



Bio-Rad Protein Assay Applications Bibliography

General Discussion

The following references contain discussions about the Bradford (Bio-Rad) Protein Assay. These include analyses of the dye-binding method, and adaptations of the assay methodology which increase its versatility.

Bradford, M., "A Rapid and Sensitive Method for the Quantitation of Microgram Quantities of Protein Utilizing the Principle of Protein-Dye Binding," *Anal. Biochem.*, **72**, 248 (1976).

Bearden, J., "Quantitation of Sub-Microgram Quantities of Protein by an Improved Protein-Dye Binding Assay," *Biochim. Biophys. Acta*, **533**, 525 (1978).

Brogdon, W. and Dickinson, C., "A Microassay System for Measuring Esterase Activity and Protein Concentration in Small Samples and in High-Pressure Liquid Chromatography Eluate Fractions," *Anal. Biochem.*, **131**, 499 (1983).

Compton, S. J. and Jones, C. G., "Mechanism of Dye Response and Interference in the Bradford Protein Assay," *Anal. Biochem.*, **151**, 369 (1985).

Duhamel, R., Meezan, E. and Brendel, K., "The Addition of SDS to the Bradford Dye-Binding Protein Assay, A Modification with Increased Sensitivity to Collagen," *J. Biochem. Biophys. Methods*, **5**, 67 (1981).

Fanger, B., "Adaptation of the Bradford Protein Assay to Membrane-Bound Proteins by Solubilizing in Glucopyranoside Detergents," *Anal. Biochem.*, **162**, 11 (1987).

Gogstad, G. and Krutnes, M., "Measurement of Protein in Cell Suspensions Using the Coomassie Brilliant Blue Dye-Binding Assay," *Anal. Biochem.*, **126**, 355 (1982).

Macart, M. and Gerbaut, L., "An Improvement of the Coomassie Blue Dye Binding Method Allowing an Equal Sensitivity to Various Proteins: Application to Cerebrospinal Fluid," *Clinica Chimica Acta*, **122**, 93 (1982).

Marshall, T. and Williams, K., "Phenol Addition to the Bradford Dye Binding Assay Improves Sensitivity and Gives a Characteristic Response with Different Proteins," *J. Biochem. Biophys. Methods*, **13**, 145 (1986).

McIntosh, J., "Application of a Dye-Binding Method to the Determination of Protein in Urine and CSF," *Clin. Chem.*, **23**, 1939 (1977).

Nesbitt, J., Sauer, D. and Culpepper, R., "Use of the Bradford Dye-Binding Method for CSF and Urine Total Protein Determination," *J. Amer. Med. Technol.*, **40**, 278 (1978).

Pierce, J. and Suelter, C., "An Evaluation of the Coomassie Brilliant Blue G-250 Dye Binding Method for Quantitative Protein Determination," *Anal. Biochem.*, **81**, 478 (1977).

Pilgrim, L., "Protein Measurement," *U. Georgia Research Reporter*, **9**, 6 (1977).

Pollard, H. B., Menard, R., Brandt, H. A., Pazoles, C. J., Creutz, C. E. and Ramu, A., "Application of Bradford Protein Assay to Adrenal Gland Subcellular Fractions," *Anal. Biochem.*, **86**, 761 (1978).

Ramagli, L. and Rodriguez, L., "Quantitation of Microgram Amounts of Protein in Two-Dimensional Polyacrylamide Gel Electrophoresis Sample Buffer," *Electrophoresis*, **6**, 559 (1985).

Rubin, R. and Warren, R., "Quantitation of Microgram Amounts of Protein in SDS-Mercaptoethanol-Tris Electrophoresis Sample Buffer," *Anal. Biochem.*, **83**, 773 (1977).

Simpson, I. and Sonne, O., "A Simple, Rapid, and Sensitive Method for Measuring Protein Concentration in Subcellular Membrane Fractions Prepared by Sucrose Density Ultracentrifugation," *Anal. Biochem.*, **119**, 424 (1982).

Snyder, J. and Desborou, S., "Rapid Estimation of Potato Tuber Total Protein-Content with Coomassie Brilliant Blue G-250," *Theor. Appl. Genet.*, **52**, 135 (1978).

