

Solvent Products

Solvent Packaging

Solvent Delivery
Systems



Solvents Application Guide

Table of Contents

Purity Chart.....	4	GC Resolv.....	13	Electronic Grade.....	24
Solvent Selection Guide.....	5	HPLC.....	13-15	Histological.....	24-25
Innovative Packaging.....	6-7	Certified ACS Plus.....	16	Laboratory Grade.....	26-27
High Volume Solvent		Certified ACS.....	16-20	Technical Grade.....	27
Delivery Systems.....	8-9	Certified.....	20-22	Reagent Grade.....	28
Solvents by Grade.....	10-30	Pesticide Grade.....	22-23	Multicompendial.....	28-29
Optima LC/MS.....	10-11	Spectranalyzed.....	23-24	Special Solvent Grades.....	29-30
Optima.....	12	Scintanalyzed.....	24	Solvents by Alphabetical Order ...	31-42



- 1 The bar code is an alphanumeric code widely used by industry for inventory and tracking control. This "three-of-nine" code meets the Health Industry Bar Code (HIBC) Supplier Labeling Standard as prepared by the Health Industry Business Communication Council (HIBCC).
- 2 This area contains the appropriate precautionary signal word, i.e., DANGER, WARNING, CAUTION, and highlights the potential health hazards, both acute and chronic, of the product.
- 3 In this area, the molecular formula and formula weight are indicated along with the product Chemical Abstract Service (CAS) Registry Number. The Date Received is blank so that the customer can write in the date the bottle was received to assist in inventory control.
- 4 The Maximum Limits of Impurities are indicated.
- 5 The full Product Name and Purity Grade are listed.
- 6 **Fisher Chemical® ChemAlert® Guide** A three-part guide that provides the Safety Code, the NFPA Code and a Storage Code.
- 7 **NFPA Code:** The National Fire Protection Association (NFPA) ratings provide information relating to the severity of health, flammability, reactivity, and related hazards that may be presented by short-term, acute exposure to material during handling under emergency conditions such as spills, leaks, and fires. Each section displays a numerical severity rating from four (most severe) to zero (least severe). The fourth section, as stipulated by NFPA, is left blank and is reserved for indicating any unusual reactivity with water or oxidizing properties. (For more information, see the Technical References section.)
- 8 **Storage Code:** This color-coded bar provides an instant guide to storage. The storage code is also spelled out for additional clarification. The five storage colors and their descriptions are listed on p. 3.
- 9 The U.S. Department of Transportation (DOT) shipping information includes the DOT Proper Shipping Name and the UN/NA Number identifies the United Nations/North American numerical designation for transportation hazards.
- 10 The catalog number is shown for identification and reference.
- 11 The address and telephone number of the manufacturer, Fisher Chemical, is listed.
- 12 **Package Size:** The amount of material per unit or case.
- 13 Lot Number

NOTE: ChemAlert is an instant guide only. It should be supplemented by reading the rest of the label (which provides detailed instructions in the event of accidental exposure, spill, or fire, and applicable OSHA, DOT, and ANSI® data) and also by the appropriate Material Safety Data Sheet (MSDS) and standard references.

Acetonitrile		HPLC		
Also meet 6 ACS specifications 7				
F	Quantity	Packaging	Cat. No.	Each Case of
	1L	Amber Glass	A990-1	93.24 6/397.15
	1L	Amber Glass/Safe-Cote*	A990SX-1	100.47 6/453.48
	2.5L	Amber Glass	A990-212	191.24 4/680.73
	4L	Amber Glass	A990-4	228.23 4/648.19
	4L	Amber Glass/Safe-Cote*	A990SX-4	240.74 4/683.67
	19L	NOWPak* 1 (Plastic Exterior)	A998N1-19	773.72
Also available in recycled FisherPak* and NOWPak* containers.				
Methyl Cyanide, Ethanenitrile				
Melting point: -94.3°C				
CH ₃ CN; C ₂ H ₃ N; F.W. 41.05				
Submicron Filtered, [75-05-8]				
Product Specifications				
Actual Lot Analysis is reported on label.				
	Assay (CH ₃ CN)			>99.0%
	Color (APHA)			≤10
Optical Absorbance:				
	at 190nm			≤1.00
	at 195nm			≤0.15
	at 200nm			≤0.07
	at 205nm			≤0.05
	at 210nm			≤0.04
	at 220nm			≤0.02
	at 254nm			≤0.01
Fluorescence Background (as Quinine Sulfate) To pass test				
LC Gradient Suitability To pass test				
	Refractive Index at 25°C			1.3405-1.3425
	Residue after Evaporation			≤1ppm
	Water			≤0.01%
	Titration Acid			≤0.008mEq/g
	Titration Base			≤0.006mEq/g
	Density at 25°C			0.775-0.780g/mL

- Product name
- Company logo
- Container size
- Synonym(s), Formula, Formula Weight, CAS Registry Number. Also: physical data; special application notes.
- Product specifications
- Type of primary container
- Catalog number
- Purity grade. Grades of FisherChemicals are described on next page.
- Price
- Storage code color. ChemAlert* storage codes are described at right.
- Spill cleanup kit symbol. See right for more information.

ACCEPTED NOMENCLATURE

- Chemical salts and compounds are listed under element name: "Sodium Benzoate," not "Benzoate of Sodium."
- Prefixes for organic compounds—such as alpha, iso, meta, secondary, and abbreviations thereof—do not affect alphabetical order. (Exceptions: Cases where prefix is considered part of the name.)
- Numerical prefixes, such as tri and tetra, are treated as an integral part of the name. The prefix mono is not used.

PACKAGING FOR SAFETY, CONVENIENCE, AND PRODUCT QUALITY

FisherChemical* solvents come in a wide variety of innovative packaging designed for safety, environmental protection, convenient handling and storage, and preservation of product integrity. Packaging also meets current government regulations. Shipping cartons covered by the U.S. Department of Transportation (DOT) regulations also meet international hazardous materials regulations and standards. The primary container is included in the product description of most chemicals in this guide.

Primary containers include:

- Plastic and glass bottles, jars
- Tin can
- SafeCote* PVC-coated bottles
- SafeTin* cans
- Aluminum cans
- Square poly bottles
- Steel Drum/Pail
- PolyPac* containers
- Fisher Pak

CHEM ALERT STORAGE CODES PROVIDE GUIDELINES



RED (R): Flammable. Store in an area segregated for flammable reagents.



BLUE (B): Health hazard. Toxic if inhaled, ingested, or absorbed through skin. Store in secure area.



YELLOW (Y): Reactive and oxidizing reagents. May react violently with air, water, or other substances. Store away from flammable and combustible materials.



WHITE (W): Corrosive. May harm skin, eyes, mucous membranes. Store away from red-, yellow-, and blue-coded reagents.



GRAY (G): Presents no more than moderate hazard in any of the categories above. For general chemical storage.

EXCEPTION: Reagent incompatible with other reagents of the same color bar. Store separately.

SPILL CLEANUP KIT SYMBOLS

To simplify ordering the appropriate spill cleanup kits, the following symbols have been used to identify chemicals that require special cleanup materials and the type of cleanup kit required. These are one-time-use kits containing everything necessary for spill cleanup.



Mercury Warning. This product contains Mercury.



Acid Spill Emergency Cleanup Kit for neutralizing and absorbing up to 1 liter of acid. **Cat. No. 18-061A.**



Caustic Spill Emergency Cleanup Kit for neutralizing and absorbing up to 1 liter of caustic. **Cat. No. 18-061C.**



Solvent Spill Emergency Cleanup Kit for absorbing up to 4 liters of solvent. **Cat. No. 18-061B.**



Styrofoam®-free EcoSafPak™ packaging. We now deliver all 4 x 4L, 1 x 4L, 6 x 1L, 6 x 500mL, 1 x 2.5L, 1 x 1L, 1 x 500mL, 2 x 4L, 2 x 1L and 2 x 500mL glass bottles (including the Safe-Cote™ glass bottles) of your favorite products in eco-friendly EcoSafPak packaging manufactured by an SFI certified manufacturer.



FIBER USED IN THIS PRODUCT LINE MEETS THE SOURCING REQUIREMENTS OF THE SFI PROGRAM
WWW.SFIPROGRAM.ORG



EcoSafPak packaging minimizes the use of non-recyclable material through the use of the revolutionary Hexacomb design inserts.

The Hexacomb is:

- Fully Recyclable
- Sturdy and Durable
- Versatile

Improved safety & handling

- EcoSafPak is the only package of its kind to pass the most demanding ISTA (International Safe Transport Association) "3A" test, consisting of:
 - 17 Drops to simulate real-world parcel shipment handling
 - Shock testing
 - Vibration testing
- Corrugated material is shock absorbing
- Greater stability in the frame of the box
- Staggered handholds for ease of transport

The EcoSafPak is now manufactured by an SFI-certified manufacturer.

The Sustainable Forestry Initiative® (SFI) program:

- is based on the premise that responsible environmental behavior and sound business decisions can co-exist
- is a comprehensive system of principles, objectives, and performance measures developed by professional foresters, conservationists, and scientists
- promotes the perpetual growing and harvesting of trees with the long-term protection of wildlife, plants, soil, and water

SFI's Nine Objectives:

- Sustainable Forestry
- Responsible Practices
- Reforestation and Productive Capacity
- Forest Health and Productivity
- Long-Term Forest and Soil Productivity
- Protection of Water Resources
- Protection of Special Sites and Biological Diversity
- Legal Compliance
- Continual Improvement

Fisher Chemical Products Purity Grades

GRADE	DEFINITION	APPLICATION	CERTIFICATE OF ANALYSIS
OPTIMA* LC/MS	Solvents and additives of exceptionally high purity specially designed and use-tested on LC/MS instruments. Solvents impurity levels in ppb.	LC/MS, HPLC, GC and spectrophotometry	Provided with each shipment
OPTIMA*	Acids and solvents of extremely high purity. Acids are analyzed for 65 metals by ICP/MS; impurity levels in ppt. Solvent impurity levels in ppm. UV absorbance curves and sample chromatograms available on request. For OPTIMA acids, a typical lot analysis is given in the catalog. Impurity levels in an actual lot received may vary from the amount listed.	HPLC, GC, plasma/ICP, spectrophotometry, and pesticide residue analysis	Provided with each shipment
GC <i>Resolv</i> *	Solvents with the highest purity and lot-to-lot consistency. Free of contaminants to the ppb level, including those listed in Contract Laboratory Program Target Compound List. Meet ACS specifications. Chromatogram available on request.	Gas Chromatography (GC)	Provided with each shipment
HPLC	Solvents manufactured specifically for use with HPLC instruments. Meet all ACS specifications. Submicron filtered.	HPLC and spectrophotometry procedures	Available on request
Certified ACS Plus	Acids that, in addition to meeting or exceeding the latest specifications of the ACS, are analyzed for more than 16 metals. Actual lot analysis on label.	Analytical applications with tighter metal specifications	Available on request
Certified ACS	Reagent chemicals that meet or exceed the latest ACS Specifications. Actual lot analysis on label.	Analytical applications requiring tight specifications	Available on request
Certified	Reagent chemicals for which the purity standard is established by Fisher Scientific. Purity is guaranteed to meet published maximum limits of impurities.	General analytical procedures	Available on request
Plasma Grade	Solvents manufactured for use with plasma/ICP instruments. Impurity levels in ppt. Packaged in acid-cleaned polyethylene bottles.	Plasma/ICP, environmental testing, trace-metal analysis	Available on request
Environmental Grade	Solvents for use in HPLC analysis, trace-organic analysis, and environmental testing. Packaged in pre-cleaned glass bottles in M3.5 (Class 100) cleanroom. Shipped with Certificate of Analysis.	HPLC, trace-organic analysis, environmental testing	Provided with each shipment
Pesticide	Solvents for use in analysis of pesticide residue. Meet or exceed ACS standards of purity for pesticide residue analysis.	GC with electron capture detector (ECD), pesticide residue analysis	Available on request
USP/NF/FCC/EP/BP/JP	Reagent chemicals that meet or surpass specifications of the United States Pharmacopeia (USP), the National Formulary (NF), the Food Chemicals Codex (FCC), the European Pharmacopeia (EP), the British Pharmacopeia (BP), and/or the Japanese Pharmacopeia (JP).	Food and drug laboratories, biological testing	20L or greater; 10kg or greater
Spectranalyzed*	Solvents for use in spectrophotometry. Also meet ACS specifications. Actual lot analysis on label.	Ultraviolet and visible wavelength detectors (UV-Vis)	Available on request
Biotechnology	Solvents and reagents that have been specially purified and assayed for biotechnology applications.	Electrophoresis, molecular biology, sequencing, and peptide and oligonucleotide synthesis	Available on request
Scintanalyzed*	Solvents, fluors, and prepared cocktails for liquid scintillation counting. Includes nonflammable, nontoxic, biodegradable ScintiSafe* cocktails.	Liquid scintillation counting	Available on request
Electronic	Solvents manufactured to ensure low levels of metal contamination. Meet Semiconductor Equipment and Materials Institute (SEMI) requirements. Actual lot analysis on label.	Electronics and circuit board manufacturing	Available on request
TraceMetal	Acids manufactured to achieve low metal contamination measurable in ppm to ppb range. Each lot is analyzed for more than 55 metals by ICP/MS. For TraceMetal acids, a typical lot analysis is given in the catalog. Impurity levels in an actual lot received may vary from amounts listed.	Primarily used in digestion of samples prior to instrument (ICP) analysis	Provided with each shipment
Histology	Solvents and products that are specially prepared for use in the histology laboratory setting. Solvents are filtered for tissue processing applications.	Tissue processing, clinical or histology procedures	Available on request
Laboratory and Technical	Chemicals of reasonable quality and purity for use in procedures where no official standards are required.	Manufacturing and general laboratory use	Available on request

