J.-P. *et al.* (1990) *J. Neurochem.* **54**, 1812.
45. Barthel, F. *et al.* (1993) *DNA and Cell Biology* **12**, 553.
46. Remy, J.S. *et al.* (1994) *Bioconj. Chem.* **5**, 647.
47. Keogh, M.-C. *et al.* (1997) *Gene Therapy* **4**, 162.
48. Demeneix, B.A. *et al.* (1994) *BioTechniques* **16**, 496.
49. Tsukamoto, M. *et al.* (1995) *Nature Gen.* **9**, 243.
50. Douar, A.-M. *et al.* (1996) *Gene Therapy* **3**, 789.
51. Gao, X., and Huang, L. (1996) *Biochemistry* **35** (3), 1027.
52. Foley, B.T., Moehring, J.M. and Moehring, T.J. (1995) *J. Biol. Chem.* **270**, 23218.
53. Harada, N. *et al.* (1997) *J. Biol. Chem.* **272**, 15232.
54. Smit, M.J. *et al.* (1996) *J. Biol. Chem.* **271**, 7574.
55. Zhang, L., David, G. and Esko, J.D. (1995) *J. Biol. Chem.* **270**, 27127.
56. Hinz, M., Moore, M.J. and Bindereif, A. (1996) *J. Biol. Chem.* **271**, 19001.
57. Nibbs, R.J.B. *et al.* (1997) *J. Biol. Chem.* **272**, 12495.
58. Venkatakrisnan, G. and Exton, J.H. (1996) *J. Biol. Chem.* **271**, 5066.
59. Fisher, E.A. *et al.* (1997) *J. Biol. Chem.* **272**, 20427.
60. Jeannin, P. *et al.* (1997) *J. Biol. Chem.* **272**, 15613.
61. Klafki, H.-W. *et al.* (1996) *J. Biol. Chem.* **271**, 28655.
62. Olofsson, A. *et al.* (1995) *J. Biol. Chem.* **270**, 31294.
63. Von Krempelhuber, A., Muller, F. and Fuhrmann, U. (1994) *J. Steroid Biochem. Mol. Biol.* **48** (5-6), 511.
64. Epstein, J.A. *et al.* (1995) *J. Biol. Chem.* **270**, 11719.
65. Choi, Y.-C. and Chae, C.-B. (1993) *Mol. Cell. Biol.* **13** (9), 5538.
66. Kizer, N., Guo, X.-L. and Hruska, K. (1997) *Proc. Natl. Acad. Sci. USA* **94**, 1013.

67. Loeffler, J.-P. and Behr, J.-P. (1993) *Meth. Enzymol.* **217**, 599.
68. Capaccioli, S. *et al.* (1993) *Biochem. Biophys. Res. Comm.* **197** (2), 818.
69. Nemoto, Y. *et al.* (1996) *J. Biol. Chem.* **271**, 13542.
70. Gaiddon, C. *et al.* (1994) *J. Biol. Chem.* **269** (36), 22663.
71. Ichijo, H. *et al.* (1997) *Science* **275**, 90.
72. Döhr, O. *et al.* (1997) *Mol. Pharmacol.* **51**, 703.
73. Lesch, K.-P. *et al.* (1996) *Science* **274**, 1527.
74. Demeneix, B.A. *et al.* (1994) *BioTech.* **16** (3), 496.
75. Rosoff, M.L., Wei, J. and Nathanson, N.M. (1996) *Proc. Natl. Acad. Sci. USA* **93**, 14889.
76. Boukhnikachvili, T. *et al.* (1997) *FEBS Lett.* **409** (2), 188.
77. Loyter, A., Scangos, G.A. and Ruddle, F.H. (1982) *Proc. Natl. Acad. Sci. USA* **79**, 422.
78. Frost, E. and Williams, J. (1978) *Virology* **91**, 39.
79. Lopata, M.A., Cleveland, D.W. and Sollner-Webb, B. (1984) *Nucl. Acids Res.* **12**, 5707.
80. Lowy, D.R., Rands, E. and Scolnick, E.M. (1978) *J. Virology* **26**, 291.
81. Lewis, W.H. *et al.* (1980) *Somat. Cell Genet.* **6**, 333.
82. Luthman, H. and Magnusson, G. (1983) *Nucl. Acids Res.* **11**, 1295.
83. Gorman, C.M., Howard, B.H. and Reeves, R. (1983) *Nucl. Acids Res.* **11**, 7631.
84. Brash, D.E. *et al.* (1987) *Mol. Cell Biol.* **7**, 2031.
85. Wilson, S.P., and Smith, A.L. (1996) *Anal. Biochem.* **246**, 148.
86. Al Molish, M.I. and Dubes, G.R. (1973) *J. Gen. Virol.* **18**, 189.
87. Freshney, R.I., (1987) *Culture of Animal Cells* A.R. Liss, Inc., NY.
88. Alam, J. and Cook, J.L. (1990) *Anal. Biochem.* **188**, 245.
89. Rosenthal, N. (1987) *Meth. Enzymol.* **152**, 704.
90. Groskreutz, D. and Schenborn, E. (1996) *Reporter Systems, In: Recombinant Proteins: Detection and Isolation.* Tuan, R., ed., Humana Press, Clifton, NJ.
91. Wood, K.V. (1995) *Curr. Opin. Biotech.* **6**, 50.
92. Hollon, T. and Yoshimura, F. K. (1989) *Anal. Biochem.* **182**, 411.
93. DeWet, J.R. *et al.* (1985) *Proc. Natl. Acad. Sci. USA* **82**, 7870.
94. DeWet, J.R. *et al.* (1987) *Cell. Biol.* **7**, 725.
95. Thompson, J.F. *et al.* (1991) *Gene* **103**, 171.
96. Pazzagli, M. *et al.* (1992) *Anal. Biochem.* **204**, 315.
97. Wood, K.V. (1991) In: *Bioluminescence and Chemiluminescence: Current Status.* Stanley, P.E. and Kricka, L.J., eds., John Wiley and Sons, NY.
98. Matthews, J.C. *et al.* (1977) *Biochemistry* **16**, 85.
99. Alton, N.K. and Vapnek, D. (1979) *Nature* **282**, 864.
100. Leslie, A.G.W. *et al.* (1988) *Proc. Natl. Acad. Sci. USA* **85**, 4133.
101. Seed, B. and Sheen, J.-Y. (1988) *Gene* **67**, 271.
102. Neumann, J.R. *et al.* (1987) *BioTechniques* **5**, 444.
103. Young, D.C. (1993) *Anal. Biochem.* **215**, 24.
104. Groskreutz, D.J. *et al.* (1995) *Promega Notes* **50**, 2.
105. Lorenz, W.W. *et al.* (1991) *Proc. Natl. Acad. Sci. USA* **88**, 4438.
106. Farr, A. and Roman, A. (1991) *Nucl. Acids Res.* **20**, 920.
107. Groskreutz, D.J. *et al.* (1996) *Promega Notes* **55**, 2.
108. Shaw, W.V. (1975) *Meth. Enzymol.* **43**, 737.
109. Cullen, B. *et al.* (1987) *Meth. Enzymol.* **152**, 687.
110. Hall, C.V. *et al.* (1983) *J. Molec. Applied Gen.* **2**, 101.
111. Silhavy, T. *et al.* (1972) In: *Experiments in Molecular Genetics*, Miller, J.H., ed., Cold Spring Harbor Laboratory, Cold Spring Harbor, NY.