

periodic table of the elements

Tabla periódica de los elementos

1 1 H 1.0079 -259.1 -252.76 2.20 -1, 1 1s ¹	2 4 Be 9.0122 1287 2471 1.47											18 2 He 4.0026 -272.2 -268.93 1s ²					
3 2 Li 6.941 180.5 1342 0.97	4 2 Be 9.0122 1287 2471 1.47											10 2 Ne 20.1797 -248.609 -246.053 [He] 2s ² 2p ⁶					
11 3 Na 22.990 97.80 883 1.01	12 2 Mg 24.305 650 1090 1.23											18 2 Ar 39.948 -185.36 -185.85 [Ne] 3s ² 3p ⁶					
19 4 K 39.098 63.5 759 0.91	20 2 Ca 40.078 842 1484 1.04	21 3 Sc 44.956 1541 2836 1.20	22 4 Ti 47.867 1668 3407 1.32	23 5 V 50.942 1910 3407 1.45	24 6 Cr 51.996 1907 2671 1.56	25 7 Mn 54.938 1246 2061 1.60	26 8 Fe 55.845 1538 2861 1.64	27 9 Co 58.933 1495 2927 1.70	28 10 Ni 58.693 1455 2913 1.75	29 11 Cu 63.546 1084.62 2562 1.75	30 12 Zn 65.409 419.53 907 1.66	31 3 Ga 69.723 29.76 2204 1.82	32 4 Ge 72.64 938.25 2833 2.02	33 5 As 74.922 subl. 616 2204 2.20	34 6 Se 78.96 221 685 2.48	35 7 Br 79.904 -7.2 58.8 2.74	36 2, 4 Kr 83.798 -157.36 -153.34 [Ar] 3d ¹⁰ 4s ² 4p ⁶
37 5 Rb 85.468 39.3 688 0.89	38 2 Sr 87.62 777 1377 0.99	39 3 Y 88.906 1522 3345 1.11	40 4 Zr 91.224 1855 4409 1.22	41 5 Nb 92.906 2623 4744 1.23	42 6 Mo 95.94 2477 4639 1.30	43 7 Tc *97.907 2623 4262 1.36	44 8 Ru 101.07 2334 4150 1.42	45 9 Rh 102.91 1964 3695 1.45	46 10 Pd 106.42 1554.8 2963 1.30	47 11 Ag 107.87 961.78 2162 1.42	48 12 Cd 112.41 321.07 767 1.46	49 3 In 114.82 156.6 2072 1.49	50 4 Sn 118.71 231.93 2602 1.72	51 5 Sb 121.76 630.63 1587 1.82	52 6 Te 127.60 449.51 988 2.01	53 7 I 126.90 113.7 184.4 2.21	54 2, 4, 6 Xe 131.29 -111.74 -108.09 [Kr] 4d ¹⁰ 5s ² 5p ⁶
55 6 Cs 132.91 28.44 671 0.86	56 2 Ba 137.33 727 1897 0.97	57-71 Lanthanides Lantánidos	72 4 Hf 178.49 2233 4603 1.23	73 5 Ta 180.95 3017 5455 1.33	74 6 W 183.84 3422 5555 1.40	75 7 Re 186.21 3185 5596 1.46	76 8 Os 190.23 3033 5012 1.52	77 9 Ir 192.22 2446 4428 1.55	78 10 Pt 195.08 1768.2 3825 1.42	79 11 Au 196.97 1064.18 2856 1.42	80 12 Hg 200.59 -38.83 356.62 1.44	81 1, 3 Tl 204.38 304 1473 1.44	82 2, 4 Pb 207.2 327.46 1749 1.55	83 3, 5 Bi 208.98 271.4 1564 1.67	84 2, 4, 6 Po *208.98 254 962 1.76	85 -1, 1, 3, 5, 7 At *209.99 302 336.95 1.96	86 2 Rn *222.02 -71 -61.7 [Xe] 4f ¹⁴ 5d ¹⁰ 6s ² 6p ⁶