Fisher Chemical High Purity Solvents Selection Guide

Solvent	Available Grades*	UV Cutoff (nm)	Boiling Point (°C)	Density (g/mL, 25°C)	Refractive Index (25°C)	Melting Point (°C)	Polarity Index (P')	Elutropic Value on Silica (ε°)	Viscosity (cP, 20°C)	Flash Point (°C)	Mol. Wt.
Acetone	GOHPSEF	330	56.1	0.7857	1.3568	-94.3	5.1	0.53	0.36	20	58.08
Acetonitrile	OHP	190	81.6	0.7780	1.3415	-50.0	5.8	0.52	0.36	2	41.05
1-Butanol	HS	215	117.7	0.8098	1.3972	-88.6	3.9	-	2.98	35	74.12
Chloroform	OMHPS	245	61.7	1.4840	1.4445	-63.3	4.1	0.26	0.58	none	119.38
Cyclohexane	HPS	202	80.7	0.7740	1.4247	-6.5	0.2	0.03	0.90	-20	84.16
N,N-Dimethylformamide	S	268	153.0	0.9440	1.4280	-61.0	6.4	-	0.92	58	73.09
Dimethyl Sulfoxide	SH	262	189.0	1.1014	1.4783	18.5	7.2	-	2.24	87.8	78.13
Ethyl Acetate	OHPSPh	255	77.1	0.8940	1.3695	-83.9	4.4	0.38	0.45	-4	88.11
Ethyl Ether	HPS	218	34.6	0.7134	1.3500	-116.3	2.8	0.43	0.24	-45	74.12
Glycerol	MSPH	205	290.0	1.2613	1.4746	18.2	-	-		193	92.09
Heptane	HSO	197	98.4	0.6838	1.3855	-90.6	0.2	0.01	0.40	-4	100.20
Hexanes	GOHPSPH	195	69.0	0.6630	1.3759	-95.3	0.1	0.01	0.31	-23	86.18
Isooctane	OHPS	205	99.2	0.6919	1.3895	109.5	0.1	0.01	0.50	28	114.23
Methanol	GPtOHPSLw	205	64.7	0.7915	1.3288	-97.8	5.1	0.73	0.55	12	32.04
Methylene Chloride	GOHPS	233	39.5	1.3180	1.4215	-96.7	3.1	0.32	0.30	N/A	84.93
N-Methylpyrrolidinone	M	275	202.2	1.03	1.469	-24.4	-	-	1.67	95	99.13
Pentane	HPS	190	36.1	0.6264	1.3555	-129.7	0.0	0.00	0.22	-49	72.15
Petroleum Ether	OP	-	35-60	0.6400	1.3610	-	0.1	-	-	-18	-
2-Propanol	OHPSLwPh	205	82.3	0.7855	1.3772	-90.0	3.9	0.63	2.40	-12	60.10
Tetrahydrofuran	OHS	210	66.1	0.8892	1.4060	-108.3	4.0	0.35	0.55	-14	72.11
Toluene	OHSSc	285	110.6	0.8660	1.4940	-95.0	2.4	0.22	0.59	-4	92.14
Water	OH	-	100.0	0.9982	1.3330	0.0	10.2	-	1.00	N/A	18.02

*G = GC *Resolv*[®] Pt = Purge & Trap O = Optima[®] M = Molecular Biology Grade H = HPLC Grade P = Pesticide Grade S = Spectranalyzed[®] E = Electronic Grade
Sc = Scintanalyzed[®] Lw = Low water

Ph = Pharmacopeia, one or more of the following: USP, NF, FCC, EP, BP, JP

Excellence in Packaging through Innovation

Fisher Chemical products come in a variety of innovative packaging options designed for safety, environmental protection, convenient handling and storage, and preservation of product integrity. Our packaging is compliant with all government regulations.



Plastic Bottles—Safety and Convenience

Available when chemical properties are compatible to minimize the risk of breakage, provide lighter weight packaging and easier, more economical shipping.

Amber Glass Bottles—Quality and Reliability

Used to package photosensitive chemicals, to protect them from light.

Safe-Cote® PVC-Coated Bottles—Safety

These tough bottles provide the purity of glass and most of the benefits of plastic for the storing and dispensing of solvents and acids. If the bottle breaks, glass fragments and liquids are more likely to remain trapped.

Aluminum Cans—Quality and Reliability

Designed to contain ethers, Fisher Chemical 1L and 4L aluminum cans feature a round, seamless construction, reducing the possibility of leakage through a seam.



Lock in Quality and Safety with Fisher Chemical Products

Fisher Chemical products are now packaged with the innovative, new FisherLOCK™ Cap, designed to help maintain the highest standards in product quality and chemical safety practices.

With an exclusive color-coded design (patent pending), the FisherLOCK™ Cap is engineered to lock to the transfer bead of the bottle when it is applied during manufacturing to a filled bottle of laboratory chemicals. The cap design includes an interior ring, visible from various angles, that offers resistance until it separates from the cap when the bottle is first opened.

The presence and location of that ring imparts tamper evidence to the configuration—without the drawbacks and complications of a plastic seal over the cap. The FisherLOCK™ Cap is easy to use, and supports product quality, reliability of use, and safety in use of the chemical products you purchase.



The FisherLOCK™ Cap LOCKS in quality, safety, reliability, and convenience:

QUALITY

- Provides a tight, tamper-evident, secure seal to assure chemical contents arrive fresh and unopened
- Eliminates polyethylene glycol contamination that is possible with a plastic overseal

SAFETY

- Caps are designed to resist back-off during transport, reducing risk of leakage
- Color-coded rings indicate storage requirements and hazard categories, and enhance proper recognition, handling and storage, even before the bottle is removed from the case

RELIABILITY

- Rigorously tested for chemical compatibility
- Bottle threads are unchanged, allowing you to continue to attach the opened bottle to standard equipment

CONVENIENCE

- Cap design facilitates correct initial torque application during manufacturing, thus eliminating caps that may be hard to twist open
- Larger ridges on the exterior of the cap make it easier to open
- Caps readily reseal after initial opening



High Volume Solvent Delivery Systems

The eco-friendly option for your laboratory

High Volume Solvent Systems offer environmentally friendly solvent handling solutions for virtually unlimited applications, enhancing safety and improving productivity within your lab.

Safety and Convenience

High Volume Solvent Delivery Systems increase lab efficiency while reducing lab-operating costs by eliminating:

- Repeated solvent testing
- Multiple lots of material
- Bottle rinsing
- Disposal costs

Environmentally Friendly

- Reduce the amount of solid waste generated in your laboratory (A single reusable 200 liter FisherPak® container replaces approximately 50 four-liter bottles and packing material)
- Minimize the release of flammable or toxic solvent liquids and vapors
- Eliminate bottle rinsing—Fisher Scientific cleans and refills empty containers

Enhanced Solvent Safety

High Volume Solvent Delivery Systems incorporate safety features to protect the lab and the environment by offering a unique combination of mechanical and manual controls to prevent unwanted solvent flow. The bottle-free, closed system eliminates the potential for glass bottle breakage and makes the risk of spills and exposure to vapors negligible.

Virtually Unlimited Applications

- High-performance liquid chromatography (HPLC)
- Sample preparation for spectroscopic analysis
- Process synthesis, extractions, preparative chromatography, and high-volume gas chromatography sample preparation

High Volume Solvent Delivery Systems from Fisher Scientific

FisherPak

The Fisher Chemical FisherPak is a refillable dispensing system that protects solvent purity and reduces solvent spills. The reusable 316 stainless-steel container is pressurized with an inert gas to deliver the solvent, so the system remains closed and the solvent remains pure.

Key Features:

- Custom solvents blended and/or tested to your specific requirements
- Customizable to your needs with our extensive list of accessories
- ASME3 approved containers available for operating pressures over 15psi
- Integrated fork channels are available for larger sizes
- Factory Mutual (FM) approved

Available Sizes: 19L, 28L, 50L, 115L, 200L, 1350L



BasicPak

The Fisher Chemical BasicPak is a returnable, reusable drum made from high-grade 304 stainless steel. This system is designed to deliver inert gas safely to the drum with 7-15psig without contamination.

Key Features:

- Equipped with Mico-Matic coupling technology
- Safety pin on dust cap provides tamper evidence
- Heavy-duty fork channels made of 316 stainless steel are available

Available Sizes: 19L, 50L, 115L, 200L



DelPak

DelPak combines the strength and durability of steel with the chemical compatibility of polyethylene.

Key Features:

- Constructed in FDA-approved high density blow molded polyethylene liner
- Steel overpack (pail)
- Inner liner bottle has a two-inch buttress plug with a three-fourth inch NPT center reducer plug and handle for easy removal
- Nine-gauge wire handle with plastic grip for carrying

Available Size: 18L



NowPak I

The inner liner of NowPak I is constructed with an inert fluorocarbon polymer that is compatible with all high-purity solvents. The inner liner collapses around the dip tube during dispensing, which prevents air from entering the system. The sealed dispensing options also eliminate introduction of lab impurities into the solvent.



Key Features:

- Six different dispensing options to meet a variety of needs
- Liners are pre-cleaned in a class 10/100 cleanroom environment to provide ultimate purity
- Liner materials have no pigments, adhesives, additives or stabilizers
- Double-containment design enhances safety and meets OSHA guidelines
- HDPE overpack includes a UV-blocking additive
- Overpack can be reused or recycled

Available Size: 19L

NowPak II

The NowPak II precleaned liner technology is offered in a pressurizable, stainless-steel overpack. The system permits ultra-clean chemical dispensing without direct contact of the chemical with the drive gas. Drive gas pressure is applied to the outside of the liner, minimizing gas entrainment or micro-bubble formation. As the fill material is dispensed, the inner liner collapses. After emptying the system, the liner is simply removed and the overpack is ready for insertion of a new, precleaned liner. The overpack, closer, and dip tube are designed for reuse.



