



Magnetic Materials

Charge Transfer Complexes	182	Superconductor Precursors	187
Acceptors	182	Lanthanides	187
Donors	183	Other Metal Compounds	188
Other Donors & Acceptors	184	Anhydrous Metal Halides:	
Inorganic Magnetic Materials	185	<i>See Materials for Thin Films/</i>	
Metals	185	<i>Microelectronics</i>	255
Oxides	186	Molecular Based Superconductors:	
Salts	186	<i>See Charge Transfer Complexes</i>	182

An introductory overview of barrier materials, high κ dielectric materials, and atomic layer deposition (ALD) is provided in the Technical Appendix section.

Superconductors

Superconductors are materials which exhibit no electrical resistance below a certain temperature defined as the critical temperature (T_c). Prior to 1986, the highest T_c reported was 20 K for Nb_3Ge and Nb_3Sn .¹ In 1986–87, however, a group led by Johannes Bednorz and Karl Müller reported the ceramic oxides $YBa_2Cu_3O_7$ (Aldrich product **32,862-6**) and $La_{2-x}Ba_xCuO_{4-x}$ which superconduct above the boiling point of nitrogen (77K).^{2,3} Materials whose T_c s are greater than the boiling point of nitrogen (a common, readily available cryogenic coolant) are referred to as high-temperature superconductors (HTS). For their work, Bednorz and Müller were awarded the Nobel Prize in Physics in 1987.⁴

Other more exotic compounds such as fullerides have also exhibited superconducting properties. Fullerides of the formula A_xC_{60} ($A = K, Rb, Cs$) are reported to have superconducting character.⁵ Although superconductive compounds have been known for nearly a century, the relatively mundane compound magnesium boride has only recently been demonstrated to exhibit superconductivities. Magnesium boride, MgB_2 (Aldrich product **55,391-3**) is not only superconductive but its critical temperature is surprisingly high for a simple ceramic material ($T_c = 39$ K).⁶ Figure 1 shows an image of a MgB_2 wire segment with a tungsten boride core. The wire is formed by reaction of magnesium vapor with a boron filament. The grain structure seen in this image is visible under polarized light.⁷

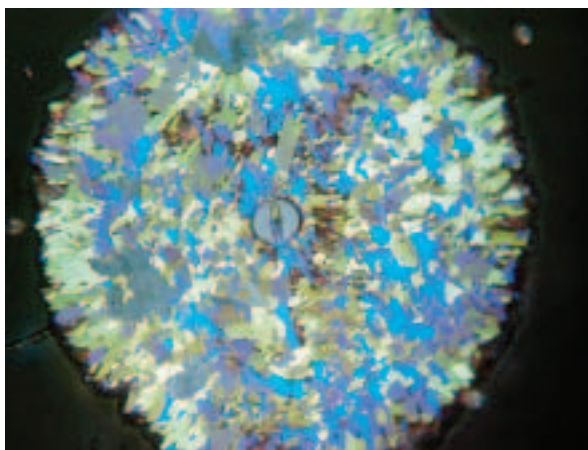


Figure 1. The cross section of a MgB_2 wire segment. (Image courtesy of D.K. Finnemore, S.L. Bud'ko, P.C. Canfield, Ames Laboratory, Iowa State University.)

Spintronics

Spintronics (short for spin-based electronics), sometimes called magnetoelectronics, is the term given to microelectronic devices that function by exploiting the spin of electrons. The most common use of spintronics today is in computer hard drives. Here memory storage is based on giant magnetoresistance (GMR), a spintronic effect. There is current research focussing on bringing magnetic random-access memory (MRAM) to market. Spintronic based MRAMs should rival the speed and rewritability of conventional RAM and retain their state (and thus memory) even when the power is turned off. Motorola has recently

developed a 256-kb MRAM (see Figure 2) based on a single magnetic tunnel junction and a single transistor. This MRAM has read/write cycles of less than 50 nanoseconds.

