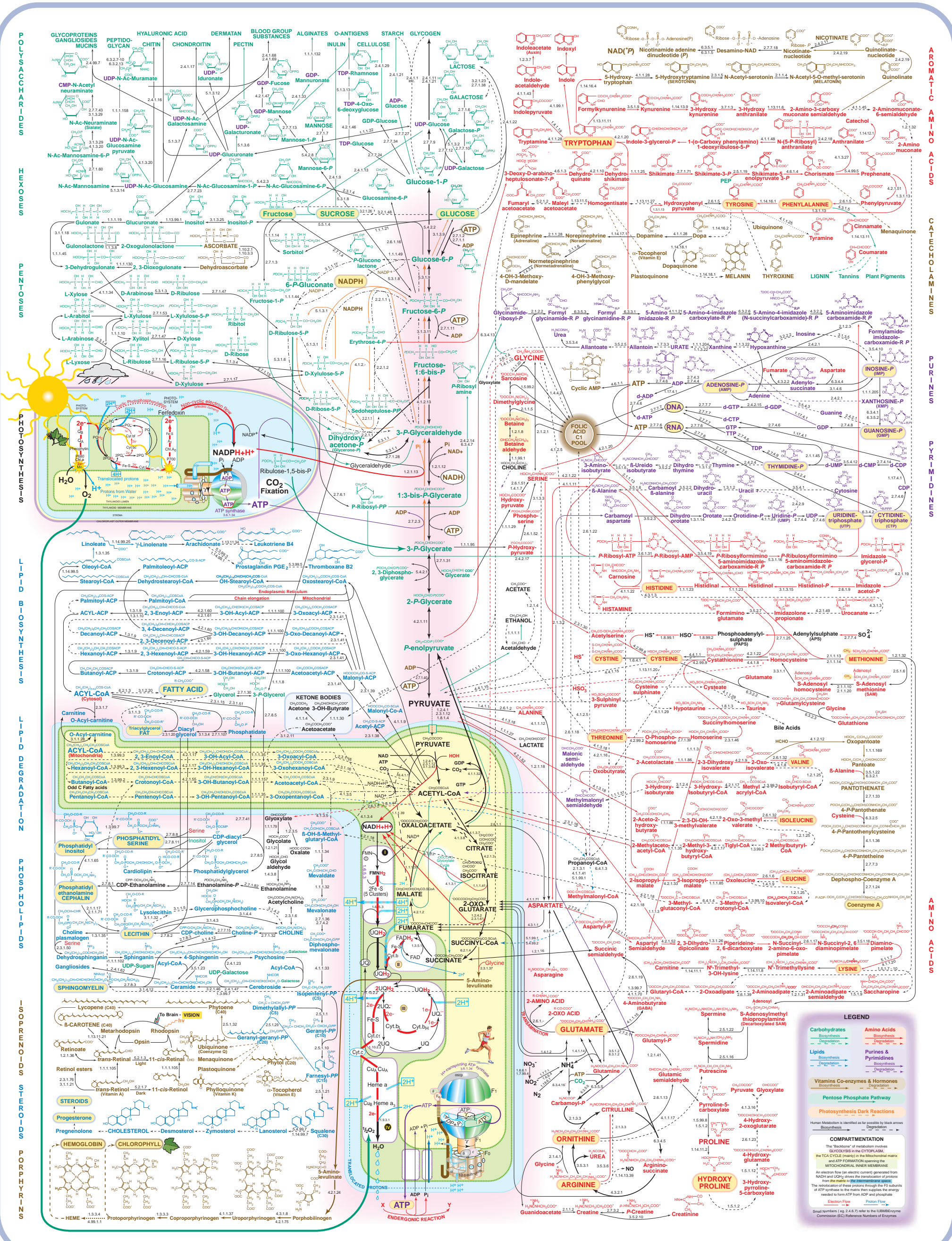


# Metabolic Pathways



© 2003 International Union of Biochemistry and Molecular Biology www.iubmb.org

22<sup>nd</sup> Edition Designed by Donald E. Nicholson, D.Sc., The University of Leeds, England – and Sigma-Aldrich