Key Features:

- Reusable stainless-steel (304) overpack with 15psi maximum operating pressure
- Pressure or pump dispense connector options
- Fluoropolymer or co-extruded polyethylene liner options
- Breakseal or thread on/off closure options

Available Sizes: 10L, 19L

Optima® LC/MS Solvents



Acetonitrile, Methanol, 2-Propanol (IPA) and Water

- Lower contamination from plasticizer peaks
- Exceptionally low metal ion content
- Fewer background peaks
- Higher Signal intensity
- Low LC/UV response

The joining of liquid chromatography (LC) with mass spectrometry (MS) has become an indispensable tool for various fields of research. The value of LC/MS derives from its ability to combine separation chemistry with selective mass ion detection. As instrumentation advances lead to ever-lower analyte detection limits, it is crucial for the chromatographer to consider the level of purity when selecting appropriate solvents for use in the LC/MS mobile phase. The Fisher Chemical product line offers superior high-purity solvents designed to meet the required purity level of advanced LC/MS systems.

For more information, including a white paper titled 'Optimizing Mobile Phase Solvent Purity for LC/MS', go to www.FisherLCMS.com.

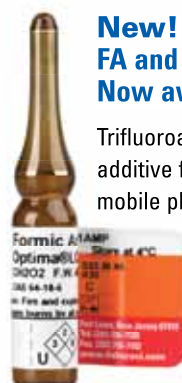
New! FA and TFA Mobile Phase Blends, OPTIMA LC/MS Grade

High purity solvents are pre-blended with modifiers such as formic acid (FA) or trifluoroacetic acid (TFA) to provide ready-to-use mobile phase for LC/MS and LC/UV applications.



New! FA and TFA, OPTIMA LC/MS Grade - Now available in ampules packaging

Trifluoroacetic Acid is an ultrapure reagent used as an additive for the formulation of solvent blends for the mobile phase in LC/MS applications.



Optima LC/MS Solvents

Optima LC/MS Solvents are manufactured and tested specifically to meet the stringent purity requirements of advanced Optima LC/MS systems.

- Lower contamination from plasticizer peaks
- Exceptionally low metal ion content
- Fewer background peaks
- Higher signal intensity
- Low LC/UV response



Optima LC/MS

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetonitrile	Optima LC/MS	500mL	Amber Glass	A955-500
Acetonitrile	Optima LC/MS	1L	Amber Glass	A955-1
Acetonitrile	Optima LC/MS	2.5L	Amber Glass	A955-212
Acetonitrile	Optima LC/MS	4L	Amber Glass	A955-4
Methanol	Optima LC/MS	500mL	Amber Glass	A456-500
Methanol	Optima LC/MS	1L	Amber Glass	A456-1
Methanol	Optima LC/MS	2.5L	Amber Glass	A456-212
Methanol	Optima LC/MS	4L	Amber Glass	A456-4
2-Propanol	Optima LC/MS	500mL	Amber Glass	A461-500
2-Propanol	Optima LC/MS	1L	Amber Glass	A461-1
2-Propanol	Optima LC/MS	2.5L	Amber Glass	A461-212
2-Propanol	Optima LC/MS	4L	Amber Glass	A461-4
Water	Optima LC/MS	500mL	Amber Glass	W6-500
Water	Optima LC/MS	1L	Amber Glass	W6-1
Water	Optima LC/MS	2.5L	Amber Glass	W6-212
Water	Optima LC/MS	4L	Amber Glass	W6-4

Solvent Blends

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
0.1% FA in Water	Optima LC/MS	500mL	Amber Glass	LS118-500
0.1% FA in Water	Optima LC/MS	1L	Amber Glass	LS118-1
0.1% FA in Water	Optima LC/MS	2.5L	Amber Glass	LS118-212
0.1% FA in Water	Optima LC/MS	4L	Amber Glass	LS118-4
0.1% TFA in Water	Optima LC/MS	500mL	Amber Glass	LS119-500
0.1% TFA in Water	Optima LC/MS	1L	Amber Glass	LS119-1
0.1% TFA in Water	Optima LC/MS	2.5L	Amber Glass	LS119-212
0.1% TFA in Water	Optima LC/MS	4L	Amber Glass	LS119-4
0.1% FA in Acetonitrile	Optima LC/MS	500mL	Amber Glass	LS120-500
0.1% FA in Acetonitrile	Optima LC/MS	1L	Amber Glass	LS120-1
0.1% FA in Acetonitrile	Optima LC/MS	2.5L	Amber Glass	LS120-212
0.1% FA in Acetonitrile	Optima LC/MS	4L	Amber Glass	LS120-4
0.1% TFA in Acetonitrile	Optima LC/MS	500mL	Amber Glass	LS121-500
0.1% TFA in Acetonitrile	Optima LC/MS	1L	Amber Glass	LS121-1
0.1% TFA in Acetonitrile	Optima LC/MS	2.5L	Amber Glass	LS121-212
0.1% TFA in Acetonitrile	Optima LC/MS	4L	Amber Glass	LS121-4

Solvent Additives

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Formic Acid	Optima LC/MS	50mL	Poly Bottle	A117-50
Formic Acid	Optima LC/MS	10 x 1mL	Ampule	A117-10X1AMP
Formic Acid	Optima LC/MS	1mL	Ampule	A117-1AMP
Formic Acid	Optima LC/MS	0.5mL	Ampule	A117-05AMP
Formic Acid	Optima LC/MS	2mL	Ampule	A117-2AMP
Trifluoroacetic Acid	Optima LC/MS	50mL	Poly Bottle	A116-50
Trifluoroacetic Acid	Optima LC/MS	10 x 1mL	Ampule	A116-10X1AMP
Trifluoroacetic Acid	Optima LC/MS	1mL	Ampule	A116-1AMP
Trifluoroacetic Acid	Optima LC/MS	0.5mL	Ampule	A116-05AMP
Trifluoroacetic Acid	Optima LC/MS	2mL	Ampule	A116-2AMP

Optima Solvents

Optima Solvents meet extremely high purity (ppm) levels. They are manufactured for use when contaminant-free performance is essential—HPLC, GC, plasma/ICP, spectrophotometry, environmental testing, and other analytical applications.

- Contaminant-free to ppb and ppm levels, depending on solvent
- Certificate of Analysis provided with each shipment
- Supplied in specially cleaned bottles
- Blanketed with inert gas to maintain purity
- Chromatograms available on request

Optima

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetone	Optima	1L	Amber Glass	A929-1
Acetone	Optima	4L	Amber Glass	A929-4
Acetone	Optima	4L	Glass/Safe-Cote	A929SK-4
Acetone	Optima	28L	FisherPak	A929RS28
Acetone	Optima	115L	FisherPak	A929RS115
Acetonitrile	Optima	1L	Amber Glass	A996-1
Acetonitrile	Optima	4L	Amber Glass	A996-4
Acetonitrile	Optima	4L	Amber Glass/Safe-Cote	A996SK4
Acetonitrile	Optima	19L	FisherPak	A996N219
Chloroform (Approx. 50ppm Amylene as Preservative)	Optima	4L	Amber Glass	C297-4
Ethyl Acetate	Optima	4L	Amber Glass	E196-4
Ethyl Acetate	Optima	4L	Safe-Cote	E196SK-4
Hexanes	Optima	1L	Amber Glass	H303-1
Hexanes	Optima	4L	Amber Glass	H303-4
Hexanes	Optima	4L	Amber Glass/Safe-Cote	H303SK-4
Hexanes	Optima	19L	FisherPak	H303SS19
n-Hexane, 95%	Optima	1L	Amber Glass	H306-1
n-Hexane, 95%	Optima	4L	Amber Glass	H306-4
n-Hexane, 95%	Optima	4L	Amber Glass/Safe-Cote	H306SK-4
Isooctane	Optima	4L	Amber Glass	O301-4
Methanol	Optima	1L	Amber Glass	A454-1
Methanol	Optima	4L	Amber Glass	A454-4
Methanol	Optima	4L	Amber Glass/Safe-Cote	A454SK-4
Methanol	Optima	115L	FisherPak	A454SS-115
Methanol	Optima	200L	FisherPak	A454SS-200
Methylene Chloride	Optima	1L	Amber Glass	D151-1
Methylene Chloride	Optima	4L	Amber Glass	D151-4
Methylene Chloride	Optima	4L	Amber Glass/Safe-Cote	D151SK-4
Methylene Chloride	Optima	19L	FisherPak	D151SS19
Petroleum Ether	Optima	4L	Amber Glass	E120-4
Petroleum Ether	Optima	4L	Amber Glass/Safe-Cote	E120SK-4
Petroleum Ether	Optima	28L	FisherPak	E120SS-28
Petroleum Ether	Optima	200L	FisherPak	E120SS-200
2-Propanol	Optima	4L	Amber Glass	A464-4
2-Propanol	Optima	4L	Amber Glass/Safe-Cote	A464SK-4
Tetrahydrofuran	Optima	1L	Amber Glass	T427-1
Tetrahydrofuran	Optima	4L	Amber Glass	T427-4
Tetrahydrofuran	Optima	4L	Amber Glass/Safe-Cote	T427SK-4
Toluene	Optima	4L	Amber Glass	T291-4
Toluene	Optima	4L	Safe-Cote/Amber Glass	T291SK-4
Toluene	Optima	19L	FisherPak	T291SS19
Toluene	Optima	200L	FisherPak	T291RS200
Water	Optima	4L	Amber Glass	W7-4
Water	Optima	4L	Safe-Cote/Amber Glass	W7SK-4

GC Resolv Solvents

GC Resolv Solvents exhibit the very highest purity and lot-to-lot consistency for gas chromatography applications.

- Contaminant-free to ppb levels
- Meet ACS specifications
- Certificate of Analysis provided with each shipment
- Supplied in specially cleaned bottles
- Blanketed with inert gas to maintain purity
- Chromatograms available on request

GC Resolv

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetone	GC Resolv	4L	Amber Glass	A928-4
n-Hexane	GC Resolv	4L	Amber Glass	H307-4
Methanol	GC Resolv	4L	Amber Glass	A457-4
Methylene Chloride	GC Resolv	4L	Amber Glass	D154-4
Methanol (Low Water)	GC Resolv	200L	BasicPak	A935RB200
2-Propanol	GC Resolv	200L	FisherPak	A464RS200

HPLC Grade Solvents

HPLC Grade Solvents are manufactured especially for use with HPLC instruments.