Spector, T., "Refinement of Coomassie Blue Method of Protein Quantitation — Simple and Linear Spectrophotometric Assay for Less-Than-Or-Equal-To 0.5 to 50 Microgram of Protein," *Anal. Biochem.*, **86**, 142 (1978).

Van Kley, H. and Hale, S., "Assay for Protein by Dye Binding," *Anal. Biochem.*, **81**, 485 (1977).

Specific Applications

The following references cite the dye-binding assay as a means of protein quantitation in specific applications.

Ahmad, F., Ahmad, P. M., Pieretti, L. and Watters, G. T., "Purification & Subunit Structure of Rat Mammary Gland Acetyl Co-Enzyme A Carboxylase," *J. Biol. Chem.*, **253**, 1733 (1978).

Altman, K., Pyliotis, N. and Mukkur, T., "A New Method for the Extraction and Purification of K99 Pili from Enterotoxigenic *E. coli* and their Characterization," *Biochem. J.*, **201**, 505 (1982).

- Antoniw, J. and Pierpoint, W., "Purification and Properties of One of b Proteins From Virus-Infected Tobacco Plants," *J. Gen. Virol.*, **39**, 343 (1978).
- Apps, D. and Glover, L., "Isolation and Characterization of Magnesium Adenosine Triphosphatase From Chromaffin Granule Membrane," *FEBS Letters*, **85**, 254 (1978).
- Babul, J., "Phosphofructokinases From *Escherichia coli*-Purification and Characterization on Non-Allosteric Isoenzyme," *J. Biol. Chem.*, **253**, 4350 (1978).
- Barnes, D. and Colowick S., "Stimulation of Sugar Uptake and Thymidine Incorporation in Mouse 3T3 Cells by Calcium-Phosphate and Other Extracellular Particles," *Proc. Nat. Acad. Sci.*, (USA), **74**, 5593 (1977).
- Benne, R., Edman, J., Traut, R. R. and Hershey, J. W. B., "Phosphorylation of Eukaryotic Protein Synthesis Initiation-Factors," *Proc. Nat. Acad. Sci.*, (USA), **75**, 108 (1978).
- Benne, R., Brownlue, M. L. and Hershey, J. W. B., "Purification and Characterization of Protein Synthesis Initiation Factors EIF-1, EIF-4C, EIF-40, and EIF-5 From Rabbit Reticulocytes," *J. Biol. Chem.*, **253**, 3070 (1978).
- Bennett, V., "Purification of an Active Proteolytic Fragment of Membrane Attachment Site for Human Erythrocyte Spectrin," *J. Biol. Chem.*, **253**, 2292 (1978).
- Boothby, J., Jasper, D., Lutz, H. and Rollins, M., "Gel Electrophoresis-Derived Enzyme-Linked Immunosorbent Assay of Serum from Cows Resistant to and Cows Susceptible to Challenge Exposure with *Mycoplasma bovis*," *Am. J. Vet. Res.*, **43**, 553 (1982).
- Brandon, C. and Wu, J., "Purification and Properties of Choline Acetyltransferase from *Torpedo Californica*," *J. Neurochem.*, **30**, 791 (1978).
- Brown, W., Vellom, D., Schnepf, E. and Greer, C., "Purification of a Surface-Bound Hexosaminidase From Spores of *Bacillus cereus* T.," *FEMS Microbiol. Letters*, **3**, 247 (1978).
- Burchell, B. and Hallinan, T., "Phospholipid Content and Activity of Pure Uridine Diphosphate Glucuronyl transferase from Rat Liver," *Biochem. J.*, **171**, 821 (1978).
- Creutz, C., Pazoles, C. J. and Pollard, H. B., "Identification and Purification of an Adrenal Medullary Protein (Synexin) That Causes Calcium-Dependent Aggregation of Isolated Chromaffin Granules," *J. Biol. Chem.*, **253**, 2858 (1978).
- Cronin, M. J., Roberts, J. M. and Weiner, R. I., "Dopamine and Dihydroergocryptine Binding to Anterior Pituitary and Other Brain Areas of Rat and Sheep," *Endocrinology*, **103**, 302 (1978).