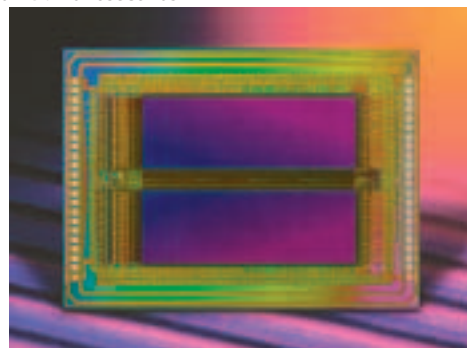


Figure 2. A 256-kb MRAM based on modern spintronics technology. (Image courtesy of Motorola Corp.)

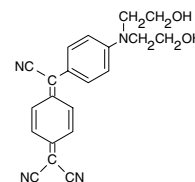
References (1) Galvalier, J.R. *App. Phys. Lett.* **1973**, *23*, 480. (2) Wu, M. K. et al. *Phys. Rev. Lett.* **1987**, *58*, 908. (3) Beno, M.A. et al. *Appl. Phys. Lett.* **1987**, *51*, 57. (4) <http://www.nobel.se/> (5) Hebard, A.F. et al. *Nature* **1991**, *350*, 600. (6) Nagamatsu, J. et al. *Nature* **2001**, *410*, 63. (7) Canfield, P.C. et al. *Phys. Res. Lett.* **2001**, *86*, 1877.

Charge Transfer Complexes

Acceptors

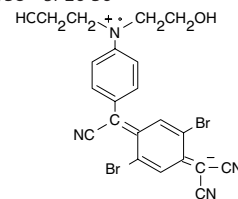
2-[4-((4-(Bis(2-hydroxyethyl)amino)phenyl)(cyano)methylene)-2,5-cyclohexadien-1-ylidene]malonitrile

57,210-1 CAS No. 287922-16-1 50 mg
 $C_{21}H_{18}N_4O_2$ FW 358.4 250 mg
98%
 R: 20/21/22-36/37/38 S: 26-36



2-[4-((4-(Bis(2-hydroxyethyl)amino)phenyl)(cyano)methylene)-2,5-dibromo-cyanohexadien-1-ylidene]malonitrile

57,650-6 CAS No. 287922-14-9 5 mg
 $C_{21}H_{16}Br_2N_4O_2$ FW 516.2 25 mg
98%
 R: 36/37/38 S: 26-36

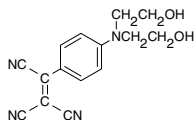


Charge Transfer Complexes

Acceptors

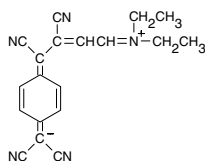
[4-[Bis(2-hydroxyethyl)amino]phenyl]-1,1,2-ethylenetri-carbonitrile

57,212-8 CAS No. 56672-91-4 50 mg
(HOCH₂CH₂)NC₆H₄C(CN)C(CN)₂ 250 mg
FW 282.3
98%
solid
 λ_{max} 514 nm
R: 36/37/38 S: 26-36



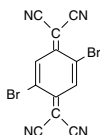
N-[3-Cyano-3-[4-(dicyanomethyl)phenyl]-2-propenylidene]-N-ethyl-ethaniminium inner salt

57,213-6 CAS No. 174280-29-6 50 mg
C₁₇H₁₆N₄ FW 276.3 250 mg
 λ_{max} 698 nm (lit.)
R: 36/37/38 S: 26-36



2,5-Dibromo-7,7,8,8-tetracyanoquinodimethane

57,209-8 CAS No. 56403-70-4 5 mg
C₁₂H₂Br₂N₄ FW 362.0 25 mg
98%
R: 36/37/38 S: 26-36



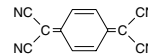
Tetracyanoethylene

T880-9 (Percyanoethylene; Ethylenetetracarbonitrile; 1 g
TCNE) 5 g
CAS No. 670-54-2 10 g
(NC)₂C:C(CN)₂ FW 128.1 25 g
98%
mp 197-199 (lit.)
Licensed under U.S. Patent 3 166 584
R: 25 S: 36/37/39-45