- Meet ACS specifications
- Submicron filtered
- Actual lot analysis on label
- Supplied in specially cleaned bottles
- Blanketed with inert gas to maintain purity
- Certificate of Analysis available on request

HPLC

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetone	HPLC	1L	Amber Glass	A949-1
Acetone	HPLC	1L	Amber Glass/Safe-Cote	A949SK-1
Acetone	HPLC	4L	Amber Glass	A949-4
Acetone	HPLC	4L	Amber Glass/Safe-Cote	A949SK-4
Acetone	HPLC	19L	NOWPak I	A949N1-19
Acetone	HPLC	115L	FisherPak	A949SS-115
Acetonitrile	HPLC	1L	Amber Glass	A998-1
Acetonitrile	HPLC	1L	Amber Glass/Safe-Cote	A998SK-1
Acetonitrile	HPLC	2.5L	Amber Glass	A998212
Acetonitrile	HPLC	4L	Amber Glass	A998-4
Acetonitrile	HPLC	4L	Amber Glass/Safe-Cote	A998SK-4
Acetonitrile	HPLC	19L	NOWPak I	A998N1-19
Alcohol	HPLC	4L	Amber Glass	A995-4
Alcohol	HPLC	200L	FisherPak	A995RS200
1-Butanol	HPLC	1L	Amber Glass	A383-1
1-Butanol	HPLC	1L	Amber Glass/Safe-Cote	A383SK-1
1-Butanol	HPLC	4L	Amber Glass	A383-4
1-Butanol	HPLC	4L	Amber Glass/Safe-Cote	A383SK-4

HPLC (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
n-Butyl Chloride	HPLC	4L	Amber Glass	B429-4
n-Butyl Chloride	HPLC	4L	Amber Glass/Safe-Cote	B429SK-4
Chloroform (Approx. 50ppm Pentene as Preservative)	HPLC	1L	Amber Glass	C607-1
Chloroform (Approx. 50ppm Pentene as Preservative)	HPLC	1L	Amber Glass/Safe-Cote	C607SK-1
Chloroform (Approx. 50ppm Pentene as Preservative)	HPLC	4L	Amber Glass	C607-4
Chloroform (Approx. 50ppm Pentene as Preservative)	HPLC	4L	Amber Glass/Safe-Cote	C607SK-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	HPLC	1L	Amber Glass	C606-1
Chloroform (Approx. 0.75% Ethanol as Preservative)	HPLC	1L	Amber Glass/Safe-Cote	C606SK-1
Chloroform (Approx. 0.75% Ethanol as Preservative)	HPLC	4L	Amber Glass	C606-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	HPLC	4L	Amber Glass/Safe-Cote	C606SK-4
Cyclohexane	HPLC	1L	Amber Glass	C620-1
Cyclohexane	HPLC	1L	Amber Glass/Safe-Cote	C620SK-1
Cyclohexane	HPLC	4L	Amber Glass	C620-4
Cyclohexane	HPLC	4L	Amber Glass/Safe-Cote	C620SK-4
Dimethyl Sulfoxide	HPLC	4L	Amber Glass	D159-4
Ethyl Acetate	HPLC	1L	Amber Glass	E195-1
Ethyl Acetate	HPLC	1L	Amber Glass/Safe-Cote	E195SK-1
Ethyl Acetate	HPLC	4L	Amber Glass	E195-4
Ethyl Acetate	HPLC	4L	Amber Glass/Safe-Cote	E195SK-4
Ethyl Acetate	HPLC	19L	NOWPak I	E195N1-19
Ethyl Ether Anhydrous (Stabilized)	HPLC	4L	Tin Can	E198-4
Ethyl Ether Anhydrous (Stabilized)	HPLC	19L	FisherPak	E198SS19
Heptane	HPLC	1L	Amber Glass	H350-1
Heptane	HPLC	1L	Amber Glass/Safe-Cote	H350SK-1
Heptane	HPLC	4L	Amber Glass	H350-4
Heptane	HPLC	4L	Amber Glass/Safe-Cote	H350SK-4
Heptane	HPLC	19L	FisherPak	H350RS19
Heptane	HPLC	200L	FisherPak	H350RS200
Hexanes	HPLC	1L	Amber Glass	H302-1
Hexanes	HPLC	1L	Amber Glass/Safe-Cote	H302SK-1
Hexanes	HPLC	4L	Amber Glass	H302-4
Hexanes	HPLC	4L	Amber Glass/Safe-Cote	H302SK-4
Hexanes	HPLC	50L	FisherPak	H302SS-50
Hexanes	HPLC	115L	FisherPak	H302SS-115
Hexanes	HPLC	200L	FisherPak	H302SS-200
Isooctane	HPLC	1L	Amber Glass	O296-1
Isooctane	HPLC	1L	Amber Glass/Safe-Cote	O296SK-1
Isooctane	HPLC	4L	Amber Glass	O296-4
Isooctane	HPLC	4L	Amber Glass/Safe-Cote	O296SK-4
Methanol	HPLC	1L	Amber Glass	A452-1
Methanol	HPLC	1L	Amber Glass/Safe-Cote	A452SK-1
Methanol	HPLC	4L	Amber Glass	A452-4
Methanol	HPLC	4L	Amber Glass/Safe-Cote	A452SK-4
Methanol	HPLC	19L	NOWPak I	A452N1-19

HPLC (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Methyl tert-Butyl Ether	HPLC	4L	Amber Glass	E127-4
Methyl tert-Butyl Ether	HPLC	200L	FisherPak	E127RS200
Methylene Chloride (Not Stabilized)	HPLC	1L	Amber Glass	D150-1
Methylene Chloride (Not Stabilized)	HPLC	1L	Amber Glass/Safe-Cote	D150SK-1
Methylene Chloride (Not Stabilized)	HPLC	4L	Amber Glass	D150-4
Methylene Chloride (Not Stabilized)	HPLC	4L	Amber Glass/Safe-Cote	D150SK-4
Methylene Chloride (Stabilized)	HPLC	1L	Amber Glass	D143-1
Methylene Chloride (Stabilized)	HPLC	1L	Amber Glass/Safe-Cote	D143SK-1
Methylene Chloride (Stabilized)	HPLC	4L	Amber Glass	D143-4
Methylene Chloride (Stabilized)	HPLC	4L	Amber Glass/Safe-Cote	D143SK-4
Methylene Chloride (Stabilized)	HPLC	50L	FisherPak	D143SS-50
Methylene Chloride (With Cyclohexene Preservative)	HPLC	1L	Amber Glass	D138-1
Methylene Chloride (With Cyclohexene Preservative)	HPLC	4L	Amber Glass	D138-4
Methylene Chloride (With Cyclohexene Preservative)	HPLC	4L	Amber Glass/Safe-Cote	D138SK-4
Pentane	HPLC	1L	Amber Glass	P399-1
Pentane	HPLC	1L	Amber Glass/Safe-Cote	P399SK-1
Pentane	HPLC	4L	Amber Glass	P399-4
Pentane	HPLC	4L	Amber Glass/Safe-Cote	P399SK-4
Pentane	HPLC	19L	Steel Drum	P399RS19
2-Propanol	HPLC	1L	Amber Glass	A451-1
2-Propanol	HPLC	1L	Amber Glass/Safe-Cote	A451SK-1
2-Propanol	HPLC	4L	Amber Glass	A451-4
2-Propanol	HPLC	4L	Amber Glass/Safe-Cote	A451SK-4
Tetrahydrofuran	HPLC	1L	Amber Glass	T425-1
Tetrahydrofuran	HPLC	1L	Amber Glass/Safe-Cote	T425SK-1
Tetrahydrofuran	HPLC	4L	Amber Glass	T425-4
Tetrahydrofuran	HPLC	4L	Amber Glass/Safe-Cote	T425SK-4
Tetrahydrofuran	HPLC	50L	FisherPak	T425SS-50
Toluene	HPLC	1L	Amber Glass	T290-1
Toluene	HPLC	1L	Amber Glass/Safe-Cote	T290SK-1
Toluene	HPLC	4L	Amber Glass	T290-4
Toluene	HPLC	4L	Amber Glass/Safe-Cote	T290SK-4
1,2,4-Trichlorobenzene	HPLC	4L	Amber Glass	O4846-4
Triethylamine	HPLC	100mL	Amber Glass	O4884-100
Vinyl acetate	HPLC	115L	BasicPak	O5057FB115
Water	HPLC	1L	Amber Glass	W5-1
Water	HPLC	1L	Amber Glass/Safe-Cote	W5SK-1
Water	HPLC	4L	Amber Glass	W5-4
Water	HPLC	4L	Amber Glass/Safe-Cote	W5SK-4
Water	HPLC	19L	NOWPak I	W5-N119

Certified ACS Plus Solvents

Certified ACS Plus Solvents meet or exceed the latest specifications of the ACS and are suitable for applications with tighter metal specifications.

- Certificate of Analysis available on request

Certified ACS Plus

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
2-Propanol	Certified ACS Plus	500mL	Amber Glass	A416-500
2-Propanol	Certified ACS Plus	1L	Amber Glass	A416-1
2-Propanol	Certified ACS Plus	4L	Amber Glass	A416-4
2-Propanol	Certified ACS Plus	4L	Poly Bottle	A416P-4
2-Propanol	Certified ACS Plus	4L	SafeTin	A416S-4
2-Propanol	Certified ACS Plus	4L	Amber Glass/Safe-Cote	A416SK-4
2-Propanol	Certified ACS Plus	20L	Steel Pail	A416-20
2-Propanol	Certified ACS Plus	200L	Steel Drum	A416-200

Certified ACS Solvents

Certified ACS Solvents meet or exceed the latest ACS specifications and are suitable for analytical applications that require tight specifications.

- Actual lot analysis on label
- Certificate of Analysis available on request

Certified ACS

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetic Anhydride	Certified ACS	100mL	Amber Glass	A10-100
Acetic Anhydride	Certified ACS	500mL	Amber Glass	A10-500
Acetic Anhydride	Certified ACS	1L	Amber Glass	A10-1
Acetic Anhydride	Certified ACS	4L	Amber Glass	A10-4
Acetone	Certified ACS	500mL	Amber Glass	A18-500
Acetone	Certified ACS	1L	Amber Glass	A18-1
Acetone	Certified ACS	4L	Amber Glass	A18-4
Acetone	Certified ACS	4L	Poly Bottle	A18P-4
Acetone	Certified ACS	4L	SafeTin	A18S-4
Acetone	Certified ACS	4L	Amber Glass/Safe-Cote	A18SK-4
Acetone	Certified ACS	20L	Steel Pail	A18-20
Acetone	Certified ACS	200L	Steel Drum	A18-200
Acetone	Certified ACS	350 gal	Steel Tote	A18CU1300
Acetonitrile	Certified ACS	1L	Amber Glass	A21-1
Acetonitrile	Certified ACS	4L	Amber Glass	A21-4
Acetonitrile	Certified ACS	20L	Steel Pail	A21-20
Acetonitrile	Certified ACS	200L	Steel Drum	A21-200
Aniline	Certified ACS	500mL	Amber Glass/Poison Pack	A740I-500
Aniline	Certified ACS	4L	Amber Glass/Poison Pack	A740I-4
n-Amyl Alcohol	Certified ACS	500mL	Amber Glass	A394-500
n-Amyl Alcohol	Certified ACS	4L	Amber Glass	A394-4
1-Butanol	Certified ACS	500mL	Amber Glass	A399-500
1-Butanol	Certified ACS	1L	Amber Glass	A399-1
1-Butanol	Certified ACS	4L	Amber Glass	A399-4