- Crouch, R., Kimsey, G., Priest, D. G. and Sarda, A., "Effect of Streptozotocin on Erythrocyte and Retinal Superoxide Dismutase," *Diabetologia*, **15**, 53 (1978).
- Decker, D. and Levinson, D., "Quantitation of Rat Liver Xanthine Oxidase by Radioimmunoassay," *Arthritis and Rheumatism*, **25**, 326 (1982).
- Dehlinger, P. and Litt, M., "Ornithine Decarboxylase Induction and Nucleolar RNA Synthesis in Friend Leukemia-Cells," *Biochem. Biophys. Res. Comm.*, **81**, 1054 (1978).
- Dolhofer, R. and Wieland, O., "In vitro Glycosylation of Hemoglobins by Different Sugars and Sugar Phosphates," *FEBS Letters*, **85**, 86 (1978).
- Duhamel, R., "Differential Staining of Collagens and Non-Collagens with Coomassie Brilliant Blue G and R," *Collagen Rel. Res.*, **3**, 195 (1983).
- Dunn, S., "Identification of Altered Subunit in Inactive F1-ATPase of an *Escherichia coli*-Unca Mutant," *Biochem. Biophys. Res. Comm.*, **82**, 596 (1978).
- Farley, M., O'Brien, T., Moyer, D. and Taylor, C., "The Detection of Estrogen Receptors in Gynecologic Tumors Using Immunoperoxidase and the Dextran-Coated Charcoal Assay," *Cancer*, **49**, 2153 (1982).
- Feinstein, S. and Altman, S., "Context Effects on Nonsense Codon Suppression in *Escherichia coli*," *Genetics*, **88**, 201 (1978).
- Forrester, L. J., Barrett, J. T. and Campbell, B. J., "Red Blood Cell Lysis Induced by Venom of Brown Recluse Spider - Role of Sphingomyelinase-D," *Arch. Biochem. Biophys.*, **187**, 355 (1978).
- Fuller, D. J. M., Byus, C. V. and Russell, D. H., "Specific Regulation by Steroid Hormones of Amount of Type-1 Cyclic AMP-Dependent Protein-Kinase Holoenzyme," *Proc. Nat. Acad. Sci.*, (USA), **75**, 223 (1978).
- Ganu, V. and Alworth, W., "Epoxide Hydrase Structure-Activity-Relationships of *In vitro* Stimulators of Microsomal Activity," *Biochemistry*, **17**, 2876 (1978).
- Gast, R. and Lee, J., "Isolation of the *In vivo* Emitter in Bacterial Bioluminescence," *Proc. Nat. Acad. Sci.*, (USA), **75**, 833 (1978).
- Goldstein, L.D., Reynolds, C. P. and Perezpol, J. R., "Isolation of Human Nerve Growth Factor From Placental Tissue," *Neurochem. R.*, **3**, 175 (1978).
- Gros, C., Giros, B. and Schwartz, J., "Identification of Aminopeptidase M as an Enkephalin-Inactivating Enzyme in Rat Cerebral Membranes," *Biochemistry*, **24**, 2179 (1985).
- Halling, S. M., Burtis, K. C. and Doi, R. H., "Reconstitution Studies Show That Rifampicin Resistance is Determined by Largest Polypeptide of *Bacillus subtilis* RNA-Polymerase," *J. Biol. Chem.*, **252**, 9024 (1977).
- Hamaguchi, H., *et al.*, "Genetic Analysis of Human Lymphocyte Proteins by Two-Dimensional Electrophoresis," *Human Genetics*, **60**, 176 (1982).
- Hardie, D. and Cohen, P., "Regulation of Fatty-Acid Biosynthesis—Simple Procedure for Purification of Acetyl Co-A Carboxylase from Lactating Rabbit Mammary Gland, and its Phosphorylation by Endogenous Cyclic AMP-Dependent and AMP-Independent Protein Kinase Activities," *FEBS Letters*, **91**, 1 (1978).
- Hersh, L. B., Coe, B. and Casey, L., "Fluorometric Assay for Choline Acetyltransferase and its Use in Purification of Enzyme from Human Placenta," *J. Neurochem.*, **30**, 1077 (1978).
- Hupe, D. and Behrens, N., "A Method for Assaying Orotate Phosphoribosyltransferase and Measuring Phosphoribosylpyrophosphate," *Anal. Biochem.*, **161**, 20 (1987).

