

Bio-Beads® SM-2 Hydrophobic Adsorbents

Applications Bibliography

Bio-Beads SM adsorbents are neutral, macroporous polymeric beads of high surface area for adsorbing organics from aqueous solutions. Each bead is composed of a large number of highly crosslinked microspheres, giving it high surface area and uniform pores. Bio-Beads adsorbents can be used with a variety of solvents, including alcohols, petroleum ether, diethyl ether, hexane, and solvent mixtures, as well as with aqueous media. They have excellent physical stability and withstand temperatures to 250 °C.

Bio-Beads SM-2 nonpolar polystyrene adsorbents are particularly useful for the adsorption of nonpolar substances or surface active agents from aqueous solutions.

The following reference cite the use of Bio-Beads SM-2 adsorbents for the analysis of organic compounds, drugs, biologicals, foods, pesticides, chemicals, metals, and detergents. For more information on Bio-Beads SM adsorbents, request the Bio-Beads SM Instruction Manual.

Applications for Bio-Beads SM-2

Application	Reference
Organics	
Trace organics in water	Wigilius, B., Boren, H., Cvarlberg, G. E., Grimvall, A., Lundgren, B. V. and Savenhed, R., <i>J. Chromatog.</i> , 391 , 169-182 (1987).
	Junk, G.A., Richard, J. J., Grieser, M. D., Witiak, D., Witiak, J. L., Arguello, M. D., Vick, R., Svec, H. J., Fritz, J. S. and Calder, G. V., <i>J. Chromatogr.</i> , 99 , 745-762 (1974).
	Jahanir, L. M. and Samuelson, O., <i>J. Chromatogr.</i> , 193 , 197-206 (1980).
	Moore, R.A. and Karasek, F. W., <i>Inter. J. Environ. Anal. Chem.</i> , 17 , 187-202 (1984).
Osterroht, C., <i>J. Chromatogr.</i> , 101 , 289-298 (1974).	
Chlorinated hydrocarbons in water	Picer, N. and Picer, M., <i>J. Chromatogr.</i> , 193 , 357-369 (1980).
Separation of nitro- and chlorophenols	Grieser, M.D. and Pietrzyk, D. J., <i>Anal. Chem.</i> , 45 , 1348-1353 (1973).
Organic bases	Chu, C. and Pietrzyk, D. J., <i>Anal. Chem.</i> , 46 (3), 330-340 (1974).

Application	Reference
Drugs	
Retention and sorption of sulfas	Rotsch, T. D., Sydor, R. J. and Pietrzyk, D. J., <i>J. Chrom. Sci.</i> , 17 , 339-344 (1979).
Separation of cannabinoids from cannabis extract	Hendriks, H., Batterman, S., Bos, R., Huizing, H. J. and Malingre, Th. M., <i>J. Chromatogr.</i> , B205 , 444-450 (1981).
Determination of benzo[a]pyrene in cigarette smoke condensate	Robinson, J. L., Marshall, M. A., Draganjac, M. E. and Noggle, L. C., <i>Anal. Chim. Acta</i> , 115 , 229-238 (1980).
Narcotic antagonists in human urine	Digregorio, G.J., <i>J. Chromatogr.</i> , 101 , 424-427 (1974).
Drugs in urine	Roerig, D. L., Lewand, D., Mueller, M. and Wang, R. I. H., <i>J. Chromatogr.</i> , 110 , 349-359 (1975).
	Bastos, M. L., Jukofsly, D. and Mule, S. J., <i>J. Chromatogr.</i> , 81 , 93 (1973).
Morphine and codeine in urine	Ulrich, L. and Ruegsegger, P., <i>Arch. Toxicol.</i> , 45 , 241-248 (1980). German.
Morphine in urine	Kullberg, M. P. and Gorodetzky, C. W., <i>Clinical Chem.</i> , 20 (2), 177-183 (1974).
Methaqualone in blood plasma	Hux, R. A., Mohammed, H. Y. and Cantwell, F. F., <i>Anal. Chem.</i> , 54 , 113-117 (1982).