Certified ACS (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
1-Butanol	Certified ACS	4L	SafeTin	A399S-4
1-Butanol	Certified ACS	20L	Steel Pail	A399-20
Carbon Disulfide	Certified ACS	500mL	Amber Glass	C184-500
Carbon Disulfide	Certified ACS	2.5L	Glass	C184-212
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	500mL	Amber Glass	C298-500
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	1L	Amber Glass	C298-1
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	4L	Amber Glass	C298-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	4L	SafeTin	C298S-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	4L	Amber Glass/Safe-Cote	C298SK-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	20L	Steel Pail	C298-20
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	200L	Steel Drum	C298-200
Cyclohexane	Certified ACS	500mL	Amber Glass	C556-500
Cyclohexane	Certified ACS	1L	Amber Glass	C556-1
Cyclohexane	Certified ACS	4L	Amber Glass	C556-4
1,2-Dichloroethane	Certified ACS	500mL	Amber Glass	E175-500
1,2-Dichloroethane	Certified ACS	4L	Amber Glass	E175-4
1,2-Dichloroethane	Certified ACS	19L	FisherPak	E175RS19
1,2-Dichloroethane	Certified ACS	20L	Steel Pail	E175-20
1,2-Dichloroethane	Certified ACS	50L	FisherPak	E175RS50
N,N-Dimethylformamide	Certified ACS	500mL	Amber Glass	D119-500
N,N-Dimethylformamide	Certified ACS	1L	Amber Glass	D119-1
N,N-Dimethylformamide	Certified ACS	4L	Amber Glass	D119-4
N,N-Dimethylformamide	Certified ACS	4L	SafeTin	D119S-4
N,N-Dimethylformamide	Certified ACS	20L	Poly Pail	D119-20
N,N-Dimethylformamide	Certified ACS	200L	Steel Drum	D119200
Dimethyl Sulfoxide	Certified ACS	500mL	Amber Glass	D128-500
Dimethyl Sulfoxide	Certified ACS	1L	Amber Glass	D128-1
Dimethyl Sulfoxide	Certified ACS	4L	Amber Glass	D128-4
Dimethyl Sulfoxide	Certified ACS	50L	FisherPak	D128RS50
1,4-Dioxane	Certified ACS	500mL	Amber Glass	D111-500
1,4-Dioxane	Certified ACS	4L	Amber Glass	D111-4
Ethyl Acetate	Certified ACS	500mL	Amber Glass	E145-500
Ethyl Acetate	Certified ACS	1L	Amber Glass	E145-1
Ethyl Acetate	Certified ACS	4L	Amber Glass	E145-4
Ethyl Acetate	Certified ACS	4L	SafeTin	E145S-4
Ethyl Acetate	Certified ACS	4L	Amber Glass/Safe-Cote	E145SK-4
Ethyl Acetate	Certified ACS	20L	Steel Pail	E145-20
Ethyl Acetate	Certified ACS	200L	Steel Drum	E145-200

Certified ACS (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Ethyl Ether Anhydrous (BHT Stabilized)	Certified ACS	500mL	Aluminum Can	E138500
Ethyl Ether Anhydrous (BHT Stabilized)	Certified ACS	1L	Aluminum Can	E138-1
Ethyl Ether Anhydrous (BHT Stabilized)	Certified ACS	4L	Aluminum Can	E138-4
Ethyl Ether Anhydrous (BHT Stabilized)	Certified ACS	20L	Steel Pail	E138-20
Formaldehyde, 37% by Weight (With Preservative)	Certified ACS	500mL	Amber Glass	F79-500
Formaldehyde, 37% by Weight (With Preservative)	Certified ACS	1L	Amber Glass	F79-1
Formaldehyde, 37% by Weight (With Preservative)	Certified ACS	4L	Amber Glass	F79-4
Formaldehyde, 37% by Weight (With Preservative)	Certified ACS	4L	Poly Bottle	F79P-4
Formaldehyde, 37% by Weight (With Preservative)	Certified ACS	20L	Rigid Poly Bottle/Corrugated Box	F79-20
Formaldehyde, 37% by Weight (With Preservative)	Certified ACS	20L	PolyPac	F79P-20
Formaldehyde, 37% by Weight (With Preservative)	Certified ACS	200L	Poly Drum	F79-200
Formamide	Certified ACS	1L	Amber Glass	F84-1
Glycerol	Certified ACS	500mL	Amber Glass	G33-500
Glycerol	Certified ACS	1L	Amber Glass	G33-1
Glycerol	Certified ACS	4L	Amber Glass	G33-4
Glycerol	Certified ACS	20L	Rigid Poly Bottle/Corrugated Box	G33-20
Glycerol	Certified ACS	200L	Steel Drum	G33-200
Hexanes	Certified ACS	500mL	Amber Glass	H292-500
Hexanes	Certified ACS	1L	Amber Glass	H292-1
Hexanes	Certified ACS	4L	Amber Glass	H292-4
Hexanes	Certified ACS	4L	Amber Glass/Safe-Cote	H292SK-4
Hexanes	Certified ACS	20L	Steel Pail	H292-20
Hexanes	Certified ACS	28L	FisherPak	H292SS-28
Hexanes	Certified ACS	50L	FisherPak	H292SS-50
Hexanes	Certified ACS	115L	FisherPak	H292SS-115
Hexanes	Certified ACS	200L	Steel Drum	H292-200
Hexanes	Certified ACS	200L	FisherPak	H292SS-200
Isobutyl Alcohol	Certified ACS	1L	Amber Glass	A397-1
Isobutyl Alcohol	Certified ACS	4L	Amber Glass	A397-4
Isooctane	Certified ACS	1L	Amber Glass	O299-1
Isooctane	Certified ACS	4L	Amber Glass	O299-4
Isooctane	Certified ACS	115L	FisherPak	O299RS115
Methanol	Certified ACS	500mL	Amber Glass	A412-500
Methanol	Certified ACS	1L	Amber Glass	A412-1
Methanol	Certified ACS	4L	Amber Glass	A412-4

Certified ACS (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Methanol	Certified ACS	4L	Poly Bottle	A412P-4
Methanol	Certified ACS	4L	Amber Glass/Safe-Cote	A412SK-4
Methanol	Certified ACS	20L	Unlined Steel Pail	A412-20
Methanol	Certified ACS	20L	Unichrome Steel Pail	A434-20
Methanol	Certified ACS	200L	Steel Drum	A412-200
Methyl Ethyl Ketone	Certified ACS	500mL	Amber Glass	M209-500
Methyl Ethyl Ketone	Certified ACS	1L	Amber Glass	M209-1
Methyl Ethyl Ketone	Certified ACS	4L	Amber Glass	M209-4
Methyl Ethyl Ketone	Certified ACS	4L	SafeTin	M209S-4
Methyl Ethyl Ketone	Certified ACS	20L	Steel Pail	M209-20
Methyl Ethyl Ketone	Certified ACS	200L	Steel Drum	M209-200
Methyl iso-Butyl Ketone	Certified ACS	1L	Amber Glass	M213-1
Methyl iso-Butyl Ketone	Certified ACS	4L	Amber Glass	M213-4
Methyl iso-Butyl Ketone	Certified ACS	20L	Steel Pail	M213-20
Methyl iso-Butyl Ketone	Certified ACS	200L	Steel Drum	M213-200
Methylene Chloride (Stabilized)	Certified ACS	500mL	Amber Glass	D37-500
Methylene Chloride (Stabilized)	Certified ACS	1L	Amber Glass	D37-1
Methylene Chloride (Stabilized)	Certified ACS	4L	Amber Glass	D37-4
Methylene Chloride (Stabilized)	Certified ACS	4L	Amber Glass/Safe-Cote	D37SK-4
Methylene Chloride (Stabilized)	Certified ACS	20L	Steel Pail	D37-20
Methylene Chloride (Stabilized)	Certified ACS	200L	Steel Drum	D37-200
Morpholine	Certified ACS	1L	Amber Glass	M263-1
Nitrobenzene	Certified ACS	500mL	Amber Glass/Poison Pack	N911-500
Nitrobenzene	Certified ACS	4L	Amber Glass/Poison Pack	N911-4
Nitromethane	Certified ACS	500mL	Amber Glass	N98-500
Petroleum Ether	Certified ACS	500mL	Amber Glass	E139-500
Petroleum Ether	Certified ACS	1L	Amber Glass	E139-1
Petroleum Ether	Certified ACS	4L	Amber Glass	E139-4
Petroleum Ether	Certified ACS	4L	SafeTin	E139S-4
Petroleum Ether	Certified ACS	4L	Amber Glass/Safe-Cote	E139SK-4
Petroleum Ether	Certified ACS	20L	Steel Pail	E139-20
Petroleum Ether	Certified ACS	200L	Steel Drum	E139-200
Pyridine	Certified ACS	500mL	Amber Glass	P368-500
Pyridine	Certified ACS	1L	Amber Glass	P368-1
Pyridine	Certified ACS	4L	Amber Glass	P368-4
Pyridine	Certified ACS	4L	SafeTin	P368S-4
Toluene	Certified ACS	500mL	Amber Glass	T324-500
Toluene	Certified ACS	1L	Amber Glass	T324-1
Toluene	Certified ACS	4L	Amber Glass	T324-4
Toluene	Certified ACS	4L	SafeTin	T324S-4

Certified ACS (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Toluene	Certified ACS	4L	Amber Glass/Safe-Cote	T324SK-4
Toluene	Certified ACS	20L	Steel Pail	T324-20
Toluene	Certified ACS	200L	Steel Drum	T324-200
Trichloroethylene (Stabilized)	Certified ACS	500mL	Amber Glass	T341-500
Trichloroethylene (Stabilized)	Certified ACS	4L	Amber Glass	T341-4
Trichloroethylene (Stabilized)	Certified ACS	20L	Steel Pail	T341-20
Xylene	Certified ACS	500mL	Amber Glass	X5-500
Xylene	Certified ACS	1L	Amber Glass	X5-1
Xylene	Certified ACS	1 gal.	Poly Bottle	X5P-1GAL
Xylene	Certified ACS	4L	Amber Glass	X5-4
Xylene	Certified ACS	4L	SafeTin	X5S-4
Xylene	Certified ACS	4L	Amber Glass/Safe-Cote	X5SK-4
Xylene	Certified ACS	20L	Steel Pail	X5-20
Xylene	Certified ACS	200L	Steel Drum	X5-200

Certified Solvents

Certified Solvents meet the purity standard established by Fisher Scientific and are manufactured for general analytical applications.

- Purity is guaranteed to meet published maximum limits on impurities
- Actual lot analysis on label
- Certificate of Analysis available on request

Certified

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetonitrile	Certified	500mL	Amber Glass	O1034-500
Acetophenone	Certified	500mL	Amber Glass	A22-500
AcetylAcetone	Certified	500mL	Amber Glass	A25-500
tert-AmylAlcohol	Certified	1L	Amber Glass	A730-1
Benzyl Alcohol	Certified	500mL	Amber Glass	A396-500
Benzyl Alcohol	Certified	4L	Amber Glass	A396-4
Benzyl Alcohol	Certified	200L	Steel Drum	A396-200
sec-Butanol	Certified	1L	Amber Glass	O1664-1
sec-Butanol	Certified	4L	Amber Glass	O1664-4
tert-Butanol	Certified	500mL	Poly Bottle	A401-500
tert-Butanol	Certified	1L	Poly Bottle	A401-1
n-Butyl Bromide	Certified	500mL	Amber Glass	B400-500
n-Butyl Chloride	Certified	1L	Amber Glass	B416-1
n-Butyl Chloride	Certified	4L	Amber Glass	B416-4
n-Butylamine	Certified	500mL	Amber Glass	B415-500
Chlorobenzene	Certified	500mL	Amber Glass	B255-500
Chlorobenzene	Certified	1L	Amber Glass	B255-1
Chlorobenzene	Certified	200L	FisherPak	B254RS200
Cyclohexanone	Certified	1L	Amber Glass	O2109-1

