

Energy-Dispersive XRF

Excitation and Filter Guide

1	2	3	4	5	6	7	8	9	10
H HYDROGEN 1.008	He HELIUM 4.003	Li LITHIUM 6.941	Be BERYLLIUM 9.012	B BORON 10.81	C CARBON 12.011	N NITROGEN 14.01	O OXYGEN 16.00	F FLUORINE 19.00	Ne NEON 20.18
11 Na SODIUM 22.99	12 Mg MAGNESIUM 24.31	13 Al ALUMINUM 26.98	14 Si SILICON 28.09	15 P PHOSPHOROUS 30.97	16 S SULFUR 32.06	17 Cl CHLORINE 35.45	18 Ar ARGON 39.95	19 K POTASSIUM 39.10	20 Ca CALCIUM 40.08
21 Sc SCANDIUM 44.96	22 Ti TITANIUM 47.90	23 V VANADIUM 50.94	24 Cr CHROMIUM 52.00	25 Mn MANGANESE 54.94	26 Fe IRON 55.85	27 Co COBALT 58.93	28 Ni NICKEL 58.70	29 Cu COPPER 63.55	30 Zn ZINC 65.38
31 Ga GALLIUM 69.72	32 Ge GERMANIUM 72.59	33 As ARSENIC 74.92	34 Se SELENIUM 78.96	35 Br BROMINE 79.90	36 Kr KRYPTON 83.80	37 Rb RUBIDIUM 85.47	38 Sr STRONTIUM 87.62	39 Y YTTORIUM 88.91	40 Zr ZIRCONIUM 91.22
41 Nb NIOBIUM 92.91	42 Mo MOLYBDENUM 95.94	43 Tc TECHNETIUM (98)	44 Ru RUTHENIUM 101.07	45 Rh RHODIUM 102.91	46 Pd PALLADIUM 106.4	47 Ag SILVER 107.87	48 Cd CADMIUM 112.41	49 In INDIUM 114.82	50 Sn TIN 118.69
51 Sb ANTIMONY 121.75	52 Te TELLURIUM 127.6	53 I IODINE 126.90	54 Xe XENON 131.3	55 Cs CESIUM 132.91	56 Ba BARIUM 137.33	57 La LANTHANUM 138.91	58 Ce CERIUM 140.12	59 Pr PRASEODYMIUM 140.91	60 Nd NEODYMIUM 144.24
61 Pm PROMETHIUM (145)	62 Sm SAMARIUM 150.4	63 Eu EUROPIUM 151.96	64 Gd GADOLINIUM 157.25	65 Tb TERBIUM 158.93	66 Dy DYSPROSIUM 162.5	67 Ho HOLMIUM 164.93	68 Er ERBIUM 167.26	69 Tm THULIUM 168.93	70 Yb YTTERIUM 173.04
71 Lu LUTETIUM 174.96	72 Hf HAFNIUM 178.49	73 Ta TANTALUM 180.95	74 W TUNGSTEN 183.85	75 Re RHENIUM 186.21	76 Os OSMIUM 190.2	77 Ir IRIDIUM 192.22	78 Pt PLATINUM 195.09	79 Au GOLD 196.97	80 Hg MERCURY 200.59
81 Tl THALLIUM 204.37	82 Pb LEAD 207.2	83 Bi BISMUTH 208.98	84 Po POLONIUM (209)	85 At ASTATINE (210)	86 Rn RADON (222)	87 Fr FRANCIUM (223)	88 Ra RADIUM 226.03	89 Ac ACTINIUM (227)	90 Th THORIUM 232.04
91 Pa PROTACTINIUM 231.04	92 U URANIUM 238.03	93 Np NEPTUNIUM 237.05	94 Pu PLUTONIUM (244)	95 Am AMERICIUM (243)	96 Cm CURIUM (247)	97 Bk BERKELIUM (247)	98 Cf CALIFORNIUM (251)	99 Es EINSTEINIUM (254)	100 Fm FERMIUM (257)
101 Md MENDELEVIUM (258)	102 No NOBELIUM (259)	103 Lr LAWRENCIUM (260)	104 Rf RUFORDIUM (261)	105 Db DUBNIUM (262)	106 Sg SEABORGIUM (263)	107 Bh BOHRIUM (264)	108 Hs HASSIUM (265)	109 Mt MOSCOVIUM (266)	110 Ds DARMSTADTIUM (267)
111 Cn COCHIN (268)	112 Fl FLEROVIUM (269)	113 Mc MOSCOVIUM (270)	114 Lv LIVERMORIUM (271)	115 Ts TENNESSIUM (272)	116 Og OGANESSIUM (273)	117 Bh BOHRIUM (274)	118 Hs HASSIUM (275)	119 Mt MOSCOVIUM (276)	120 Ds DARMSTADTIUM (277)