Cat. No. M3782



<b>Argentina</b> SIGMA-ALDRICH DE ARGENTINA S.A. Free Tel: 0811 888 7446 Tel: (+54) 11 4556 1472 Fax: (+54) 11 4552 1698	<b>Belgium</b> SIGMA-ALDRICH N.V.S.A. Free Tel: 0800 14747 Free Fax: 0800 14745 Tel: (+32) 3 899 13 01 Fax: (+32) 3 899 13 11	<b>Chile</b> SIGMA-ALDRICH QUIMICA LIMITADA Free Tel: 800 174 725 Tel: (+56) 2 495 7395 Fax: (+56) 2 495 7396	<b>Denmark</b> SIGMA-ALDRICH DENMARK A/S Tel: (+45) 43 56 59 00 Fax: (+45) 43 56 59 05	<b>Germany</b> SIGMA-ALDRICH CHEMIE GMBH Free Tel: 0800 51 55 000 Free Fax: 0800 64 90 000 Tel: (+49) 89 6513 10 Fax: (+49) 89 6513 1160	<b>India</b> SIGMA-ALDRICH CHEMICALS PRIVATE LIMITED Bangalore: (+91) 80 6621 9400 New Delhi: (+91) 11 4358 8000 Mumbai: (+91) 22 2570 2364 Hyderabad: (+91) 40 4015 5488 Kolkata: (+91) 33 4013 8003 Bangalore: (+91) 80 6621 9650 New Delhi: (+91) 11 4358 8001 Mumbai: (+91) 22 2579 7589 Hyderabad: (+91) 40 4015 5466 Kolkata: (+91) 33 4013 8016	<b>Ireland</b> SIGMA-ALDRICH IRELAND LTD. Free Tel: 1800 200 888 Free Fax: 1800 602 222 Tel: (+353) 402 20370 Fax: (+353) 402 20375	<b>Japan</b> SIGMA-ALDRICH JAPAN K.K. Tel: (+81) 3 5796 7300 Tel: (+81) 3 5796 7315	<b>Korea</b> SIGMA-ALDRICH KOREA Free Tel: (+82) 80 023 7111 Free Fax: (+82) 80 023 8111 Tel: (+82) 31 329 9000 Fax: (+82) 31 329 9000	<b>Israel</b> SIGMA-ALDRICH ISRAEL LTD. Free Tel: 1 800 70 2222 Tel: (+972) 8 948 4100 Fax: (+972) 8 948 4200	<b>Italy</b> SIGMA-ALDRICH S.r.l. Numero Verde: 800 827018 Tel: (+39) 02 3341 7310 Fax: (+39) 02 3801 0737	<b>Mexico</b> SIGMA-ALDRICH QUIMICA, S.A. de C.V. Free Tel: 01 800 007 5300 Free Fax: 01 800 712 9920 Tel: (+52) 722 276 1600 Fax: (+52) 722 276 1601	<b>The Netherlands</b> SIGMA-ALDRICH CHEMIE BV Free Tel: 0800 022 9089 Free Fax: 0800 022 9088 Tel: (+31) 78 620 5411 Fax: (+31) 78 620 5421	<b>New Zealand</b> SIGMA-ALDRICH NEW ZEALAND LTD. Free Tel: 0800 936 666 Free Fax: 0800 937 777 Tel: (+61) 9841 0555 Fax: (+61) 9841 0500	<b>Norway</b> SIGMA-ALDRICH NORWAY AS Tel: (+47) 23 17 60 00 Fax: (+47) 23 17 60 10	<b>Palau</b> SIGMA-ALDRICH SPC. z.o.o. Tel: (+48) 61 829 01 20 Fax: (+48) 61 829 01 20	<b>Portugal</b> SIGMA-ALDRICH QUIMICA, S.A. Free Tel: 800 202 178 Free Fax: 800 202 178 Tel: (+351) 21 924 3555 Fax: (+351) 21 924 2610	<b>Russia</b> SIGMA-ALDRICH RUS, LLC Tel: (+7) 495 621 5979 Fax: (+7) 495 621 5923	<b>Singapore</b> SIGMA-ALDRICH PTE. LTD. Tel: (+65) 6779 1200 Free Fax: (+65) 6779 1822 Tel: (+34) 91 661 96 97 Fax: (+34) 91 661 96 42	<b>Spain</b> SIGMA-ALDRICH QUIMICA, S.A. Free Tel: 900 101 376 Free Fax: 900 102 028 Tel: (+34) 91 661 96 97 Fax: (+34) 91 661 96 42	<b>Sweden</b> SIGMA-ALDRICH SWEDEN AB Tel: (+46) 8 742 4200 Fax: (+46) 8 742 4243	<b>South Africa</b> SIGMA-ALDRICH (PTY) LTD. Free Tel: 0800 100 800 Free Fax: 0800 100 800 Tel: (+27) 11 979 1188 Fax: (+27) 11 979 1189	<b>Switzerland</b> SIGMA-ALDRICH CHEMIE GMBH Free Tel: 0800 80 00 80 Free Fax: 0800 80 00 80 Tel: (+41) 81 755 2828 Fax: (+41) 81 755 2815	<b>United Kingdom</b> SIGMA-ALDRICH COMPANY LTD. Free Tel: 0800 771 181 Free Fax: 0800 378 785 Tel: (+44) 1747 833 000 Fax: (+44) 1747 833 313	<b>United States</b> SIGMA-ALDRICH Toll-Free: 800 325 5052 Tel: (+1) 314 771 5765 Fax: (+1) 314 771 5765	<b>Vietnam</b> SIGMA-ALDRICH PTE. LTD. VN R.O. Tel: (+84) 316 2810 Fax: (+84) 6258 4238
--	--	---	---	---	---	--	--	---	---	--	--	---	--	--	---	--	---	--	---	--	---	---	---	--	--