

Spark optical emission spectrometry (OES) periodic chart

Elements and matrices analyzable by OES

1	<table border="1"> <tr> <td>1</td><td>H HYDROGEN</td><td colspan="14"></td><td>2</td> </tr> <tr> <td>2</td><td>Li LITHIUM</td><td>Be BERYLLIUM</td><td colspan="12"></td><td>B BORON</td><td>C CARBON</td><td>N NITROGEN</td><td>O OXYGEN</td><td>F FLUORINE</td><td>Ne NEON</td> </tr> <tr> <td>3</td><td>Na SODIUM</td><td>Mg MAGNESIUM</td><td colspan="12"></td><td>Al ALUMINUM</td><td>Si SILICON</td><td>P PHOSPHORUS</td><td>S SULFUR</td><td>Cl CHLORINE</td><td>Ar ARGON</td> </tr> <tr> <td>4</td><td>K POTASSIUM</td><td>Ca CALCIUM</td><td>Sc SCANDIUM</td><td>Ti TITANIUM</td><td>V VANADIUM</td><td>Cr CHROMIUM</td><td>Mn MANGANESE</td><td>Fe IRON</td><td>Co COBALT</td><td>Ni NICKEL</td><td>Cu COPPER</td><td>Zn ZINC</td><td>Ga GALLIUM</td><td>Ge GERMANIUM</td><td>As ARSENIC</td><td>Se SELENIUM</td><td>Br BROMINE</td><td>Kr KRYPTON</td> </tr> <tr> <td>5</td><td>Rb RUBIDIUM</td><td>Sr STRONTIUM</td><td>Y YTTORIUM</td><td>Zr ZIRCONIUM</td><td>Nb NIOBIUM</td><td>Mo MOLYBDENUM</td><td>Tc TECHNETIUM</td><td>Ru RUTHENIUM</td><td>Rh RHODIUM</td><td>Pd PALLADIUM</td><td>Ag SILVER</td><td>Cd CADMIUM</td><td>In INDIUM</td><td>Sn TIN</td><td>Sb ANTIMONY</td><td>Te TELLURIUM</td><td>I IODINE</td><td>Xe XENON</td> </tr> <tr> <td>6</td><td>Cs CESIUM</td><td>Ba BARIUM</td><td>La LANTHANUM</td><td>Hf HAFNIUM</td><td>Ta TANTALUM</td><td>W TUNGSTEN</td><td>Re RHENIUM</td><td>Os OSMIUM</td><td>Ir IRIDIUM</td><td>Pt PLATINUM</td><td>Au GOLD</td><td>Hg MERCURY</td><td>Tl THALLIUM</td><td>Pb LEAD</td><td>Bi BISMUTH</td><td>Po POLONIUM</td><td>At ASTATINE</td><td>Rn RADON</td> </tr> <tr> <td>7</td><td>Fr FRANCIUM</td><td>Ra RADIUM</td><td>Ac ACTINIUM</td><td colspan="14"></td> </tr> </table>																1	H HYDROGEN															2	2	Li LITHIUM	Be BERYLLIUM													B BORON	C CARBON	N NITROGEN	O OXYGEN	F FLUORINE	Ne NEON	3	Na SODIUM	Mg MAGNESIUM													Al ALUMINUM	Si SILICON	P PHOSPHORUS	S SULFUR	Cl CHLORINE	Ar ARGON	4	K POTASSIUM	Ca CALCIUM	Sc SCANDIUM	Ti TITANIUM	V VANADIUM	Cr CHROMIUM	Mn MANGANESE	Fe IRON	Co COBALT	Ni NICKEL	Cu COPPER	Zn ZINC	Ga GALLIUM	Ge GERMANIUM	As ARSENIC	Se SELENIUM	Br BROMINE	Kr KRYPTON	5	Rb RUBIDIUM	Sr STRONTIUM	Y YTTORIUM	Zr ZIRCONIUM	Nb NIOBIUM	Mo MOLYBDENUM	Tc TECHNETIUM	Ru RUTHENIUM	Rh RHODIUM	Pd PALLADIUM	Ag SILVER	Cd CADMIUM	In INDIUM	Sn TIN	Sb ANTIMONY	Te TELLURIUM	I IODINE	Xe XENON	6	Cs CESIUM	Ba BARIUM	La LANTHANUM	Hf HAFNIUM	Ta TANTALUM	W TUNGSTEN	Re RHENIUM	Os OSMIUM	Ir IRIDIUM	Pt PLATINUM	Au GOLD	Hg MERCURY	Tl THALLIUM	Pb LEAD	Bi BISMUTH	Po POLONIUM	At ASTATINE	Rn RADON	7	Fr FRANCIUM	Ra RADIUM	Ac ACTINIUM															1
1	H HYDROGEN															2																																																																																																																																							
2	Li LITHIUM	Be BERYLLIUM													B BORON	C CARBON	N NITROGEN	O OXYGEN	F FLUORINE	Ne NEON																																																																																																																																			
3	Na SODIUM	Mg MAGNESIUM													Al ALUMINUM	Si SILICON	P PHOSPHORUS	S SULFUR	Cl CHLORINE	Ar ARGON																																																																																																																																			