- Ikegami, M. and Fraenkel-Conrat, H., "RNA-Dependent RNA-Polymerase of Tobacco Plants," *Proc. Nat. Acad. Sci.*, (USA), **75**, 2122 (1978).
- Iqbal, M. J., Ford, P. and Johnson, M. W., "Purification of Sex Hormone-Binding Globulin Using Affinity Matrix in an Electrophoretic Cell System," *FEBS Letters*, **87**, 235 (1978).
- Jablonski, E. and DeLuca, M., "Purification and Properties of the NADH and NADPH-Specific FMN Oxidoreductases from *Beneckea harveyi*," *Biochemistry*, **16**, 2932 (1977).
- Jacquot, J., Benali, R., Zahm, J. and Puchelle, E., "Identification of Different Molecular Forms of Human Airway Lysozyme," *Anal. Biochem.*, **160**, 227 (1987).
- Jarrett, H. and Penniston, J., "Purification of Ca^{2+} -Stimulated ATPase Activator from Human Erythrocytes — its Membership in Class of Ca^{2+} -Binding Modulator Proteins," *J. Biol. Chem.*, **253**, 4676 (1978).
- Johnson, W. and Kramer, T., "Effect of Copper Deficiency on Erythrocyte Membrane Protein of Rats," *J. Nutrition*, **117**, 1085 (1987).
- Kasahara, M. and Penefsky, H., "High Affinity Binding of Mono Valent P(I) by Beef Heart Mitochondrial Adenosine Triphosphatase," *J. Biol. Chem.*, **253**, 4180 (1978).
- Kobiler, D., Beyer, E. C. and Barondes, S. H., "Developmentally Regulated Lectins from Chick, Muscle, Brain and Liver Have Similar Chemical and Immunological Properties," *Develop. Bio.*, **64**, 265 (1978).
- Kolakofsky, D. and Altman, S., "Endo-Ribonuclease Activity Associated with Animal RNA Viruses," *J. Virology*, **25**, 274 (1978).
- Kramer, A., Frick, M. and Keller, W., "Separation of Multiple Components of HeLa Cell Nuclear Extracts Required for Pre-messenger RNA Splicing," *J. Biol. Chem.*, **262**, 17630 (1978).
- Kuhn, B., *et al.*, "Genetic Control of Chalcone-Flavanone Isomerase Activity in *Callistephus chinensis*," *Planta*, **138**, 199 (1978).
- Kuret, J. and Schulman, H., "Purification and Characterization of a Ca^{2+} /Calmodulin-Dependent Protein Kinase from Rat Brain," *Biochemistry*, **23**, 5495 (1984).
- La Porte, D. and Storm, D., "Detection of Calcium-Dependent Regulatory Protein-Binding Components Using I^{125} -Labeled Calcium-Dependent Regulatory Protein," *J. Biol. Chem.*, **253**, 3374 (1978).
- Liu, L. and Wang, J., "*Micrococcus luteus* Gyrase-Active Components and a Model for Its Supercoiling of DNA," *Proc. Nat. Acad. Sci.*, (USA), **75**, 2098 (1978).
- Mahan, K., Hoffmann, P. and Hoffmann, C., "A Reproducible Method for the Optimum Infection of Human Mononuclear Cells and Lymphoid Cell Lines By Herpes Simplex Virus Type I," *J. Virological Methods*, **4**, 63 (1982).
- McCain, D., *et al.*, "Assay for Early Cytoplasmic Effects of the SRC Gene Product of Rous Sarcoma Virus," *Proc. Nat. Acad. Sci.*, (USA), **75**, 2750 (1978).