Certified (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Cyclohexanone	Certified	4L	Amber Glass	02109-4
Cyclohexanone	Certified	200L	Stainless Steel Drum	02109SS200
2-Chloro-2-methylpropane	Certified	500mL	Amber Glass	01875-500
Decane	Certified	500mL	Amber Glass	02128-500
o-Dichlorobenzene	Certified	1L	Amber Glass	02231-1
Diisopropylamine	Certified	4L	Amber Glass	02412-4
1,2-Dimethoxyethane	Certified	1L	Amber Glass	02430-1
1,2-Dimethoxyethane	Certified	4L	Amber Glass	02430-4
N,N-Dimethylacetamide	Certified	500mL	Amber Glass	D135-500
Ethylbenzene	Certified	1L	AmberGlass	02751-1
Ethylene Glycol	Certified	500mL	Amber Glass	E178-500
Ethylene Glycol	Certified	1L	Amber Glass	E178-1
Ethylene Glycol	Certified	4L	Amber Glass	E178-4
Ethylene Glycol	Certified	200L	Steel Drum	E178-200
Ethylene Glycol Monomethyl Ether	Certified	500mL	Amber Glass	E182-500
Ethylene Glycol Monomethyl Ether	Certified	4L	Amber Glass	E182-4
Ethylene Glycol Monomethyl Ether	Certified	20L	Steel Pail	E182-20
Formalin, Buffered, 10% (Acetate Buffer)	Certified	4L	Poly Bottle	SF99-4
Formalin, Buffered, 10% (Acetate Buffer)	Certified	20L	PolyPac	SF99-20
Heptane	Certified	1L	Amber Glass	03008-1
Heptane	Certified	4L	Amber Glass	03008-4
Heptane	Certified	115L	FisherPak	03008RS115
Hexanes	Certified	500mL	Amber Glass	H291-500
Hexanes	Certified	4L	Amber Glass	H291-4
Hexanes	Certified	4L	SafeTin	H291S-4
Hexanes	Certified	20L	Steel Pail	H291-20
Hexanes	Certified	200L	Steel Drum	H291-200
Isoamyl Alcohol	Certified	500mL	Amber Glass	A393-500
Isoamyl Alcohol	Certified	4L	Amber Glass	A393-4
Isopropanol, 10% (v/v)	Certified	1 gal.	Poly Bottle	A460-1GAL
Isopropyl Acetate	Certified	1L	Poly Bottle	061111
Isopropyl Acetate	Certified	2.5L	Poly Bottle	06111-212
Isopropyl Ether	Certified	500mL	Amber Glass	E141-500
Isopropyl Ether	Certified	4L	Amber Glass	E141-4
Methylal	Certified	500mL	Amber Glass	M222-500
Methyl Acetate	Certified	500mL	Amber Glass	M203-500
Methyl Acetate	Certified	4L	Amber Glass	M203-4
2-Methylbutane	Certified	4L	Amber Glass	03551-4
1-Octanol	Certified	500mL	Amber Glass	A402-500

Certified (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
1-Octanol	Certified	4L	Amber Glass	A402-4
Pentane	Certified	4L	Amber Glass	O4062-4
Pentane	Certified	19L	FisherPak	O4062RS19
Pentane	Certified	20L	Steel Pail	O4062-20
Petroleum Ether	Certified	115L	FisherPak	E120SS-115
1-Propanol	Certified	500mL	Amber Glass	A414-500
1-Propanol	Certified	1L	Amber Glass	A414-1
1-Propanol	Certified	4L	Amber Glass	A414-4
1-Propanol	Certified	4L	SafeTin	A414S-4
1-Propanol	Certified	20L	Steel Pail	A414-20
1-Propanol	Certified	50L	BasicPak	A414RB50
2-Propanol	Certified	1L	Amber Glass	A417-1
2-Propanol	Certified	4L	Amber Glass	A417-4
1,1,2,2-Tetrabromoethane	Certified	100mL	Amber Glass	A29-100
Tetrahydrofuran	Certified	500mL	Amber Glass	T397-500
Tetrahydrofuran	Certified	1L	Amber Glass	T397-1
Tetrahydrofuran	Certified	4L	Amber Glass	T397-4
Tetrahydrofuran	Certified	4L	Amber Glass/Safe-Cote	T397SK-4
Tetrahydrofuran	Certified	19L	BasicPak	T397RB19
Tetrahydrofuran	Certified	20L	Steel Pail	T397-20
Tetrahydrofuran	Certified	50L	FisherPak	T397RS50
Tetrahydrofuran	Certified	200L	Steel Drum	T397-200
Tetrahydrofuran	Certified	200L	BasicPak	T397RB200
Tetrahydrofuran	Certified	200L	FisherPak	T397RS200
Triethanolamine	Certified	500mL	Amber Glass	T407-500
Triethanolamine	Certified	1L	Amber Glass	T407-1
Triethanolamine	Certified	4L	Amber Glass	T407-4
o-Xylene	Certified	4L	Amber Glass	O5081-4
p-Xylene	Certified	500mL	Amber Glass	O5082-500
p-Xylene	Certified	4L	Amber Glass	O5082-4

Pesticide Grade Solvents

Pesticide Grade Solvents are used for analysis of pesticide residue and in GC with electron capture detector.

- Meet or exceed ACS standards of purity for pesticide residue analysis
- Actual lot analysis on label

Pesticide

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetone	Pesticide	4L	Amber Glass	A40-4
Acetonitrile	Pesticide	4L	Amber Glass	A999-4
Chloroform (With Amylene Preservative)	Pesticide	4L	Amber Glass	C603-4
Cyclohexane	Pesticide	4L	Amber Glass	C553-4

Pesticide (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Ethyl Acetate	Pesticide	4L	Amber Glass	E191-4
Ethyl Ether	Pesticide	4L	Amber Glass	E199-4
Hexanes	Pesticide	4L	Amber Glass	H300-4
Isooctane	Pesticide	4L	Amber Glass	O297-4
Methanol	Pesticide	4L	Amber Glass	A450-4
Methylene Chloride	Pesticide	4L	Amber Glass	D142-4
Pentane	Pesticide	4L	Amber Glass	P400-4
Petroleum Ether	Pesticide	4L	Amber Glass	P480-4
2-Propanol	Pesticide	4L	Amber Glass	A519-4

Spectranalyzed Solvents

Spectranalyzed Solvents are for use in spectrophotometry—ultraviolet and visible wavelength detectors.

- Meet all ACS specifications
- Actual lot analysis on label
- Certificate of Analysis available on request

Spectranalyzed

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetone	Spectranalyzed	1L	Amber Glass	A19-1
Acetone	Spectranalyzed	4L	Amber Glass	A19-4
Acetone	Spectranalyzed	115L	FisherPak	A19RS115
1-Butanol	Spectranalyzed	4L	Amber Glass	A400-4
Carbon Disulfide	Spectranalyzed	500mL	Amber Glass	C573-500
Chloroform (Approx. 0.75% Ethanol Preservative)	Spectranalyzed	1L	Amber Glass	C574-1
Chloroform (Approx. 0.75% Ethanol Preservative)	Spectranalyzed	4L	Amber Glass	C574-4
Chloroform (Approx. 0.75% Ethanol Preservative)	Spectranalyzed	4L	Amber Glass/Safe-Cote	C574SK-4
Cyclohexane	Spectranalyzed	1L	Amber Glass	C555-1
Cyclohexane	Spectranalyzed	4L	Amber Glass	C555-4
1,2-Dichloroethane	Spectranalyzed	4L	Amber Glass	E190-4
N,N-Dimethylformamide	Spectranalyzed	1L	Amber Glass	D131-1
N,N-Dimethylformamide	Spectranalyzed	4L	Amber Glass	D131-4
Dimethyl Sulfoxide	Spectranalyzed	1L	Amber Glass	D136-1
Ethyl Acetate	Spectranalyzed	4L	Amber Glass	E189-4
Ethyl Ether	Spectranalyzed	1L	Tin Can	E197-1
Ethyl Ether	Spectranalyzed	4L	Tin Can	E197-4
Glycerol	Spectranalyzed	1L	Amber Glass	G153-1
Glycerol	Spectranalyzed	4L	Amber Glass	G153-4
Heptane	Spectranalyzed	4L	Amber Glass	H340-4
Hexanes	Spectranalyzed	1L	Amber Glass	H334-1
Hexanes	Spectranalyzed	4L	Amber Glass	H334-4
Isooctane	Spectranalyzed	1L	Amber Glass	O300-1
Isooctane	Spectranalyzed	4L	Amber Glass	O300-4

Spectranalyzed (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Methanol	Spectranalyzed	1L	Amber Glass	A408-1
Methanol	Spectranalyzed	4L	Amber Glass	A408-4
Methanol	Spectranalyzed	4L	Amber Glass/Safe-Cote	A408SK-4
Methylene Chloride (Stabilized)	Spectranalyzed	1L	Amber Glass	D35-1
Methylene Chloride (Stabilized)	Spectranalyzed	4L	Amber Glass	D35-4
Pentane	Spectranalyzed	1L	Amber Glass	P393-1
2-Propanol	Spectranalyzed	1L	Amber Glass	A419-1
2-Propanol	Spectranalyzed	4L	Amber Glass	A419-4
Tetrahydrofuran	Spectranalyzed	4L	Amber Glass	T424-4
Toluene	Spectranalyzed	4L	Amber Glass	T330-4

Scintanalyzed Solvents

Scintanalyzed Solvents are specially formulated and tested for liquid scintillation counting.

- Certificate of Analysis available on request

Scintanalyzed

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
1,4-Dioxane	Scintanalyzed	4L	Amber Glass	D116-4
1,4-Dioxane	Scintanalyzed	200L	Steel Drum	D116-200
Toluene	Scintanalyzed	4L	Amber Glass	T313-4
Toluene	Scintanalyzed	4L	Amber Glass/Safe-Cote	T313SK-4
Xylene	Scintanalyzed	4L	Amber Glass	X16-4

Electronic Grade Solvents

Electronic Grade Solvents are manufactured to ensure low levels of metal contamination and are suitable for use in electronics and circuit board manufacturing.

- Certificate of Analysis available on request

Electronic

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetone	Electronic	4L	Amber Glass	A946-4
Methanol	Electronic	4L	Poly Bottle	A947-4
Trichloroethylene	Electronic	4L	Amber Glass	T403-4

Histological Grade Solvents

Histological Grade Solvents are specially prepared for use in the histology laboratory setting and are filtered for tissue processing applications.

- Certificate of Analysis available on request

Histological

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetone	Histological	1 gal.	Poly Bottle	A16F1GAL
Acetone	Histological	4L	Poly Bottle	A16P-4
Acetone	Histological	4L	SafeTin	A16S-4

Histological (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetone	Histological	20L	Steel Pail	A16S-20
Alcohol	Histological	1 gal.	Rectangular Poly Bottle	A962F1GAL
Alcohol	Histological	4L	Amber Glass	A962-4
Alcohol	Histological	4L	Poly Bottle	A962P-4
Alcohol	Histological	4L	SafeTin	A962S-4
Alcohol	Histological	200L	Steel Drum	A962-200
Alcohol	Histological	200L	BasicPak	A962RB200
Alcohol	Histological	200L	FisherPak	A992RS200
1,4-Dioxane	Histological	4L	SafeTin	D56S-4
Ethanol, Anhydrous	Histological	1 gal.	Poly Bottle	A405F1GAL
Ethanol, Anhydrous	Histological	4L	Poly Bottle	A405P-4
Ethanol, Anhydrous	Histological	20L	Steel Pail	A405-20
Ethanol, CDA19	Histological	1 gal.	Poly Bottle	A406F1GAL
Ethanol, CDA19	Histological	4L	Poly Bottle	A406P-4
Ethanol, CDA19	Histological	20L	Steel Pail	A406-20
Formaldehyde, 37% by Weight	Histological	1 gal.	Poly Bottle	F75F1GAL
Formaldehyde, 37% by Weight	Histological	1 gal.	Poly Bottle	F75P1GAL
Formaldehyde, 37% by Weight	Histological	20L	PolyPac	F75P-20
Formalin Solution, 10%	Histological	4L	Poly Bottle	SF98-4
Formalin Solution, 10%	Histological	20L	PolyPac	SF98-20
Methanol	Histological	1 gal.	Poly Bottle	A433F1GAL
Methanol	Histological	4L	Poly Bottle	A433P-4
Methanol	Histological	4L	SafeTin	A433S-4
Methanol	Histological	20L	Steel Pail	A433S-20
Methanol	Histological	200L	Steel Drum	A433S-200
2-Propanol	Histological	1 gal.	Poly Bottle	A426F1GAL
2-Propanol	Histological	4L	Poly Bottle	A426P-4
2-Propanol	Histological	4L	SafeTin	A426S-4
2-Propanol	Histological	20L	Steel Pail	A426S-20
2-Propanol	Histological	200L	Steel Drum	A426S-200
Tetrahydrofuran	Histological	500mL	Amber Glass	T400-500
Tetrahydrofuran	Histological	4L	Amber Glass	T400-4
Tetrahydrofuran	Histological	20L	Steel Pail	T400-20
Toluene	Histological	1 gal.	Poly Bottle	T326F1GAL
Toluene	Histological	4L	Poly Bottle	T326P-4
Toluene	Histological	20L	Steel Pail	T326S20
Xylene	Histological	1 gal.	Poly	X3F1GAL
Xylene	Histological	1 gal.	Poly Bottle	X3P-1GAL
Xylene	Histological	4L	SafeTin	X3S-4
Xylene	Histological	20L	Steel Pail	X3S-20
Xylene	Histological	50L	BasicPak	X3RB50
Xylene	Histological	200L	Steel Drum	X3S-200

Laboratory Grade Solvents

Laboratory Grade Solvents are of reasonable quality and purity for use in procedures where no official standards are required.

