

参考資料 4

原子量累積表 (2009)

*参考資料3における原子量表を使用し、小数点3位までで表した(以下を四捨五入)

原子数	C	H	N
1	12.011	1.008	14.007
2	24.021	2.016	28.013
3	36.032	3.024	42.020
4	48.043	4.032	56.027
5	60.054	5.040	70.034
6	72.064	6.048	84.040
7	84.075	7.056	98.047
8	96.086	8.064	112.054
9	108.096	9.071	126.060
10	120.107	10.079	140.067
11	132.118	11.087	154.074
12	144.128	12.095	168.080
13	156.139	13.103	182.087
14	168.150	14.111	196.094
15	180.161	15.119	210.101
16	192.171	16.127	224.107
17	204.182	17.135	238.114
18	216.193	18.143	252.121
19	228.203	19.151	266.127
20	240.214	20.159	280.134
21	252.225	21.167	294.141
22	264.235	22.175	308.147
23	276.246	23.183	322.154
24	288.257	24.191	336.161
25	300.268	25.199	350.168
26	312.278	26.206	364.174
27	324.289	27.214	378.181
28	336.300	28.222	392.188
29	348.310	29.230	406.194
30	360.321	30.238	420.201
31	372.332	31.246	434.208
32	384.342	32.254	448.214
33	396.353	33.262	462.221
34	408.364	34.270	476.228
35	420.375	35.278	490.235
36	432.385	36.286	504.241
37	444.396	37.294	518.248
38	456.407	38.302	532.255
39	468.417	39.310	546.261
40	480.428	40.318	560.268
41	492.439	41.326	574.275
42	504.449	42.333	588.281
43	516.460	43.341	602.288
44	528.471	44.349	616.295
45	540.482	45.357	630.302
46	552.492	46.365	644.308
47	564.503	47.373	658.315
48	576.514	48.381	672.322
49	588.524	49.389	686.328
50	600.535	50.397	700.335

原子数	O	P	S
1	15.999	30.974	32.065
2	31.999	61.948	64.130
3	47.998	92.921	96.195
4	63.998	123.895	128.260
5	79.997	154.869	160.325
6	95.996	185.843	192.390
7	111.996	216.816	224.455
8	127.995	247.790	256.520
9	143.995	278.764	288.585
10	159.994	309.738	320.650
11	175.993	340.711	352.715
12	191.993	371.685	384.780
13	207.992	402.659	416.845
14	223.992	433.633	448.910
15	239.991	464.606	480.975

原子数	Br	Cl	F
1	79.904	35.453	18.998
2	159.808	70.906	37.997
3	239.712	106.359	56.995
4	319.616	141.812	75.994
5	399.520	177.265	94.992
6	479.424	212.718	113.990
7	559.328	248.171	132.989
8	639.232	283.624	151.987
9	719.136	319.077	170.986
10	799.040	354.530	189.984
11	878.944	389.983	208.982
12	958.848	425.436	227.981
13	1038.752	460.889	246.979
14	1118.656	496.342	265.978
15	1198.560	531.795	284.976

原子数	Na	K	H ₂ O
1	22.990	39.098	18.015
2	45.980	78.197	36.031
3	68.969	117.295	54.046
4	91.959	156.393	72.061
5	114.949	195.492	90.076
6	137.939	234.590	108.092
7	160.928	273.688	126.107
8	183.918	312.786	144.122
9	206.908	351.885	162.138
10	229.898	390.983	180.153
11	252.887	430.081	198.168
12	275.877	469.180	216.183
13	298.867	508.278	234.199
14	321.857	547.376	252.214
15	344.847	586.475	270.229