- McLemore, T. L., Martin, R. R., Pickard, L. R., Springer, R. R., Wray, N. P., Toppell, K. L., Mattox, K. L., Guinn, G. A., Cantrell, E. T. and Busbee, D. L., "Analysis of Arylhydrocarbon Hydroxylase Activity in Human Lung Tissue, Pulmonary Macrophages, and Blood Lymphocytes," *Cancer*, **41**, 2292 (1978).
- Millay, R., *et al.*, "The Reduced NADH-Activated Phosphoenol Pyruvate Carboxylase from *Pseudomonas MA*," *J. Biol. Chem.*, **253**, 1371 (1978).
- Ogunyemi, E., *et al.*, "Studies on Biosynthesis of Cyclitols .36. Purification of Myo-Inositol-1-Phosphate Synthase of Duckweed, *Lemna gibba*, to Homogeneity by Affinity Chromatography on NAD-Sepharose Molecular and Catalytic Properties of Enzymes," *H-S Z. Physiol. Chem.*, **359**, 613 (1978).
- Panasenko, S.M., Alazard, R. J. and Lehman, I. R., "Simple, 3-Step Procedure for Large Scale Purification of DNA-Ligase from a Hybrid Lambda Lysogen Constructed *In vitro*," *J. Biol. Chem.*, **253**, 4590 (1978).
- Pettigrew, D. W. and Frieden, C., "Rabbit Muscle Phosphofructokinase — Modification of Molecular and Regulatory Kinetic Properties with Affinity Label 5' -*p*-(Fluorosulfonyl) Benzoyl Adenosine," *J. Biol. Chem.*, **253**, 3623 (1978).
- Phillips, J. and Allison, Y., "Proton Translocation by the Bovine Chromaffin Granule Membrane," *Biochem. J.*, **170**, 661 (1978).
- Powell, J. and Morrison, J., "Purification and Properties of Aspartate-Aminotransferase and Aromatic-Amino-Acid Aminotransferase from *Escherichia coli*," *Eur. J. Biochem.*, **87**, 391 (1978).
- Reczek, P. and Vilee, C., "A Purification of Microsomal Glucose-6-Phosphatase from Human Tissue," *Biochem. Biophys. Res. Com.*, **107**, 1158 (1982).
- Riquelme, P., Fox, R. W. and Kemp, R. G., "Mouse Muscle Phosphofructokinase is Partially Phosphorylated," *Biochem. Biophys. Res. Comm.*, **81**, 864 (1978).
- Rosen, B. P., Brey, R. N. and Hasan, S. M., "Energy Transduction in *Escherichia coli* — New Mutation Affecting FO Portion of ATP-Synthetase Complex," *J. Bact.*, **134**, 1030 (1978).
- Roskoski, R., Ngan, P., Mettenburg, R. and Lund, D. D., "Isoproterenol Injection Alters Protein Kinase DEAE Cellulose Profiles in Selected Rat Tissues," *Biochem. Biophys. Res. Comm.*, **82**, 641 (1978).
- Rubin, R. W., Warren, R. H., Lukeman, D. S. and Clements, E., "Actin Content and Organization in Normal and Transformed Cells in Culture," *J. Cell. Bio.*, **78**, 28 (1978).
- Scherr, R. P. and Thauer, R., "Purification and Properties of Reduced Ferredoxin CO_2 Oxidoreductase from *Clostridium pasteurianum*, a Molybdenum Iron-Sulfur-Protein," *Eur. J. Biochem.*, **85**, 125 (1978).
- Sharma, R. K., Desai, R., Thompson, T. R. and Wang, J. H., "Purification of Heat-Stable Inhibitor Protein of Ca^{2+} -Activated Cyclic Nucleotide Phosphodiesterase by Affinity Chromatography," *Can. J. Biochem.*, **56**, 598 (1978).