47 **K L**

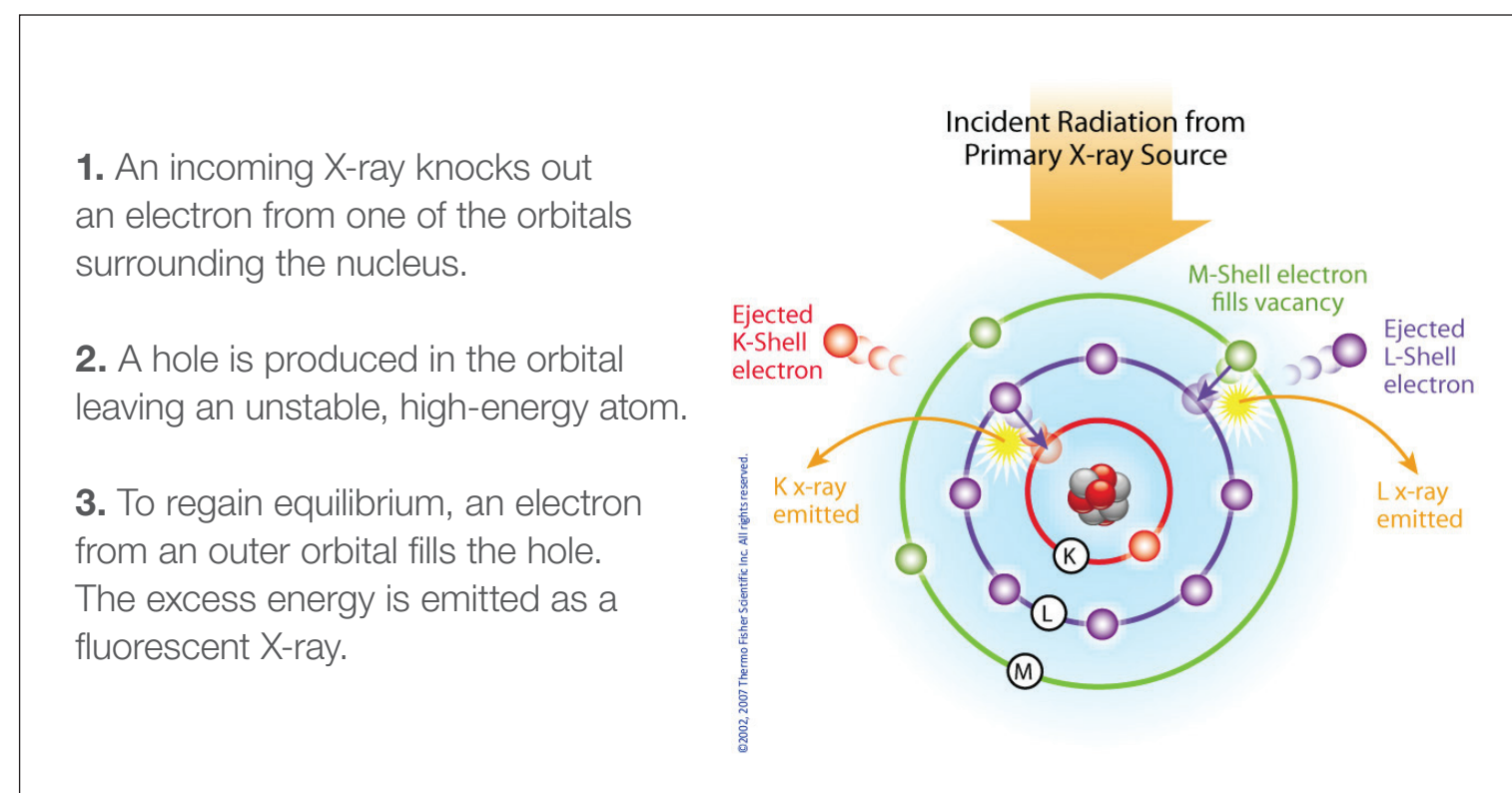
Symbol: **K**
Element Name: **L**
Atomic Weight: **107.87**

K α wtd. avg. (keV): **22.104**
K β wtd. avg. (keV): **24.987**
K abs. edge (keV): **25.517**