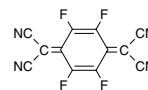
7,7,8,8-Tetracyanoquinodimethane

15,763-5 ((2,5-Cyclohexadiene-1,4-diylidene)- 1 g
dimalononitrile; TCNQ) 5 g
CAS No. 1518-16-7 10 g
C₁₂H₄N₄ FW 204.2
98%
Electron-acceptor molecule used to form charge-transfer superconductors. Williams, J.M. et al. *Organic Superconductors (Including Fullerenes): Synthesis, Structure, Properties, and Theory* Prentice-Hall: Englewood Cliffs, NJ, **1992** (Z24,430-9).
mp 287-289 °C (dec.) (lit.)
R: 22 S: 22-24/25



2,3,5,6-Tetrafluoro-7,7,8,8-tetracyanoquinodimethane

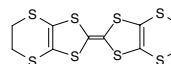
37,677-9 ((2,3,5,6-Tetrafluoro-2,5-cyclohexadiene- 5 mg
1,4-diylidene)dimalononitrile; 7,7,8,8- 25 mg
Tetracyano-2,3,5,6-
tetrafluoroquinodimethane)
CAS No. 29261-33-4
C₁₂F₄N₄ FW 276.1
97%
mp 291 °C (dec.) (lit.)
S: 22-24/25



Donors

Bis(ethylenedithio)tetrathiafulvalene

36,202-6 (Bi[5,6-dihydro-1,3-dithiolo[4,5-b][1,4]- 100 mg
dithiine-2-ylidene]; BEDT-TTF) 500 mg
CAS No. 66946-48-3
C₁₀H₈S₈ FW 384.7
98%
Used as a component for organic superconductors.
Aldrichimica Acta **1990**, 23, 54.
mp 244 °C (dec.) (lit.)
S: 22-24/25



Charge Transfer Complexes

Donors

Tetrathiafulvalene

18,318-0 (TTF; $\Delta 2,2'$ -Bi-1,3-dithiole) 100 mg
 CAS No. 31366-25-3 250 mg
 $C_6H_4S_4$ FW 204.4 1 g
97%
 Electron donor for supramolecular synthesis,¹ charge-transfer complex synthesis² with 7,7,8,8-tetracyanoquinodimethane (15,763-5), and for electron transfer to diazonium salts.³ (1) *Chem. Soc. Rev.* **1994**, 23, 41. (2) *Aldrichimica Acta* **1985**, 18, 73. (3) *Tetrahedron Lett.* **1994**, 35, 8675.
 mp 120-123 °C (lit.)
 R: 43 S: 36/37



Other Donors & Acceptors

p-Benzoquinone

B1,035-8 (Quinone) 5 g
 CAS No. 106-51-4 100 g
 $C_6H_4O_2$ FW 108.1 500 g
98% 1 kg
 Free radical inhibitor
 Vapor density 3.73 (vs air)
 Vapor pressure 0.1 mm Hg (25 °C)
 mp 113-115 °C (lit.)
 R: 23/25-36/37/38-50 S: 26-28-45-61

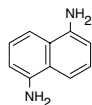


Chloranil

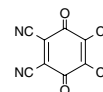
23280 CAS No. 118-75-2 10 g
 $C_6Cl_4O_2$ FW 245.9 50 g
puriss., ≥99.0% (AT) 250 g
 R: 36/38-50/53 S: 37-60-61

1,5-Diaminonaphthalene

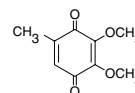
D2,120-0 (1,5-Naphthalenediamine) 25 g
 CAS No. 2243-62-1 100 g
 $C_{10}H_{10}N_2$ FW 158.2
97%
 mp 185-187 °C (lit.)
 R: 40-50/53 S: 36/37-60-61