87 7 Fr *223.02 27 677 0.86	88 2 Ra *226.03 696 1140 0.97	89-103 Actinides Actínidos	104 3 Rf *261.11	105 4, 5 Db *262.11	106 5, 6 Sg *266.12	107 6, 7 Bh *264.12	108 7, 8 Hs *277	109 8, 9 Mt *268.14	110 9, 10 Ds *281	111 10, 11 Rg *280	112 11, 12 Cn *285	113 *284 Uut [Rn] 5f ¹⁴ 6d ¹⁰ 7s ² 7p ¹	114 *287 Fl [Rn] 5f ¹⁴ 6d ¹⁰ 7s ² 7p ²	115 *288 Uup [Rn] 5f ¹⁴ 6d ¹⁰ 7s ² 7p ³	116 *291 Lv [Rn] 5f ¹⁴ 6d ¹⁰ 7s ² 7p ⁴	117 *294 Uus [Rn] 5f ¹⁴ 6d ¹⁰ 7s ² 7p ⁵	118 *294 Uuo [Rn] 5f ¹⁴ 6d ¹⁰ 7s ² 7p ⁶
		Lanthanides Lantánidos	57 3 La 138.91 920 3464 1.08	58 3, 4 Ce 140.12 799 3443 1.08	59 3, 4 Pr 140.91 931 3520 1.07	60 3 Nd 144.24 1016 3074 1.07	61 3 Pm *144.91 1042 3000 1.07	62 2, 3 Sm 150.36 1072 1794 1.07	63 2, 3 Eu 151.96 822 1596 1.01	64 3 Gd 157.25 1313 3273 1.11	65 3, 4 Tb 158.93 1359 3230 1.10	66 3 Dy 162.50 1412 2700 1.10	67 3 Ho 164.93 1472 2700 1.10	68 3 Er 167.26 1529 2868 1.11	69 2, 3 Tm 168.93 1545 1950 1.11	70 2, 3 Yb 173.04 824 1196 1.06	71 3 Lu 174.97 1663 3402 1.14
		Actinides Actínidos	89 3 Ac *227.03 1050 3198 1.00	90 3 Th *232.04 1750 4788 1.11	91 4, 5 Pa *231.04 1572	92 3, 4, 5, 6 U *238.03 1135 4131 1.22	93 3, 4, 5, 6 Np *237.05 644 3902 1.22	94 3, 4, 5, 6 Pu *244.06 640 3228 1.22	95 3, 4, 5, 6 Am *243.06 1176 2011 1.20	96 3, 4 Cm *247.07 1345	97 3, 4 Bk *247.07 996	98 3, 4 Cf *251.08 900	99 3 Es *252.08 860	100 3 Fm *257.095 1527	101 3 Md *258.10 827	102 2, 3 No *259.10	103 3 Lr *262.11

1 45
Rh
102.91
1964
3695
1.45

0, 1, 2, 3, 4, 5
[Kr] 4d⁵5s¹

- 1 Atomic number
- 2 Element symbol
- 3 Relative atomic mass
- 4 Melting point in °C
- 5 Boiling point in °C
- 6 Electronegativity (Allred, Rochow)
- 7 Oxidations states
- 8 Electron configuration

- Nonmetals, Metaloides
 - Halogens, Halógenos
 - Inert gases, Gases nobles
 - Alkali metals, Metales alcalinos
 - Transition metals, Metales de transición
 - Lanthanides, Lantánidos
 - Actinides, Actínidos
 - Other metals, Otros metales
 - Semi metals, Semi metales
- * most stable isotope, isótopo más estable



The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

MilliporeSigma and the Vibrant M are trademarks of Merck KGaA, Darmstadt, Germany. Copyright © 2017 EMD Millipore Corporation. All Rights Reserved. 2017 - 05493 06/2017