- Certificate of Analysis available on request

Laboratory

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Amyl Acetate	Laboratory	500mL	Amber Glass	A718-500
Amyl Acetate	Laboratory	4L	Amber Glass	A718-4
n-Butyl Phosphate	Laboratory	500mL	Amber Glass	B404-500
n-Butyl Phosphate	Laboratory	4L	Amber Glass	B404-4
Carbon Disulfide	Laboratory	2.5L	Glass	C183-212
Chlorobenzene	Laboratory	4L	Amber Glass	B254-4
Chlorobenzene	Laboratory	20L	Steel Pail	B254-20
Chlorobenzene	Laboratory	200L	FisherPak	B254RS200
Cyclohexanone	Laboratory	4L	Amber Glass	C550-4
Ethanolamine	Laboratory	1L	Amber Glass	M251-1
Ethanolamine	Laboratory	4L	Amber Glass	M251-4
Ethylene Glycol Monoethyl Ether Acetate	Laboratory	4L	Amber Glass	E181-4
Ethylene Glycol Monoethyl Ether	Laboratory	1L	Amber Glass	E180-1
Ethylene Glycol Monoethyl Ether	Laboratory	4L	Amber Glass	E180-4
Ethylene Glycol Monoethyl Ether	Laboratory	20L	Steel Pail	E180-20
Ethyl Ether	Laboratory	1L	Aluminum Can	E134-1
Ethyl Ether	Laboratory	4L	Tin Can	E134-4
Ethyl Ether	Laboratory	20L	Steel Pail	E134-20
Ethyl Ether, Fat Extraction	Laboratory	4L	Tin Can	E492-4
Ethyl Ether, Fat Extraction	Laboratory	20L	Steel Pail	E492-20
Ethylene Glycol	Laboratory	4L	Amber Glass	E177-4
Ethylene Glycol	Laboratory	20L	Steel Pail	E177-20
Ligroine, 100% Heptane	Laboratory	4L	Amber Glass	O3387-4
Ligroine, Hexane	Laboratory	20L	Steel Pail	O3386-20
Methanol	Laboratory	4L	Amber Glass	A411-4
Methanol	Laboratory	20L	Steel Pail	A411-20
2-Octanol	Laboratory	500mL	Amber Glass	O269-500
Petroleum Ether	Laboratory	115L	FisherPak	E120SS-115
1-Propanol	Laboratory	50L	BasicPak	A414RB50
2-Propanol	Laboratory	4L	Amber Glass	A415-4
2-Propanol	Laboratory	20L	Steel Pail	A415-20
1,1,2,2-Tetrabromoethane	Laboratory	500mL	Amber Glass	A33-500
1,1,2,2-Tetrabromoethane	Laboratory	4L	Amber Glass	A33-4
Tetrahydrofuran	Laboratory	19L	BasicPak	T397RB19
Tetrahydrofuran	Laboratory	50L	FisherPak	T397RS50
Tetrahydrofuran	Laboratory	200L	BasicPak	T397RB200

Laboratory (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Tetrahydrofuran	Laboratory	200L	FisherPak	T397RS200
Toluene	Laboratory	4L	Amber Glass	T323-4
Toluene	Laboratory	20L	Steel Pail	T323-20
Triethylene Glycol	Laboratory	4L	Amber Glass	T346-4
Xylene	Laboratory	4L	Amber Glass	X4-4
Xylene	Laboratory	20L	Steel Pail	X4-20

Technical Grade Solvents

Technical Grade Solvents are of reasonable quality and purity for use in procedures where no official standards are required.

- Certificate of Analysis available on request

Technical

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
1-Butanol	Technical	4L	Amber Glass	A398-4
Chlorobenzene	Technical	200L	FisherPak	B254RS200
Chloroform (Approx. 0.75% Ethanol as Preservative)	Technical	4L	Amber Glass	C295-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Technical	4L	SafeTin	C295S-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Technical	20L	Steel Pail	C295-20
Diacetone Alcohol	Technical	4L	Amber Glass	D17-4
Diethylenetriamine	Technical	500mL	Amber Glass	D126-500
Heptane	Technical	20L	Steel Pail	H20-20
Heptane	Technical	200L	Steel Drum	H20-200
Hexanes	Technical	4L	SafeTin	N3S-4
Hexanes	Technical	20L	Steel Drum	N3-20
Hexanes	Technical	200L	Steel Drum	N3-200
Methyl Ethyl Ketone	Technical	1L	Amber Glass	M208-1
Methyl Ethyl Ketone	Technical	4L	Amber Glass	M208-4
Methyl Ethyl Ketone	Technical	20L	Steel Pail	M208-20
n-Butyl Acetate	Technical	4L	Amber Glass	B395-4
Petroleum Ether	Technical	115L	FisherPak	E120SS-115
1-Propanol	Technical	50L	BasicPak	A414RB50
Pyridine	Technical	4L	Amber Glass	P369-4
Tetrachloroethylene	Technical	4L	Amber Glass	C182-4
Tetrachloroethylene	Technical	20L	Steel Pail	C182-20
Tetrahydrofuran	Technical	19L	BasicPak	T397RB19
Tetrahydrofuran	Technical	50L	FisherPak	T397RS50
Tetrahydrofuran	Technical	200L	BasicPak	T397RB200
Tetrahydrofuran	Technical	200L	FisherPak	T397RS200
Trichloroethylene	Technical	4L	Amber Glass	T340-4
Triethylenetetramine	Technical	500mL	Amber Glass	T410-500
Triethylenetetramine	Technical	1L	Amber Glass	T410-1

Reagent Grade Solvents

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
n-Butyl Acetate	Reagent	1L	Amber Glass	B396-1
n-Butyl Acetate	Reagent	4L	Amber Glass	B396-4
n-Butyl Phthalate	Reagent	500mL	Amber Glass	D30-500
n-Butyl Phthalate	Reagent	4L	Amber Glass	D30-4
Chlorobenzene	Reagent	200L	FisherPak	B254RS200
Cyclohexane	Reagent	4L	Amber Glass	O2093-4
Cyclohexane	Reagent	20L	Steel Pail	O2093-20
Cyclohexanol	Reagent	500mL	Amber Glass	C558-500
Cyclohexene (Contains 0.01% tert-Butylcresol as an Inhibitor)	Reagent	1L	Amber Glass	O2111-1
Diethylene Glycol	Reagent	1L	Amber Glass	D49-1
Ethyl Acetoacetate	Reagent	500mL	Amber Glass	E146-500
Octane	Reagent	1L	Amber Glass	O3980-1
Petroleum Ether	Reagent	115L	FisherPak	E120SS-115
1-Propanol	Reagent	50L	BasicPak	A414RB50
Propylene Oxide	Reagent	1L	Amber Glass	O4332-1
Tetrabutylammonium Iodide (White Crystals)	Reagent	100g	Poly Bottle	O6109-100
Tetrachloroethylene (Stabilized)	Reagent	4L	Amber Glass	O4586-4
Tetrahydrofuran	Reagent	19L	BasicPak	T397RB19
Tetrahydrofuran	Reagent	50L	FisherPak	T397RS50
Tetrahydrofuran	Reagent	200L	BasicPak	T397RB200
Tetrahydrofuran	Reagent	200L	FisherPak	T397RS200
Triethylamine	Reagent	1L	Amber Glass	O4885-1
Triethylamine	Reagent	4L	Amber Glass	O4885-4
Triethylamine	Reagent	20L	Steel Pail	O488520

Multicompendial Solvents

Multicompendial Solvents meet or surpass specifications of the United States Pharmacopeia (USP), the National Formulary (NF), the Food Chemicals Codex (FCC), the European Pharmacopeia (EP), the British Pharmacopeia (BP), and/or the Japanese Pharmacopeia (JP).

Multicompendial

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetone	NF/FCC	1L	Amber Glass	A11-1
Acetone	NF/FCC	4L	Amber Glass	A11-4
Acetone	NF/FCC	4L	SafeTin	A11S-4
Acetone	NF/FCC	20L	Steel Pail	A11-20
Acetone	NF/FCC	200L	Steel Drum	A11-200
Acetone	NF/FCC/EP	4L	Amber Glass	A94
Acetone	NF/FCC/EP	20L	Steel Pail	A920
Acetone	NF/FCC/EP	200L	Steel Drum	A9200
Benzyl Alcohol	NF	4L	Amber Glass	A395-4
Benzyl Alcohol	NF/EP/BP	4L	Amber Glass	A392-4
Ethyl Acetate	NF	4L	Amber Glass	E124-4

Multicompendial (continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Ethyl Acetate	NF	20L	Steel Pail	E124-20
Ethyl Acetate	NF/FCC	200L	FisherPak	E124RS200
Formaldehyde (40% by Volume)	USP	4L	Poly Bottle	F77P-4
Formaldehyde (40% by Volume)	USP	20L	Rigid Poly Bottle/Corrugated Box	F77-20
Formaldehyde (40% by Volume)	USP	20L	PolyPac	F77P-20
Formaldehyde (40% by Volume)	USP	200L	Steel Drum/Poly Liner	F77-200
Glycerol	USP/FCC	500mL	Amber Glass	G31-500
Glycerol	USP/FCC	1L	Amber Glass	G31-1
Glycerol	USP/FCC	4L	Amber Glass	G31-4
Glycerol	USP/FCC	20L	Rigid Poly Bottle	G31-20
Glycerol	USP/FCC	200L	Poly Drum	G31-200
Glycerol	USP/FCC/EP/BP	4L	Amber Glass	G30-4
Glycerol	USP/FCC/EP/BP	20L	Rigid Poly Bottle/Corrugated Box	G30-20
Glycerol	USP/FCC/EP/BP	200L	Steel Drum	G30200
Methanol	NF	500mL	Amber Glass	A413500
Methanol	NF	4L	Amber Glass	A4134
Methanol	NF	20L	Steel Pail	A41320
Methanol	NF	200L	Steel Drum	A413-200
Pentane	Multicompendial	28L	FisherPak	P399RS28
2-Propanol	USP	500mL	Amber Glass	A516500
2-Propanol	USP	4L	Amber Glass	A516-4
2-Propanol	USP	20L	Steel Pail	A516-20
2-Propanol	USP	200L	Steel Drum	A516-200
Propylene Glycol	USP/FCC	1L	Amber Glass	P355-1
Propylene Glycol	USP/FCC	4L	Amber Glass	P355-4
Propylene Glycol	USP/FCC	20L	Steel Pail	P355-20
Propylene Glycol	USP/FCC	200L	Steel Drum	P355-200
Triethanolamine	NF	500mL	Amber Glass	T350-500
Triethanolamine	NF	4L	Amber Glass	T350-4