4	K POTASSIUM	Ca CALCIUM	Sc SCANDIUM	Ti TITANIUM	V VANADIUM	Cr CHROMIUM	Mn MANGANESE	Fe IRON	Co COBALT	Ni NICKEL	Cu COPPER	Zn ZINC	Ga GALLIUM	Ge GERMANIUM	As ARSENIC	Se SELENIUM	Br BROMINE	Kr KRYPTON																																																																																																																																					
5	Rb RUBIDIUM	Sr STRONTIUM	Y YTTORIUM	Zr ZIRCONIUM	Nb NIOBIUM	Mo MOLYBDENUM	Tc TECHNETIUM	Ru RUTHENIUM	Rh RHODIUM	Pd PALLADIUM	Ag SILVER	Cd CADMIUM	In INDIUM	Sn TIN	Sb ANTIMONY	Te TELLURIUM	I IODINE	Xe XENON																																																																																																																																					
6	Cs CESIUM	Ba BARIUM	La LANTHANUM	Hf HAFNIUM	Ta TANTALUM	W TUNGSTEN	Re RHENIUM	Os OSMIUM	Ir IRIDIUM	Pt PLATINUM	Au GOLD	Hg MERCURY	Tl THALLIUM	Pb LEAD	Bi BISMUTH	Po POLONIUM	At ASTATINE	Rn RADON																																																																																																																																					
7	Fr FRANCIUM	Ra RADIUM	Ac ACTINIUM																																																																																																																																																				
6	<table border="1"> <tr> <td>6</td><td>Ce CERIUM</td><td>Pr PRASEODYMIUM</td><td>Nd NEODYMIUM</td><td>Pm PROMETHIUM</td><td>Sm SAMARIUM</td><td>Eu EUROPIUM</td><td>Gd GADOLINIUM</td><td>Tb TERBIUM</td><td>Dy DYSPROSIUM</td><td>Ho HOLMIUM</td><td>Er ERBIUM</td><td>Tm THULIUM</td><td>Yb YTTERIUM</td><td>Lu LUTETIUM</td> </tr> <tr> <td>7</td><td>Th THORIUM</td><td>Pa PROTACTINIUM</td><td>U URANIUM</td><td>Np NEPTUNIUM</td><td>Pu PLUTONIUM</td><td>Am AMERICIUM</td><td>Cm CURIUM</td><td>Bk BERKELIUM</td><td>Cf CALIFORNIUM</td><td>Es EINSTEINIUM</td><td>Fm FERMIUM</td><td>Md MENDELEVIUM</td><td>No NOBELIUM</td><td>Lr LAWRENCIUM</td> </tr> </table>																6	Ce CERIUM	Pr PRASEODYMIUM	Nd NEODYMIUM	Pm PROMETHIUM	Sm SAMARIUM	Eu EUROPIUM	Gd GADOLINIUM	Tb TERBIUM	Dy DYSPROSIUM	Ho HOLMIUM	Er ERBIUM	Tm THULIUM	Yb YTTERIUM	Lu LUTETIUM	7	Th THORIUM	Pa PROTACTINIUM	U URANIUM	Np NEPTUNIUM	Pu PLUTONIUM	Am AMERICIUM	Cm CURIUM	Bk BERKELIUM	Cf CALIFORNIUM	Es EINSTEINIUM	Fm FERMIUM	Md MENDELEVIUM	No NOBELIUM	Lr LAWRENCIUM	6																																																																																																								
6	Ce CERIUM	Pr PRASEODYMIUM	Nd NEODYMIUM	Pm PROMETHIUM	Sm SAMARIUM	Eu EUROPIUM	Gd GADOLINIUM	Tb TERBIUM	Dy DYSPROSIUM	Ho HOLMIUM	Er ERBIUM	Tm THULIUM	Yb YTTERIUM	Lu LUTETIUM																																																																																																																																									
7	Th THORIUM	Pa PROTACTINIUM	U URANIUM	Np NEPTUNIUM	Pu PLUTONIUM	Am AMERICIUM	Cm CURIUM	Bk BERKELIUM	Cf CALIFORNIUM	Es EINSTEINIUM	Fm FERMIUM	Md MENDELEVIUM	No NOBELIUM	Lr LAWRENCIUM																																																																																																																																									
7																	7																																																																																																																																						