L α 1 (keV): **2.984**
L β 1 (keV): **3.151**
LII abs. edge (keV): **3.528**

OPTIMIZED FILTER: **K L**
ALTERNATE FILTER: **K L**

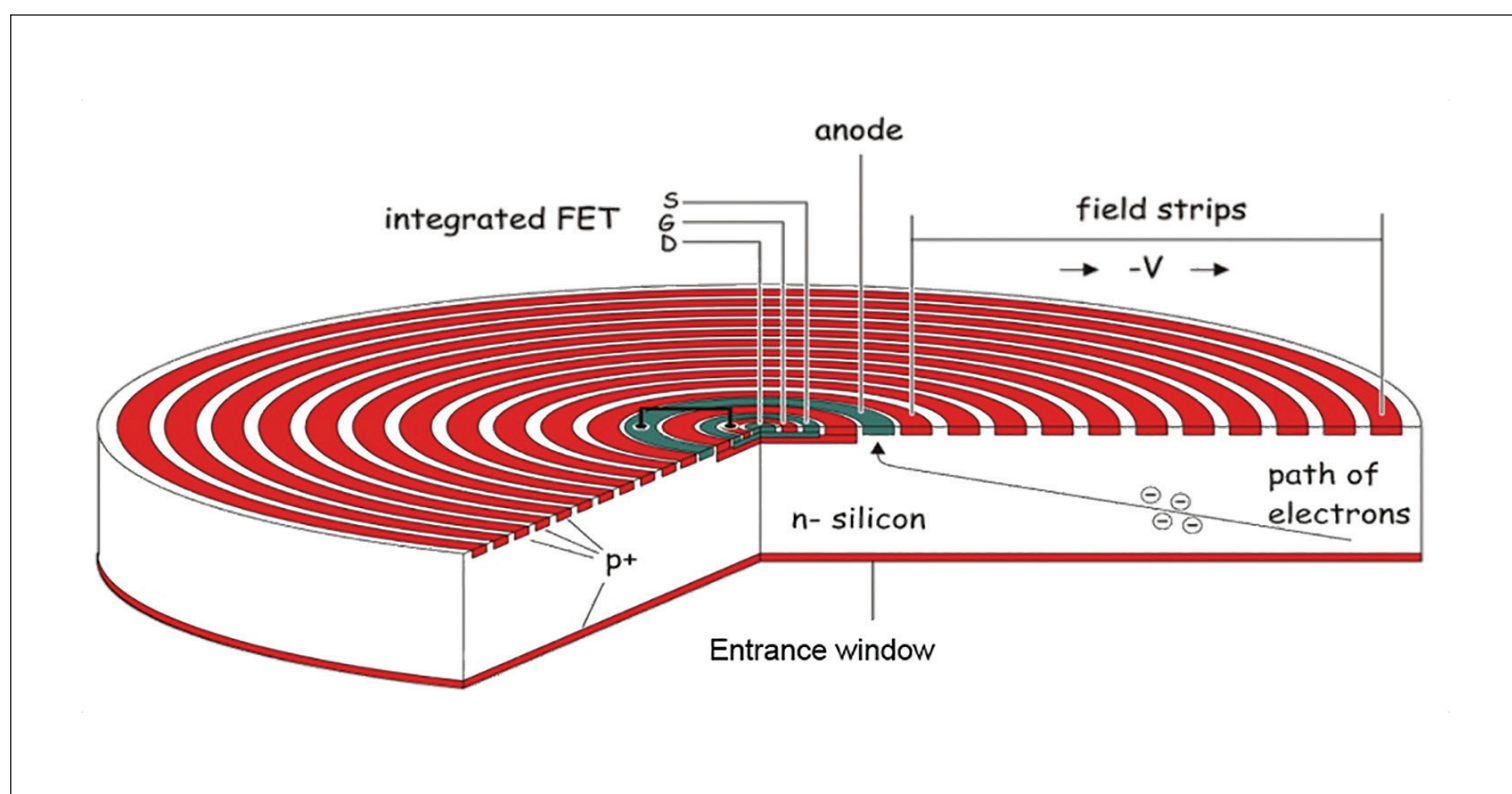
Color Code	Condition Name	Filter Number	Default kV	Optimized Atmosphere
Yellow	Low Za	0	4	Vacuum (solids) / He (liquids)
Light Yellow	Low Za II	1	8	Vacuum (solids) / He (liquids)
Orange	Low Zc	2	8	Vacuum (solids) / He (liquids)
Light Green	Low Zc	3	12	Air
Green	Mid Za	4	16	Air
Light Blue	Mid Zb	5	20	Air
Blue	Mid Zc	6	28	Air
Dark Blue	High Za	7	40	Air
Black	High Zb	8	50	Air