2,3-Dichloro-5,6-dicyano-*p*-benzoquinone

D6,040-0 (4,5-Dichloro-3,6-dioxo-1,4-cyclohexadiene-1,2-dicarbonitrile; DDQ) 5 g
 CAS No. 84-58-2 10 g
 $C_8Cl_2N_2O_2$ FW 227.0 100 g
98%
 mp 213-216 °C (lit.)
 R: 25-29 S: 22-24/25-37-45

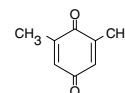
2,3-Dimethoxy-5-methyl-*p*-benzoquinone

29,956-1 (Coenzyme Q₀) 1 g
 CAS No. 605-94-7 5 g
 $C_9H_{10}O_4$ FW 182.2
99%
 mp 59-60 °C (lit.)
 R: 36/37/38 S: 26-36



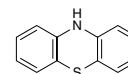
2,6-Dimethylbenzoquinone

D14,970-5 CAS No. 527-61-7 1 g
 $C_8H_8O_2$ FW 136.1 5 g
99%
 mp 71-73 °C (lit.)



Phenothiazine

P1,483-1 CAS No. 92-84-2 25 g
 $C_{12}H_9NS$ FW 199.3 500 g
98+% 1 kg
 Employed in the preparation of carbazoles and piperazines,¹ and charge-transfer semiconducting complexes.² (1) *Indian J. Chem., Sect. B: Org. Chem. Incl. Med. Chem.* **1994**, 33B, 397. (2) *Mol. Cryst. Liq. Cryst. Sci. Technol., Sect. A* **1993**, 237, 419.
 bp 371 °C/760 mm Hg (lit.)
 mp 182-187 °C (lit.)
 R: 36/37/38-43 S: 26-36

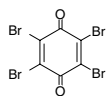


Charge Transfer Complexes

Other Donors & Acceptors

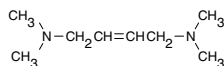
Tetrabromo-1,4-benzoquinone

34,350-1 CAS No. 488-48-2 5 g
 $C_6Br_4O_2$ FW 423.7
90%, Technical Grade
 mp 292-294 °C (lit.)
 R: 36/37/38 S: 26-37/39



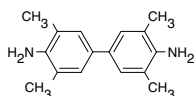
Tetrakis(dimethylamino)ethylene

23,423-0 (Octamethyl-ethylenetetramine) 25 g
 CAS No. 996-70-3
 $C_{10}H_{24}N_4$ FW 200.3
95%
 To prolong shelf-life, store under argon!
 n_D^{20} 1.48 (lit.)
 bp. 59 °C/0.9 mm Hg (lit.)
 Fp. 128 °F
 Density. 0.861 g/mL (lit.)
 R: 10-34 S: 26-36/37/39-45



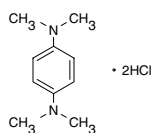
3,3',5,5'-Tetramethylbenzidine

86,033-6 (TMB) 100 mg
 CAS No. 54827-17-7 1 g
 $C_{16}H_{20}N_2$ FW 240.3 5 g
99+% 25 g
 mp 168-169 °C (lit.)
 R: 36/37/38 S: 26-36



N,N,N',N'-Tetramethyl-*p*-phenylenediamine dihydrochloride

23,469-9 (Wurster's Reagent dihydrochloride; 1 g
 Wurster's reagent) 5 g
 CAS No. 637-01-4 25 g
 $C_{10}H_{16}N_2 \cdot 2HCl$ FW 237.2
98%
 mp 222-224 °C (lit.)
 R: 36/37/38 S: 26-36



Inorganic Magnetic Materials

Metals

Cerium

CAS No. 7440-45-1
 Ce FW 140.1
 R: 20/21/22-36/37/38 S: 16-26-36/37/39

46,121-0 chips, 99.9% 10 g
 Packaged under argon 50 g
 mp 795 °C (lit.)
 Density. 6.67 g/mL (lit.)