Element

M Routinely analyzed

M Less routinely analyzed in other matrices

- Not analyzed

☉ = Radioactive

ATOMIC NUMBER

ELEMENT

Au Mg Ni Pd Pt Sn Zn
Cu In Ir Rh Ru Zr

Element

M Routinely analyzed

M Less routinely analyzed in other matrices

M

1st line: other principal matrices

2nd line: other matrices

Principal matrix element

Matrix element

Potential matrix element

Not matrix element

Physical state at room temp.

E Solid

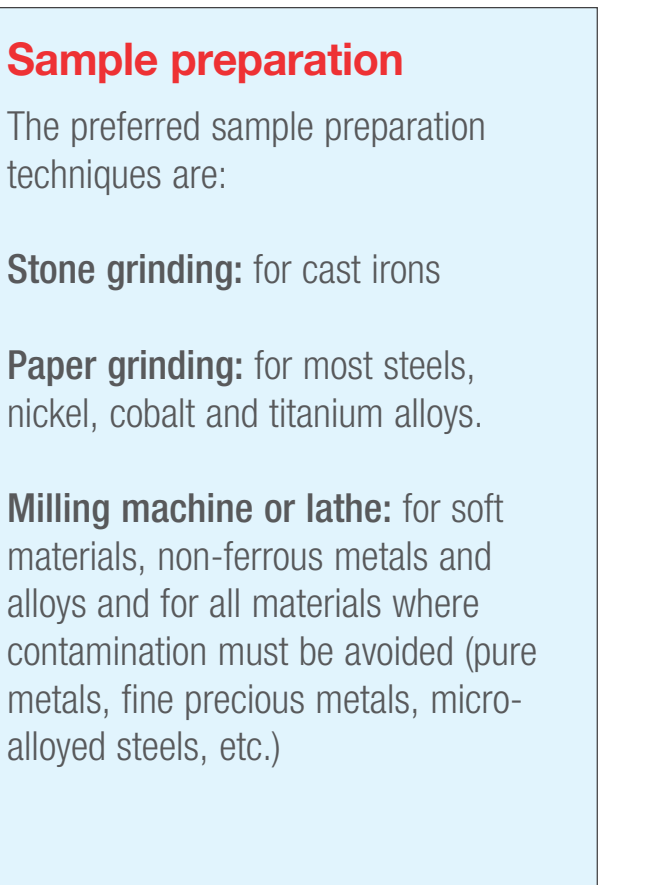
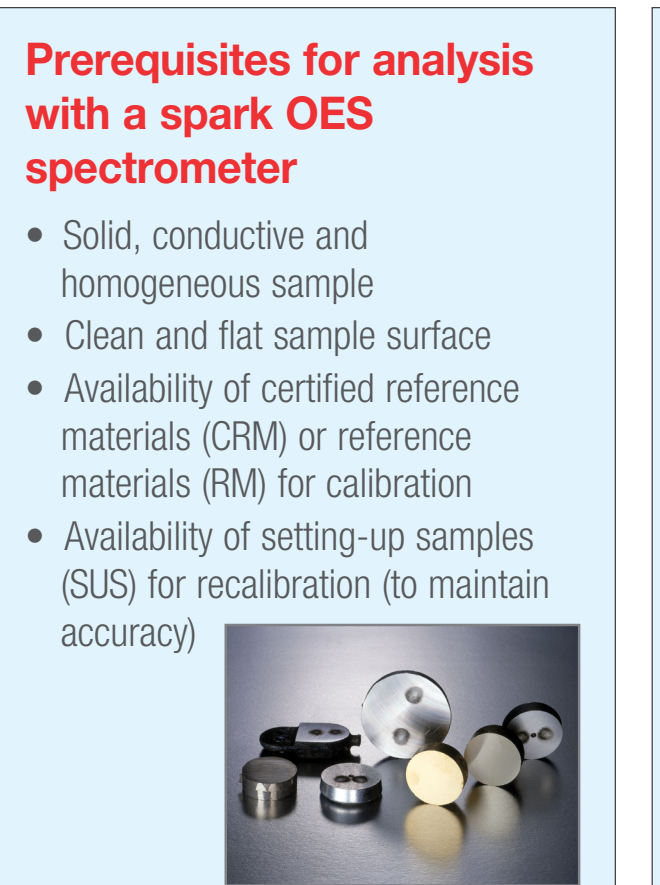
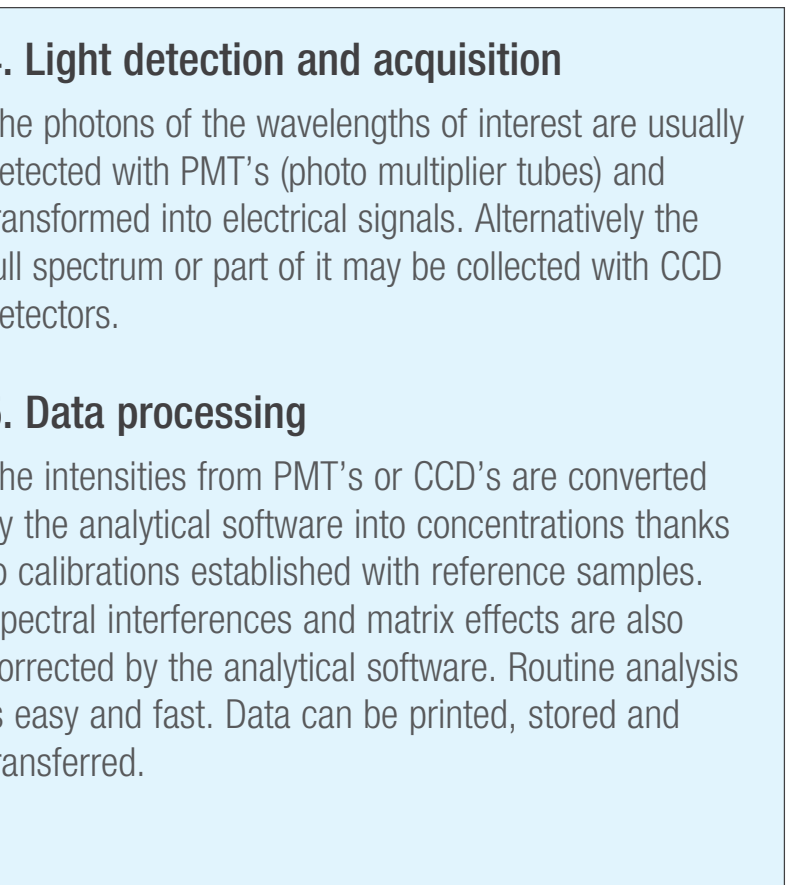
