



Malvern Panalytical

THE PERIODIC TABLE OF THE ELEMENTS

1	1 1.008 H Hydrogen																	2 4.003 He Helium	1						
2	3 6.939 Li Lithium	4 9.012 Be Beryllium																	5 10.811 B Boron	6 12.011 C Carbon	7 14.007 N Nitrogen	8 15.999 O Oxygen	9 18.998 F Fluorine	10 20.179 Ne Neon	2
3	11 22.990 Na Sodium	12 24.312 Mg Magnesium																	13 26.982 Al Aluminium	14 28.086 Si Silicon	15 30.974 P Phosphorus	16 32.064 S Sulfur	17 35.453 Cl Chlorine	18 39.948 Ar Argon	3
4	19 39.102 K Potassium	20 40.080 Ca Calcium	21 44.956 Sc Scandium	22 47.900 Ti Titanium	23 50.942 V Vanadium	24 51.996 Cr Chromium	25 54.938 Mn Manganese	26 55.847 Fe Iron	27 58.933 Co Cobalt	28 58.693 Ni Nickel	29 63.546 Cu Copper	30 65.380 Zn Zinc	31 69.720 Ga Gallium	32 72.640 Ge Germanium	33 74.922 As Arsenic	34 78.960 Se Selenium	35 79.904 Br Bromine	36 83.800 Kr Krypton	4						
5	37 85.470 Rb Rubidium	38 87.620 Sr Strontium	39 88.905 Y Yttrium	40 91.220 Zr Zirconium	41 92.906 Nb Niobium	42 95.960 Mo Molybdenum	43 98.000 Tc Technetium	44 101.070 Ru Ruthenium	45 102.905 Rh Rhodium	46 106.400 Pd Palladium	47 107.868 Ag Silver	48 112.400 Cd Cadmium	49 114.820 In Indium	50 118.690 Sn Tin	51 121.75 Sb Antimony	52 127.600 Te Tellurium	53 126.904 I Iodine	54 131.293 Xe Xenon	5						
6	55 132.905 Cs Caesium	56 137.340 Ba Barium	L 57 - 71 Lanthanides	72 178.490 Hf Hafnium	73 108.948 Ta Tantalum	74 183.850 W Tungsten	75 186.200 Re Rhenium	76 190.200 Os Osmium	77 192.200 Ir Iridium	78 195.090 Pt Platinum	79 196.967 Au Gold	80 200.590 Hg Mercury	81 204.383 Tl Thallium	82 207.190 Pb Lead	83 208.980 Bi Bismuth	84 (209) Po Polonium	85 (210) At Astatine	86 (222) Rn Radon	6						
7	87 (223) Fr Francium	88 (226) Ra Radium	A 89 - 103 Actinides	104 (267) Rf Rutherfordium	105 (270) Db Dubnium	106 (269) Sg Seaborgium	107 (270) Bh Bohrium	108 (270) Hs Hassium	109 (278) Mt Meitnerium	110 (281) Ds Darmstadtium	111 (281) Rg Roentgenium	112 (285) Cn Copernicium	113 (286) Nh Nihonium	114 (289) Fl Flerovium	115 (289) Mc Moscovium	116 (293) Lv Livermorium	117 (293) Ts Tennessee	118 (294) Og Oganesson	7						
			L	57 138.910 La Lanthanum	58 140.120 Ce Cerium	59 140.907 Pr Praseodymium	60 144.240 Nd Neodymium	61 (145) Pm Promethium	62 150.360 Sm Samarium	63 151.960 Eu Europium	64 157.250 Gd Gadolinium	65 158.92 Tb Terbium	66 162.500 Dy Dysprosium	67 164.930 Ho Holmium	68 167.260 Er Erbium	69 168.934 Tm Thulium	70 173.040 Yb Ytterbium	71 174.967 Lu Lutetium							
			A	89 (227) Ac Actinium	90 232.038 Th Thorium	91 231.000 Pa Protactinium	92 238.030 U Uranium	93 (237) Np Neptunium	94 (244) Pu Plutonium	95 (243) Am Americium	96 (247) Cm Curium	97 (247) Bk Berkelium	98 (251) Cf Californium	99 (252) Es Einsteinium	100 (257) Fm Fermium	101 (258) Md Mendelevium	102 (259) No Nobelium	103 (262) Lr Lawrencium							
	Metals	Metal-like	Lanthanides																	Non-metals	Noble gasses	Actinides			

Atomic number 29 63.546 Atomic weight
Element symbol **Cu** Element name
Copper