Application	Reference
Drugs in blood	Schlicht, H. J. and Gelbke, H. P., <i>Z. Rechtsmedizin</i> , 81 , 25-30 (1978).
Pharmaceutical drug syrups	Mohammed, H. Y. and Cantwell, F. F., <i>Anal. Chem.</i> , 50 (3), 491-496 (1978).
Extraction of drugs from human stomach fluid	Ibrahim G., Andryauskas, S. and Bastos, M. L., <i>J. Chromatogr.</i> , 108 , 107-116 (1975).
Polar and non-polar drug metabolites	Dieterle, W., Faigle, J. W. and Mory, H., <i>J. Chromatogr.</i> , 168 , 27-34 (1979).
Biological	
Purification of aminobenzyl-phosphonic acid	Landt, M., Boltz, S. C. and Butler, L. G., <i>Biochem.</i> , 17 (5), 915-919 (1978).
Free rhodamine	Spack, E. G., Jr., Packard, B., Wier, M. L. and Edidin, M., <i>Anal. Biochem.</i> , 158 , 233-237 (1986).
Crude plasma clean-up	Tamura, M., Harris, T. M. Higashimori, K., Sweetman, B. J., Blair, I. A. and Inagami, T., <i>Biochem.</i> , 26 , 2797-2806 (1987).
Leukotrienes from plasma prostaglandins	Salari, H. and Steffenrud, S., <i>J. Chromatogr.</i> , 378 , 35-44 (1986).
Prostaglandins	Hamberg, M., <i>Anal. Biochem.</i> , 55 , 368-378 (1973).
Prostaglandins adsorption from biological fluid	Leffler, C. W., Desiderio, D. M. and Wakelyn, C. E., <i>Prostaglandins</i> , 21 , 227-241 (1981).
Isolation of proline and proline from fossil bone	Stafford, T. W., Duhamel, R. C., Haynes, C. V. and Brendel, K., <i>Life Sciences</i> , 31 , 931-938 (1982).
Anabolic steroids in animal tissue and urine	Verbeke, R., <i>J. Chromatogr.</i> , 177 , 69-84 (1979).
Steroids from urine	Shackleton, C. H. L., Sjovall, J. and Wisen, O., <i>Clin. Chim. Acta</i> , 27 , 354-356 (1970).
Extraction of steroids	Bradlow, H. L., <i>Steroids</i> , 11 , 265 (1968).
Bile acids	Okuyama, S., Uemura, D. and Hirata, Y., <i>Bulletin of Chemical Society of Japan</i> , 52 (1), 124-126 (1979).
Bile acids	Schwartz, H. P., v. Bergmann, K. and Paumgartner, G., <i>Clinica Chimica Acta</i> , 50 , 197-206 (1974).