26,104-1 ingot, 99.9% 25 g
 Packaged under oil 100 g

26,300-1 99.9%, powder, -40 mesh 10 g
 Packaged under oil and nitrogen 50 g

Iron

CAS No. 7439-89-6
 Fe FW 55.85

31,035-2 granules, low in magnesium and manganese, ACS reagent 100 g
 500 g
 Mg. ≤5 ppm
 Mn. ≤0.002%
 R: 11 S: 16-33

46,037-0 foil, thickness 0.127 mm, 99.5% 2.6 g
 R: 25-36 S: 26-36/37/39-45 10.4 g

Neodymium

46,087-7 CAS No. 7440-00-8 10 g
 Nd FW 144.2 50 g
powder, -40 mesh, 99+% (Purity based on trace metals analysis)
 Packaged under argon in ampules
 Density. 7.003 g/mL (lit.)
 R: 11-36/37/38 S: 16-26-33-36/37/39

Nickel

CAS No. 7440-02-0
 Ni FW 58.69

52,035-7 flake, 20 μm 500 g
 R: 10-17-36/37/38-40-42/43 S: 16-26-33-36

26,828-3 powder, (submicron), 99.8% 25 g
 Average particle size determined by Fisher 100 g
 sub-sieve sizer
 R: 10-17-36/37/38-40-42/43 S: 16-26-33-36

26,698-1 powder, 3 μm, 99.7% 100 g
 Average particle size determined by Fisher 500 g
 sub-sieve sizer
 Density. 8.9 g/mL (lit.)
 R: 10-17-36/37/38-40-42/43 S: 16-26-33-36

26,697-3 powder, -100 mesh, 99% 100 g
 R: 10-17-36/37/38-40-42/43 S: 16-26-33-36 500 g

26,824-0 wire, diam. 0.5 mm, 99+% 35 g
 175 g of 100 m 175 g
 35 g of 20 m
 R: 40-43 S: 22-36

Ordering/Pricing: www.sigma-aldrich.com 800.558.9196 (USA) Your local Sigma-Aldrich office: See Back Cover

Magnetic Materials

ALDRICH

Inorganic Magnetic Materials

Oxides

Ferric hydroxide oxide

(Goethite)
CAS No. 20344-49-4
Fe(OH)O FW 88.85

54,626-7	crystalline powder	50 g
		250 g
37,125-4	catalyst grade, 30-50 mesh	50 g
		250 g

Iron(II,III) oxide

CAS No. 1317-61-9
Fe₃O₄ FW 231.5
R: 20/21/22-36/37/38 S: 26-36/37/39-45

51,815-8	99.99%	10 g
		50 g
31,006-9	powder, 5 μm, 98%	25 g
		500 g
		2.5 kg

Iron(II) oxide

40,086-6	CAS No. 1345-25-1	5 g
	FeO FW 71.84	25 g
	-10 mesh, 99.9%	
	Employed in the synthesis of the first example of a mixed-valence ammonium iron phosphate, NH ₄ Fe ₂ (PO ₄) ₂ . Boudin, S. Lii, K-H. <i>Inorg. Chem.</i> 1998 , <i>37</i> , 799.	
	Density.	5.7 g/mL (lit.)
	R: 11 S: 16-33-7/9	

Iron(III) oxide

(Ferric oxide)
CAS No. 1309-37-1
Fe₂O₃ FW 159.7
Density. 5.24 g/mL (lit.)
R: 36/37/38 S: 36/37/39

52,931-1	99.999%	5 g
	Fp.	>230 °F 25 g
25,572-6	99.998%	10 g
		50 g
20,351-3	99.98%	10 g
		50 g
31,005-0	powder, 5 μm, 99+%	25 g
		500 g
		2.5 kg
48,066-5	99.9%	10 g
		50 g
34,300-5	pieces, 3-12 mm, 99.8%	50 g
	Suitable for vacuum deposition. sintered	

Neodymium oxide

CAS No. 1313-97-9
Nd₂O₃ FW 336.5
Density. 7.24 g/mL (lit.)