Sharma, H. and Tihon, C., "Determination of Avidin and Streptavidin by a Modified Bradford Method," *Anal. Biochem.*, **170**, 135 (1988).

Shaw, J., Monticone, V. and Paetkau, V., "Partial Purification and Molecular Characterization of a Lymphokine (Costimulator) Required for Mitogenic Response of Mouse Thymocytes *In vitro*," *J. Immunol.*, **120**, 1967 (1978).

Sherman, M. R., Pickering, L. A., Rollwagen, F. M. and Miller, L. K., "Mero-Receptors: Proteolytic Fragments of Receptors Containing the Steroid Binding Site," *Fed. Proc.*, **37**, 167 (1978).

Sica, V., et al., "Assay of Total Estradiol Receptor in Tissue Homogenate and Tissue Fractions by Exchange with Sodium Thiocyanate at Low Temperature," *Biochemistry*, **20**, 686 (1981).

Spooner, J., Daley, L., Ware, G. and Vines, M., "Determination of Irrigation and Fertilizer Practices for Jade Plant *Crassula argentea*," *J. Am. S. Hort.*, **103**, 306 (1978).

Stripf, R. and Werner, D., "Differentiation of *Rhizobium japonicum* .2. Enzymatic-Activities in Bacteroids and Plant Cytoplasm During Development of Nodules of Glycine-Max," *Z. Naturf. C.*, **33**, 373 (1978).

Svasti, J. and Bowman, B., "Human Group-Specific Component Changes in Electrophoretic Mobility Resulting from Vitamin D Binding and From Neuraminidase Digestion," *J. Biol. Chem.*, **253**, 4188 (1978).

Thompson, H. J., Sharma, S. K. and Brown, S. A., "O-Methyltransferases of Furanocoumarin Biosynthesis," *Arch. Biochem. Biophys.*, **188**, 272 (1978).

Weiel, J., Hershey, J. W. B. and Levison, S. A., "Fluorescence Polarization Studies of Binding of Fluorescein-Labeled Initiation Factor IF3 to 30S-Ribosomal Subunits from *Escherichia coli*," *FEBS Letters*, **87**, 103 (1978).

Weisz, A., et al., "Interaction of Rat Liver Glucocorticoid Receptor with Heparin," *Biochemistry*, **23**, 5393 (1984).

Werner, D. and Stripf, R., "Differentiation of *Rhizobium japonicum* .1. Enzymatic Comparison of Nitrogenase Repressed and Derepressed Free Living Cells and of Bacteroids," *Z. Naturf. C.*, **33**, 245 (1978).

Wiedner, H., et al., "Nonessential Nature of Sulfhydryl Groups for ATPase Activity in Myosin-Cyanylation Study," *J. Biol. Chem.*, **253**, 2763 (1978).

Yagi, K., Yazawa, M., Kakiuchi, S., Ohshima, M. and Uenishi, K., "Identification of an Activator Protein for Myosin Light Chain Kinase as the Ca-dependent Modulator Protein," *J. Biol. Chem.*, **253**, 1338 (1978).

Ordering Information

Catalog Number	Product Description
500-0001	Bio-Rad Protein Assay Kit I: 450 ml Dye Reagent Concentrate and lyophilized bovine gamma globulin protein standard. Sufficient for 450 protein assay determinations with standards procedure or 2,250 assays with microassay.
500-0002	Bio-Rad Protein Assay Kit II: Same as 500-0001, except protein standard is bovine serum albumin
500-0006	Bio-Rad Protein Assay Dye Reagent Concentrate , 450 ml
500-0005	Protein Standard I: Bovine gamma globulin, lyophilized. Will provide 20 ml, 1.4 mg/ml
500-0007	Protein Standard II: Bovine serum albumin, lyophilized. Will provide 20 ml, 1.4 mg/ml

BIO-RAD

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: Regents Park, Australia, Ph. 02-9914-2800, Fx. 02-9914-2889 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. (86-10) 2046622, Fx. (86-10) 2051876 K benhavn  , 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 M nchen, 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 Veendam,
The Netherlands, Ph. 031318-540666, Fx. 0313 18-542216 Auckland, New Zealand, Ph. 09-443 3099, Fx. 09-443 3097 Kowloon, Hong Kong, Ph. 7893300,
Ph. 7891257 Singapore, Ph. (65) 272-9877, Fx. (65) 273-4835 Sundbyberg, Sweden, Ph. 46 (0) 8 627 50 00, Fx. 46 (0) 8 627 54 00 Madrid, Spain, Ph. (91) 661 70 85,
Ph. (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