Application	Reference
Cobalamins	Fenton, W. A. and Rosenberg, L. E., <i>Anal. Biochem.</i> , 90 , 119-125 (1978).
Phosphatidyl-ethanolamine from plasma	Salari, H., <i>J. Chromatogr.</i> , 419 , 103-111 (1987).
Glucuronides	Delaborde, S., Loosli, H. R., Ponelle, M., Griesser, R. and Maurer, G., <i>J. High Resolution Chromatography & Chromatography Communications</i> , 10 , 71-76 (1987).
Impregnation of resin with pentafluorobenzyl bromide microbial fatty acid analysis	Rosenfeld, J. M., Hammerberg, O. and Orvidas, M. C., <i>J. Chromatogr.</i> , 378 , 9-16 (1986).
Plant hormones	Stafford, A. E., Kunhle, J. A., Corse, J. and Hautala, E., <i>J. Chromatogr.</i> , 294 , 485-488 (1984).
Captopril (derivatized in blood)	Funke, P.T., Ivashkiv, E., Malley, M.F. and Cohen, A.I., <i>Anal. Chem.</i> , 52 , 1086-1089, (1980).
Food	
Flavonoids from aqueous fractions	Rosler, K. H. and Goodwin, R. S., <i>J. Natural Products</i> , 47 (1), 188 (1983).
Acid dyes	Uematsu, T., Kurita, T. and Hamada, A., <i>J. Chromatogr.</i> , 172 , 327-334 (1979).
Naringin and limonin from grapefruit juice	Chandler, B. V. and Johnson, R.I., <i>Proc. Int. Soc. Citriculture</i> , 2 , 885-888 (1981).
Triadimefon in grape wine	Nickless, G. and Spitzer, T., <i>J. Chromatogr.</i> , 208 , 409-413 (1981).
Environmental	
Indium treated adsorbent to concentrate humic complexes	Hiraide, M., Arima, Y. and Mizuike, A., <i>Anal. Chim. Acta</i> , 200 , 171-179 (1987).
Hydrocarbons, PCBs and fatty acids in sea water	Gomez-Bellnchon, J. I., Grimalt, J. O. and Albaiges, J., <i>Environ. Sci. Technol.</i> , 22 , 667-685 (1988).
Polynuclear hydrocarbons from water	Navratil, J. D., Sievers, R. E. and Walton, H. F., <i>Anal. Chem.</i> , 49 (14), 2260-2263 (1977).

Application	Reference
Pesticides	
Aminocarb insecticide from water	Levesque, D. and Mallet, V. N., <i>Intern. J. Environ. Anal. Chem.</i> , 16 , 139-147 (1983).
Carbamate insecticides	Sundaram, K. M. S., Szeto, S. Y. and Hindle, R., <i>J. Chromatogr.</i> , 177 , 29-34 (1979).
Chlorinated pesticides	McNeil, E. E. and Otsen, R., <i>J. Chromatogr.</i> , 132 , 277-286 (1977).
Ethyl and methyl parathion	Paschal, D. D., Bicknell, R. and Dresbach, D., <i>Anal. Chem.</i> , 49 , (11) 1551-1554 (1977).
Fenitrothion and its degradation products	Volpe, G. and Mallet, V. N., <i>Chromatographia</i> , 14 (6), 333-336 (1981).
Organophosphorous pesticides	LeBel, G. L., Williams, D. T., Griffith, G. and Benoit, F. M., <i>J. Assoc. Off. Anal. Chem.</i> , 62 , No. 2, 241-249 (1979).
Chemical	
Carboxylic acid from aqueous solutions	Jahangir, J. L. M. and Samuelson, O., <i>J. Chromatogr.</i> , 237 , 371-379 (1982).
	Hasanain, M. A. and Hines, A. L., <i>Ind. Eng. Chem. Process Des. Dev.</i> , 20 , 621-625 (1981).
Use in HPLC column as ground material	Baum, R. G., Saetre, R. and Cantwell, F. F., <i>Anal. Chem.</i> , 52 (1), 15-19 (1980).
Benzenesulfonic acids	Rotsch, T. D. and Pietryzk, D. J., <i>Anal. Chem.</i> , 52 (8), 1323-1327 (1980).
Use in TLC as stationary phase	Pietryzk, D. J., Ritsch, T. D. and Leuthauser, S. W., <i>J. Chromatogr. Sci.</i> , 17 , 555-561 (1979).
Metals	
Metal cation adsorption	Mackey, D. J., <i>J. Chromatogr.</i> , 242 , 275-287 (1982).
Metal ion adsorption using pyrocatechol loaded adsorbent	Brajter, K., Olbrych-Sleszynska, E. and Staskiewicz, M., <i>Talanta</i> , 35 (1), 65-67 (1988).
Resin impregnation with TBP for extracting actinides from fission products	Green, L. W., Elliot, N. L. and Longhurst, T. H., <i>Anal. Chem.</i> , 55 , 2394-2398 (1983).