20,385-8	99.99%	10 g
		50 g
22,865-6	99.9%	25 g
		100 g

Salts

Erbium(III) bromide hydrate

57,522-4	CAS No. 29843-93-4	5 g
	ErBr ₃ ·xH ₂ O FW 407.0	25 g
	99.999%	
	bp.	1460 °C mm Hg
	mp.	923 °C
	R: 36/37/38 S: 26-36	

Iron(II) titanate

40,087-4	CAS No. 12022-71-8	25 g
	FeO ₃ Ti FW 151.7	100 g
	-100 mesh	
	R: 20/21/22-36/37/38-40 S: 22-26-36	

Neodymium bromide hydrate

54,505-8	CAS No. 29843-90-1	10 g
	Br ₃ Nd·xH ₂ O FW 384.0	50 g
	solid	
	Density.	5.3 g/mL (lit.)
	R: 36/37/38 S: 26-36	

Neodymium chloride hexahydrate

CAS No. 13477-89-9
NdCl₃·6H₂O FW 358.7

20,383-1	99.99%	20 g
	mp.	124 °C (lit.) 100 g
	R: 36/37/38 S: 26-37/39	
28,918-3	99.9%	25 g
	R: 36/37/38 S: 26-36	100 g

Neodymium(III) phosphate hydrate

58,709-5	CAS No. 14298-32-9	5 g
	NdO ₄ P FW 239.2	
	99.99+%	
	R: 36/37/38 S: 26-36	

Neodymium(III) titanate

40,369-5	Nd ₄ O ₁₆ Ti ₅ FW 1072	100 g
	-325 mesh	500 g

Potassium hexacyanocobaltate(III)

21,864-2	(Tripotassium hexacyanocobaltate(III))	5 g
	CAS No. 13963-58-1	100 g
	K ₃ Co(CN) ₆ FW 332.3	500 g
	Applied in formation of a xerogel via a cyanogel the gel has potential use in solid-state gas sensors. Sharp, S.L. et al. <i>Chem. Mater.</i> 1998 , <i>10</i> , 880.	
	Density.	1.878 g/mL (lit.)
	R: 23/24/25 S: 28-36/37-45-7/9	

Potassium hexacyanoferrate(III)

(Potassium ferricyanide; Red prussiate)
CAS No. 13746-66-2
K₃Fe(CN)₆ FW 329.2
R: 32-36/37/38 S: 26-36

45,594-6	≥99.99%	25 g
	Density.	1.89 g/mL (lit.) 100 g

Inorganic Magnetic Materials

Salts

39,351-7	99+%, ACS reagent, powder, 10 μm	5 g
	ferro compounds.	≤0.05% 100 g
	insolubles.	≤0.005% 500 g
	Cl ⁻	≤0.01%
	SO ₄ ²⁻	≤0.01%
24,402-3	≥99.0%	5 g
	ACS reagent	100 g
	insolubles.	≤0.005% 500 g
	Density.	1.89 g/mL (lit.)
20,801-9	99%	500 g
	Density.	1.85 g/mL (lit.) 2 kg

Potassium hexacyanoruthenate(II)

37,823-2	CAS No. 314075-44-0	500 mg
	C ₆ H ₆ K ₄ N ₆ ORu FW 435.6	2.5 g
	R: 23/24/25-36/37/38 S: 22-26-36/37/39-45	

Sodium nitroferricyanide(III)

43,145-1	(Sodium pentacyanonitrosylferrate; Nitroprusside sodium; sodium nitroprusside; SNP)	10 g
	CAS No. 13755-38-9	50 g
	Na ₂ [Fe(CN) ₅ NO] · 2H ₂ O FW 297.9	
	99.99+%, ReagentPlus™	
	Meets A.C.S. reagent specifications.	
	insolubles.	≤0.01%
	total metallic impurities.	<100 ppm
	Density.	1.72 g/mL (lit.)
	Cl ⁻	≤0.02%
	SO ₄ ²⁻	Passes test
	™Trademark of Sigma-Aldrich Co.	
	R: 25-32 S: 22-37-45	