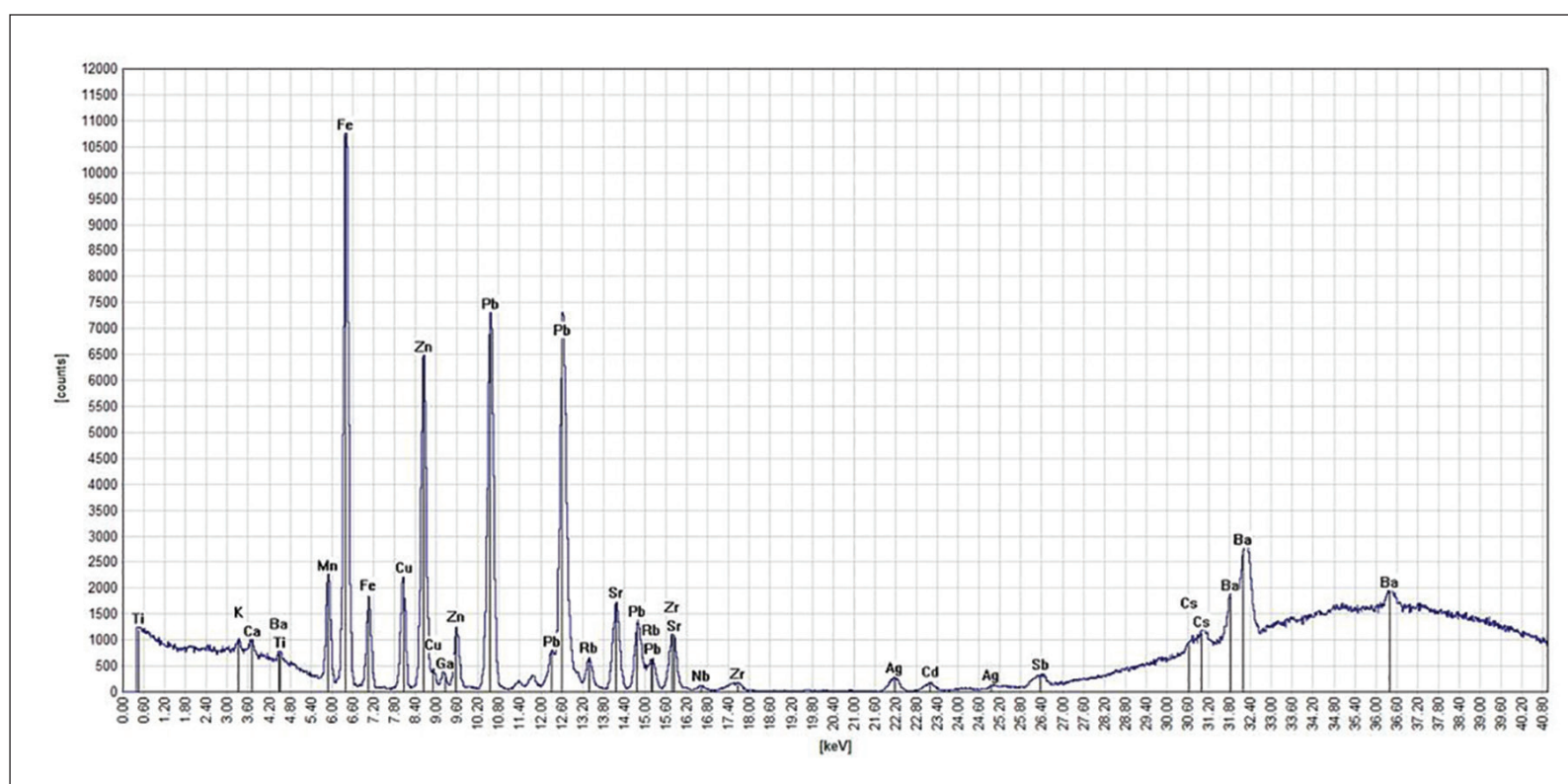
Principle of X-ray fluorescence



Different components of the EDXRF spectrometer



Detection principle of the silicon drift detector (SDD)



EDXRF detection of hazardous elements in NIST SRM 2710 Montana soil