Application	Reference
Detergents	
Triton X-100 detergent	Holloway, P. W., <i>Anal. Chem.</i> , 53 , 304-308 (1973).
	Drexler, G., Eichinger, A., Wolf, C. and Sieghart, W., <i>J. Immun. Methods</i> , 95 , 117-122 (1986).
	Bonomi, F. and Kurtz, D. M., Jr., <i>Anal. Biochem.</i> , 142 , 226-231 (1984).
	Welling, G. W., et. al., <i>J. Chromatogr.</i> , 297 , 101 (1984).
	Metsikko, K., van Meer, G. and Simons, K., <i>The EMBO Journal</i> , 5 (13), 3429-3435 (1986).
Cholate detergent	Bonomo, E. A. and Swaney, J. B., <i>J. Lipid Research</i> , 29 , 380-384 (1988).
	Jinks, D. C. and McElhaney, R. N., <i>Anal. Biochem.</i> , 164 , 331-335 (1987).
Deoxycholate	Horigome, T. and Sugano, H., <i>Anal. Biochem.</i> , 130 , 393-396 (1983).
	Lorusso, D. J. and Green, F. A., <i>Science</i> , 188 , 66 (1974).
	Shechter, I. and Bloch, K., <i>J. Biol. Chem.</i> , 246 (24), 7690-7696 (1971).
NP-40 detergent	Momoi, T., <i>Biochem. Biophys. Res. Commun.</i> , 87 , 541-549 (1979).
Emulgen 911	Gibson, G. G., and Schenkman, F. B., <i>J. Biol. Chem.</i> , 253 , 5957-5963 (1978).
	Warner, M., <i>J. Biol. Chem.</i> , 257 (21), 12995-13000 (1982).
Emulphogene BC-720	Brunch, R. C., Thotakura, N. R. and Bahl, O. P., <i>J. Biol. Chem.</i> , 261 (20), 9450-9460 (1986).
Coating microtiter plates with detergent soluble membrane proteins	Drexler, G., Eichinger, A., Wolf, C. and Sieghart, W., <i>J. Immunological Methods</i> , 95 , 117-122 (1986).



**Bio-Rad
Laboratories**

**Life Science
Group**

*Bio-Rad Laboratories Main Office, 2000 Alfred Nobel Drive, Hercules, California 94547, Ph. (510) 741-1000, Fx. (510) 741-5800
Also in: North Ryde, Australia, Ph. 02-805-5000, Fx. 02-805-1920 Wien, Austria, Ph. (1) 877 89 01, Fx. (1) 876 56 29 Nazareth, Belgium, Ph. 09-385 55 11, Fx. 09-385 65 54
Mississauga, Canada, Ph. (905) 712-2771, Fx. (905) 712-2990 Beijing, China, Ph. (01) 2046622, Fx. (01) 2051876 Copenhagen, Denmark, Ph. 39 17 9947, Fx. 39 27 1698
Espoo, Finland, Ph. 90 804 2200, Fx. 90 804 1100 Ivry sur Seine Cedex, France, Ph. (1) 49 60 68 34, Fx. (1) 46 71 24 67 München, Germany, Ph. 089 31884-0, Fx. 089 31884-100
New Delhi, India, Ph. 91-11-461-0103, Fx. 91-11-461-0765 Milano, Italy, Ph. 02-21609.1, Fx. 02-21609.399 Tokyo, Japan, Ph. 03-5811-6270 Fx. 03-5811-6272 Veenendaal,
The Netherlands, Ph. 0318-540666, Fx. 0318-542216 Auckland, New Zealand, Ph. 09-443 3099, Fx. 09-443 3097 Kowloon, Hong Kong, Ph. 7893300, Fx. 7891257 Singapore,
Ph. (65) 272-9877, Fx. (65) 273-4835 Solna, Sweden, Ph. 46 (0) 8 735 83 00, Fx 46 (0) 735 54 60 Madrid, Spain, Ph. (91) 661 70 85, Fx. (91) 661 96 98 Glattbrugg, Switzerland,
Ph. 01/809 55 55, Fx. 01/809 55 00 Hemel Hempstead, United Kingdom, Ph. 0800 181134, Fx. 01442 259118*