Magnetic Nanostructures See: Books Page 497

Superconductor Precursors

Lanthanides

Erbium(III) chloride hexahydrate

CAS No. 10025-75-9
Cl₃Er · 6H₂O FW 381.7
R: 36/37/38 S: 26-37/39

20,321-1	99.995%	20 g
		100 g
28,925-6	99.9%	25 g
		100 g

Erbium(III) nitrate pentahydrate

CAS No. 10031-51-3
ErN₃O₉ · 5H₂O FW 443.4
R: 8-37/38-41 S: 17-26-36/39

54,510-4	99.9999%	10 g
		50 g
29,816-6	99.9%	25 g
		100 g

Erbium oxide

CAS No. 12061-16-4
Er₂O₃ FW 382.5
R: 36/37/38 S: 26-36

20,323-8	99.99+%	5 g
	Density.	8.64 g/mL (lit.) 25 g
28,924-8	99.9%	5 g
		25 g
		100 g

Lanthanide oxides See: Metallic Oxides Page 453

Lanthanum oxide

20,355-6	CAS No. 1312-81-8	20 g
	La ₂ O ₃ FW 325.8	100 g
	99.999%	500 g
	Precursor to LAMOX fast ion conductors and superconductors.	
	Density.	6.51 g/mL (lit.)
	R: 36/37/38 S: 26-36	

Samarium oxide

CAS No. 12060-58-1
Sm₂O₃ FW 348.8
R: 33 S: 22-24/25

39,439-4	99.999%	5 g
	Radioactive for shipping purposes	25 g
22,867-2	99.9%	10 g
	Density.	8.347 g/mL (lit.) 100 g

Ytterbium(III) bromide hydrate

54,496-5	CAS No. 15163-03-8	5 g
	YbCl ₃ ·xH ₂ O FW 279.4	25 g
	mp	300 °C (lit.)

Yttrium See: Metals Page 346

Yttrium barium copper oxide

32,862-6	CAS No. 107539-20-8	10 g
	YBa ₂ Cu ₃ O _x , x ~ 6.7	50 g
	powder	
	R: 20/21/22-36/37/38 S: 22-26-28-36	

Yttrium chloride hexahydrate

CAS No. 10025-94-2
YCl₃ · 6H₂O FW 303.4
R: 36/37/38 S: 26-36

20,491-9	99.999%	10 g
	Density.	2.18 g/mL (lit.) 50 g
46,431-7	99.99%	25 g
	mp	100 °C (dec.) (lit.) 100 g
21,164-8	99.9%	50 g
		250 g

Yttrium hexaboride

33,326-3	CAS No. 12008-32-1	2 g
	YB ₆ FW 153.8	
	powder, -100 mesh	

Ordering/Pricing: www.sigma-aldrich.com 800.558.9196 (USA) Your local Sigma-Aldrich office: See Back Cover

Magnetic Materials

ALDRICH

Superconductor Precursors

Lanthanides

Yttrium(III) oxide

CAS No. 1314-36-9
 Y_2O_3
 Density. 5.01 g/mL (lit.)
 R: 36/37/38 S: 26-36

20,492-7	99.999%	2 g
		10 g
		50 g
20,516-8	99.99%	10 g
		50 g
		100 g
		250 g

Other Metal Compounds

Barium calcium tungsten oxide See: Metallic Oxides Page 454

Barium strontium tungsten oxide See: Metallic Oxides Page 454

Bismuth strontium calcium copper oxide See: Metallic Oxides Page 454

Lithium hydride

20,104-9	CAS No. 7580-67-8	5 g
		LiH FW 7.949
		100 g
		500 g
		mp 680 °C (lit.)
		Density. 0.82 g/mL (lit.)