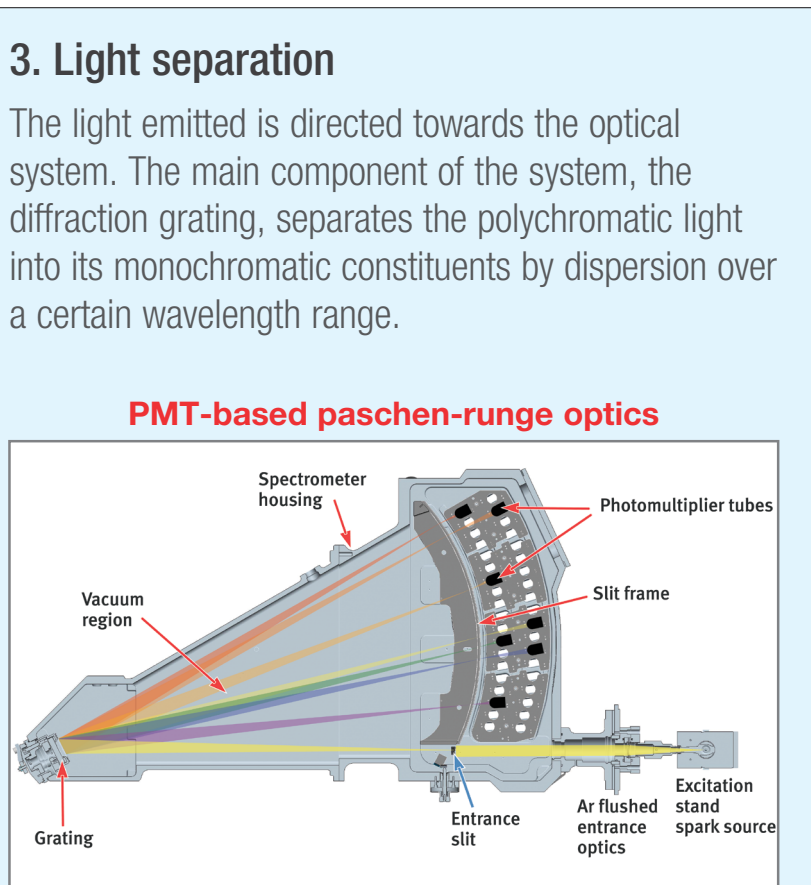
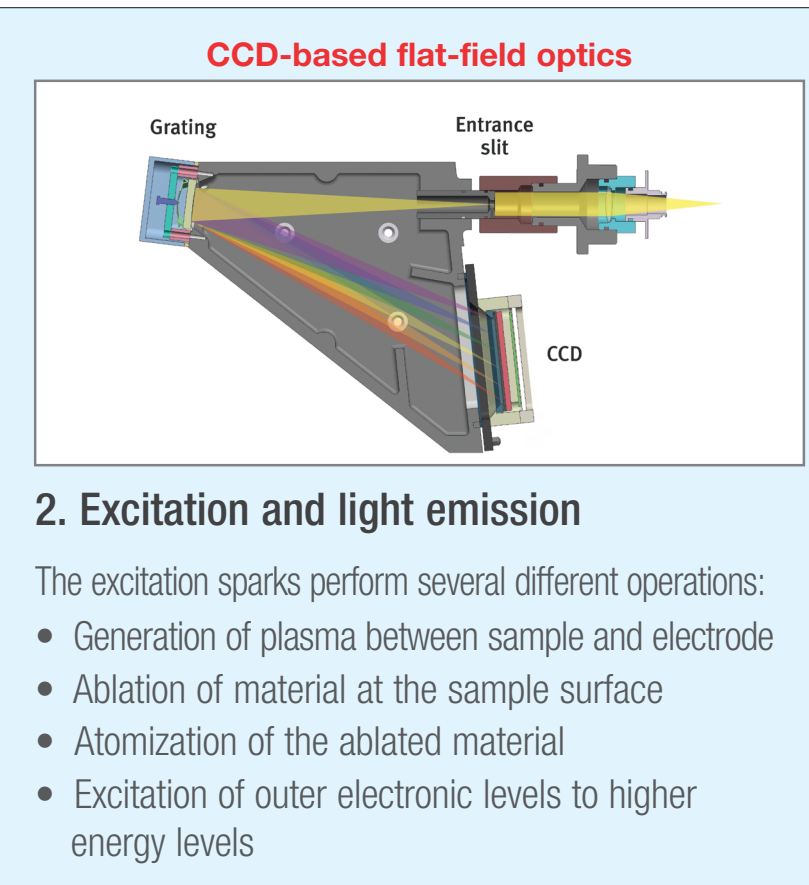
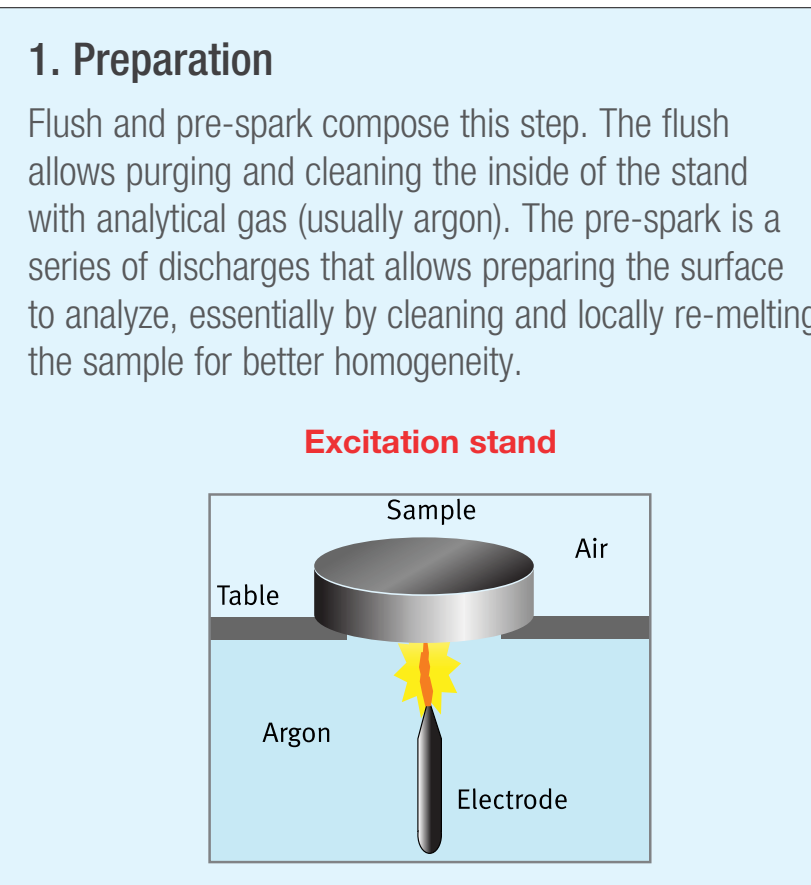
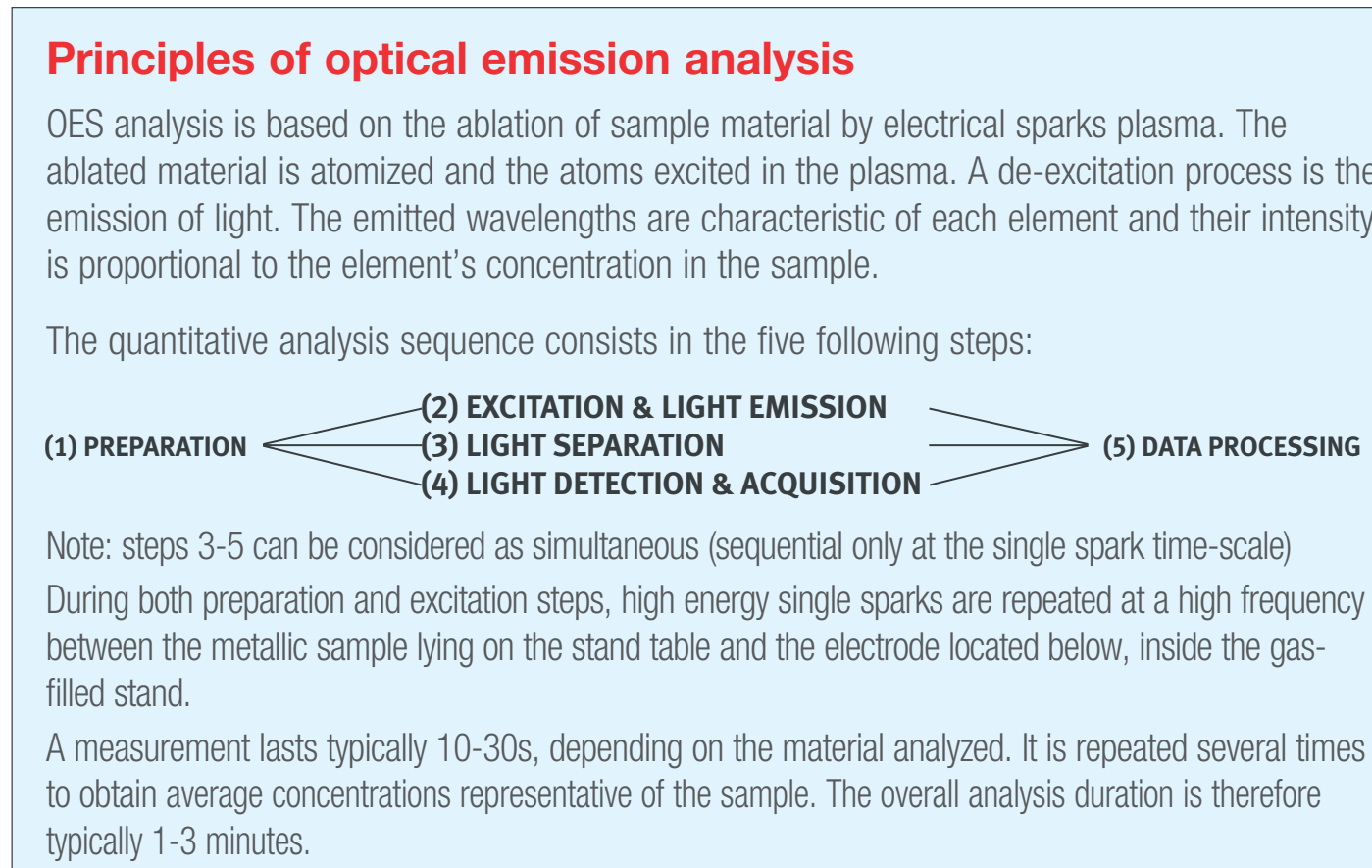
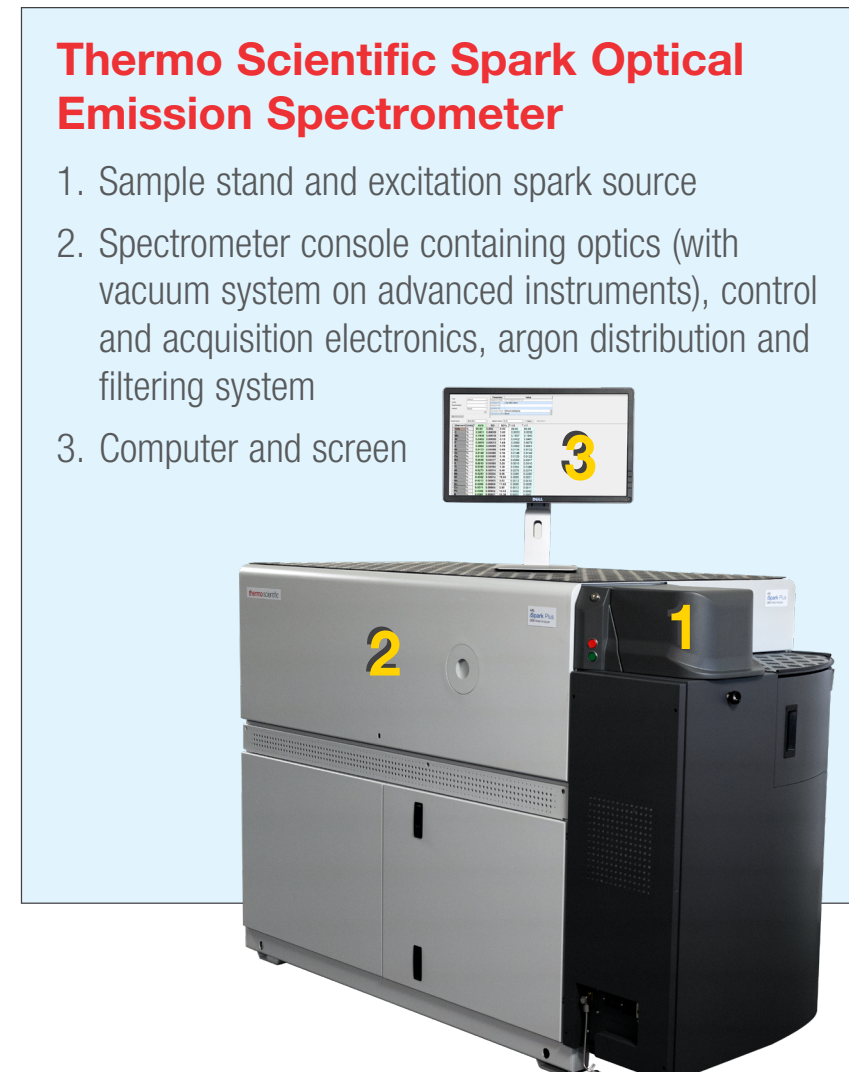
L Liquid

G Gas

☉ Non-natural (solid)

Principles of spark optical emission spectroscopy

Spark optical emission spectrometry (OES) is a form of atomic emission spectrometry (AES). OES is simple and fast, and needs little maintenance and consumables. It quantitatively analyzes most elements in solid metallic samples, from trace to percent levels. OES is therefore a preferred analysis technique in all the industries dealing with metals.



The content of this table is indicative, may depend on instrument and is subject to change. Contact your nearest Thermo Fisher Scientific sales representative for the latest and most accurate information.