R: 14-34 S: 16-26-27-36/37/39

Magnesium boride

55,391-3	CAS No. 12007-25-9	5 g
		MgB ₂ FW 45.93
		25 g
		Superconduction material with unique properties.
		References

Nature **418**, 760 (2002)
 S: 22-24/25

Molybdenum(VI) oxide

(Molybdenum trioxide)
 CAS No. 1313-27-5
 MoO_3 FW 143.9
 Precursor to LAMOX fast ion conductors and superconductors.^{1,2}
 mp 795 °C (lit.)
 Density. 4.692 g/mL (lit.)
References
 1. Lacorre, P., *Nature* **404**, 856 (2000)
 2. Arulraj, A., *Chem. Mater.* **14**, 2492 (2002)
 R: 36/37-48/20/22 S: 22-23

20,381-5	99.99%	5 g		
		Used in the solid state synthesis of a remarkable ternary, reduced molybdenum oxide, Pr ₄ Mo ₉ O ₁₈ , whose structure contains previously unknown Mo ₇ , Mo ₁₃ , and Mo ₁₉ clusters. The new cluster product is a small band gap semiconductor. Tortelier, J. Gougeon, P. <i>Inorg. Chem.</i> 1998 , <i>37</i> , 6229.		
		25 g		
		26,785-6	99.5+%, ACS reagent	100 g
		insol. dil. NH ₄ OH. ≤0.01%	500 g	
		AsO ₄ ³⁻ , PO ₄ ³⁻ and SiO ₂ ≤0.001%		
		Cl ⁻ ≤0.002%		
		NO ₃ ⁻ Passes test		
		PO ₄ ³⁻ ≤5 ppm		
		SO ₄ ²⁻ ≤0.02%		
Heavy metals. ≤0.005%				
NH ₄ ⁺ ≤0.002%				
22,181-3	10-20 μm, 99+%	100 g		
		500 g		

Potassium hydride

21,581-3	CAS No. 7693-26-7	75 g
		KH FW 40.11
		300 g

30 wt % dispersion in mineral oil
 R: 11-14/15-34 S: 16-26-27-36/37/39-45

Potassium tetrachloropalladate(II)

20,579-6	(Potassium palladium(II) chloride)	2 g
		CAS No. 10025-98-6

K₂PdCl₄ FW 326.4
 Used in the synthesis of semiconducting metal-containing polymers in which the polypyrrole backbone has a conformational energy minimum and is nearly planar. Reacts with bis(dithiolates) to metal-bis(dithiolates) with applications in laser Q-switch materials, optical CD recording media, bar code material and superconductivity.
 mp 105 °C (dec.) (lit.)
 Density. 2.67 g/mL (lit.)
References
 1. Mathis, M. et al., *Chem. Mater.* **10**, 3568 (1998)
 2. Cummings, S.D.; Eisenberg, R., *Inorg. Chim. Acta* **225** (1996)
 R: 36/37/38 S: 26-36

Potassium tetrachloroplatinate(II)

(Potassium platinum(II) chloride)
 CAS No. 10025-99-7
 K₂PtCl₄ FW 415.1
 Reacts with bis(dithiolates) to form metal-bis(dithiolates) with applications in laser Q-switch materials, liquid crystals, optical CD recording media, bar code materials and superconductivity. Reacts with capping polymers to form shape-control colloidal platinum nanoparticles for catalysis and photocatalysts.
References
 1. Ahmedi, T.S., *Science* **272**, 1924 (1996)
 2. Ahmedi, T.S., *Chem. Mater.* **8**, 1162 (1996)
 3. Cummings, S.D.; Eisenberg, R., *Inorg. Chem.* **225** (1996)
 R: 25-34-42/43 S: 22-26-36/37/39-45

32,341-1	99.99%	250 mg
		1 g
		5 g
52,085-3	Engelhard code S3020	1 g
		Pt ≥46.2%
		5 g
		25 g

Sodium hydride

22,344-1	CAS No. 7646-69-7	50 g
		NaH FW 24.00
		250 g

dry, 95% 1 kg
 Direct intercalation into C₆₀ results in the superconducting material (NaH)₄C₆₀.
 mp 800 °C (dec.) (lit.)
References
 Imaeda, K. et al., *Solid State Commun.* **99**, 479 (1996)
 R: 15 S: 24/25-43-7/8

For anhydrous metal halides See: Anhydrous Halides Page 255

For Molecular Based Superconductors see Charge Transfer Complexes Section See: Charge Transfer Complexes Page 182