Special Solvent Grades

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetonitrile, 50% (v/v)	For AOAC Method 992.22	500mL	Amber Glass	SA9125-500
Acetonitrile, 50% (v/v)	For AOAC Method 992.22	4L	Amber Glass	SA9125-4
Alcohol Reagent 20%	For AOAC Method 992.22	1L	Poly Bottle	SA9320-1
Alcohol Reagent 20%	For AOAC Method 992.22	4L	Poly Bottle	SA9320-4
Alcohol Reagent 20%	For AOAC Method 992.22	10L	Cubitainer	SA9320-10
Benzine (Petroleum Naphtha)		20L	Steel Pail	B264-20
Carbon Disulfide (Low Benzene)		500mL	Amber Glass	C185-500
Ethyl Alcohol Denatured	Proprietary Solvent	500mL	Amber Glass	A407-500
Ethyl Alcohol Denatured	Proprietary Solvent	1L	Amber Glass	A407-1
Ethyl Alcohol Denatured	Proprietary Solvent	4L	Amber Glass	A407-4

Special Solvent Grades (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Ethyl Alcohol Denatured	Proprietary Solvent	4L	Poly Bottle	A407P-4
Ethyl Alcohol Denatured	Proprietary Solvent	4L	SafeTin	A407S-4
Ethyl Alcohol Denatured	Proprietary Solvent	4L	Amber Glass/Safe-Cote	A407SK-4
Ethyl Alcohol Denatured	Proprietary Solvent	19L	BasicPak	A407RB19
Ethyl Alcohol Denatured	Proprietary Solvent	20L	Steel Pail	A407-20
Ethyl Alcohol Denatured	Proprietary Solvent	115L	BasicPak	A407RB115
Ethyl Alcohol Denatured	Proprietary Solvent	200L	Steel Drum	A407-200
Ethyl Alcohol Denatured	Proprietary Solvent	200L	BasicPak	A407RB200
IPA (Isopropyl Alcohol), 70% (v/v)		500mL	Poly Bottle	A459-500
IPA (Isopropyl Alcohol), 70% (v/v)		1L	Poly Bottle	A459-1
IPA (Isopropyl Alcohol), 70% (v/v)		4L	Poly Bottle	A459-4
IPA (Isopropyl Alcohol), 70% (v/v)		20L	PolyPac	A459-20
Kerosene (Odorless)		4L	Amber Glass	K10-4
Kerosene (Odorless)		200L	Steel Drum	K10-200
Methanol	Low Water	4L	Amber Glass	A935-4
Methanol	Purge and Trap Suitable for Volatile Organic Residue Analysis	500mL	Amber Glass	A453-500
Methanol	Purge and Trap Suitable for Volatile Organic Residue Analysis	1L	Amber Glass	A453-1
2-Propanol	Low Water	4L	Amber Glass	A520-4
Stoddard Solvent		4L	Amber Glass	S457-4
Stoddard Solvent		200L		S457200
Water	DIUF	4L	Poly Bottle	W2-4
Water	DIUF	20L	PolyPac	W2-20
Water	Environmental Grade	1L	Amber Glass	W11-1
Water	Environmental Grade	4L	Amber Glass	W11-4
Water	Plasma Grade	500mL	LDPE Bottle	W9-500
Water	Plasma Grade	1L	LDPE Bottle	W9-1
Water	Plasma Grade	2L	LDPE Bottle	W9-2

Solvents (in Alphabetical Order)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetic Anhydride	Certified ACS	100mL	Amber Glass	A10-100
Acetic Anhydride	Certified ACS	500mL	Amber Glass	A10-500
Acetic Anhydride	Certified ACS	1L	Amber Glass	A10-1
Acetic Anhydride	Certified ACS	4L	Amber Glass	A10-4
Acetone	Certified ACS	500mL	Amber Glass	A18-500
Acetone	Certified ACS	1L	Amber Glass	A18-1
Acetone	HPLC	1L	Amber Glass	A949-1
Acetone	HPLC	1L	Amber Glass/Safe-Cote	A949SK-1
Acetone	NF/FCC	1L	Amber Glass	A11-1
Acetone	Optima	1L	Amber Glass	A929-1
Acetone	Spectranalyzed	1L	Amber Glass	A19-1
Acetone	Histological	1 gal.	Poly Bottle	A16F1GAL
Acetone	Certified ACS	4L	Amber Glass	A18-4
Acetone	Certified ACS	4L	Poly Bottle	A18P-4
Acetone	Certified ACS	4L	SafeTin	A18S-4
Acetone	Certified ACS	4L	Amber Glass/Safe-Cote	A18SK-4
Acetone	Electronic Grade	4L	Amber Glass	A946-4
Acetone	GC <i>Resolv</i>	4L	Amber Glass	A928-4
Acetone	Histological	4L	Poly Bottle	A16P-4
Acetone	Histological	4L	SafeTin	A16S-4
Acetone	HPLC	4L	Amber Glass	A949-4
Acetone	HPLC	4L	Amber Glass/Safe-Cote	A949SK-4
Acetone	NF/FCC	4L	Amber Glass	A11-4
Acetone	NF/FCC	4L	SafeTin	A11S-4
Acetone	NF/FCC/EP	4L	Amber Glass	A94
Acetone	HPLC	19L	NOWPak I	A949N1-19
Acetone	Optima	4L	Amber Glass	A929-4
Acetone	Optima	4L	Amber Glass/Safe-Cote	A929SK-4
Acetone	Pesticide Grade	4L	Amber Glass	A40-4
Acetone	Spectranalyzed	4L	Amber Glass	A19-4
Acetone	Certified ACS	20L	Steel Pail	A18-20
Acetone	Histological	20L	Steel Pail	A16S-20
Acetone	NF/FCC	20L	Steel Pail	A11-20
Acetone	NF/FCC/EP	20L	Steel Pail	A920
Acetone	Certified ACS	200L	Steel Drum	A18-200
Acetone	NF/FCC	200L	Steel Drum	A11-200
Acetone	NF/FCC/EP	200L	Steel Drum	A9200
Acetone	Certified ACS	350 gal.	Steel Tote	A18CU1300
Acetonitrile	Certified	500mL	Amber Glass	01034-500
Acetonitrile	Optima LC/MS	500mL	Amber Glass	A955-500
Acetonitrile	Certified ACS	1L	Amber Glass	A21-1
Acetonitrile	HPLC	1L	Amber Glass	A998-1
Acetonitrile	HPLC	1L	Amber Glass/Safe-Cote	A998SK-1
Acetonitrile	Optima	1L	Amber Glass	A996-1
Acetonitrile	Optima LC/MS	1L	Amber Glass	A955-1
Acetonitrile	HPLC	2.5L	Amber Glass	A998212
Acetonitrile	Certified ACS	4L	Amber Glass	A21-4
Acetonitrile	HPLC	4L	Amber Glass	A998-4
Acetonitrile	HPLC	4L	Amber Glass/Safe-Cote	A998SK-4
Acetonitrile	Optima	4L	Amber Glass	A996-4
Acetonitrile	Optima	4L	Amber Glass/Safe-Cote	A996SK4
Acetonitrile	Optima LC/MS	4L	Amber Glass	A955-4

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Acetonitrile	Pesticide	4L	Amber Glass	A999-4
Acetonitrile	HPLC	19L	NOWPak I	A998N1-19
Acetonitrile	Certified ACS	20L	Steel Pail	A21-20
Acetonitrile	Certified ACS	200L	Steel Drum	A21-200
Acetonitrile, 50% (v/v)	For AOAC Method 992.22	500mL	Amber Glass	SA9125-500
Acetonitrile, 50% (v/v)	For AOAC Method 992.22	4L	Amber Glass	SA9125-4
0.1% FA in Acetonitrile	OPTIMA LC/MS	500mL	Amber Glass	LS120-500
0.1% FA in Acetonitrile	OPTIMA LC/MS	1L	Amber Glass	LS120-1
0.1% FA in Acetonitrile	OPTIMA LC/MS	2.5L	Amber Glass	LS120-212
0.1% FA in Acetonitrile	OPTIMA LC/MS	4L	Amber Glass	LS120-4
0.1% TFA in Acetonitrile	OPTIMA LC/MS	500mL	Amber Glass	LS121-500
0.1% TFA in Acetonitrile	OPTIMA LC/MS	1L	Amber Glass	LS121-1
0.1% TFA in Acetonitrile	OPTIMA LC/MS	2.5L	Amber Glass	LS121-212
0.1% TFA in Acetonitrile	OPTIMA LC/MS	4L	Amber Glass	LS121-4
Acetophenone	Certified	500mL	Amber Glass	A22-500
Acetyl Acetone	Certified	500mL	Amber Glass	A25-500
Alcohol	Histological	1 gal.	Rectangular Poly Bottle	A962F1GAL
Alcohol	Histological	4L	Amber Glass	A962-4
Alcohol	Histological	4L	Poly Bottle	A962P-4
Alcohol	Histological	4L	SafeTin	A962S-4
Alcohol	HPLC	4L	Amber Glass	A995-4
Alcohol	Histological	200L	Steel Drum	A962-200
Alcohol Reagent 20%	For AOAC Method 992.22	1L	Poly Bottle	SA9320-1
Alcohol Reagent 20%	For AOAC Method 992.22	4L	Poly Bottle	SA9320-4
Alcohol Reagent 20%	For AOAC Method 992.22	10L	Cubitainer	SA9320-10
Amyl Acetate	Laboratory	500mL	Amber Glass	A718-500
Amyl Acetate	Laboratory	4L	Amber Glass	A718-4
n-Amyl Alcohol	Certified ACS	500mL	Amber Glass	A394-500
n-Amyl Alcohol	Certified ACS	4L	Amber Glass	A394-4
tert-Amyl Alcohol	Certified	1L	Amber Glass	A730-1
Aniline	Certified ACS	500mL	Amber Glass/Poison Pack	A740I-500
Aniline	Certified ACS	4L	Amber Glass/Poison Pack	A740I-4
Benzene (Petroleum Naphtha)		20L	Steel Pail	B264-20
Benzyl Alcohol	Certified	500mL	Amber Glass	A396-500
Benzyl Alcohol	NF	500mL	Amber Glass	A395500
Benzyl Alcohol	Certified	4L	Amber Glass	A396-4
Benzyl Alcohol	NF	4L	Amber Glass	A3954
Benzyl Alcohol	NF/EP/BP	4L	Amber Glass	A392-4
Benzyl Alcohol	Certified	200L	Steel Drum	A396-200
1-Butanol	Certified ACS	500mL	Amber Glass	A399-500
1-Butanol	Certified ACS	1L	Amber Glass	A399-1
1-Butanol	HPLC	1L	Amber Glass	A383-1
1-Butanol	HPLC	1L	Amber Glass/Safe-Cote	A383SK-1
1-Butanol	Certified ACS	4L	Amber Glass	A399-4
1-Butanol	Certified ACS	4L	SafeTin	A399S-4
1-Butanol	HPLC	4L	Amber Glass	A383-4

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
1-Butanol	HPLC	4L	Amber Glass/Safe-Cote	A383SK-4
1-Butanol	Spectranalyzed	4L	Amber Glass	A400-4
1-Butanol	Technical	4L	Amber Glass	A398-4
1-Butanol	Certified ACS	20L	Steel Pail	A399-20
sec-Butanol	Certified	1L	Amber Glass	O1664-1
tert-Butanol	Certified	500mL	Poly Bottle	A401-500
n-Butyl Acetate	Reagent	1L	Amber Glass	B396-1
n-Butyl Acetate	Reagent	4L	Amber Glass	B396-4
n-Butyl Acetate	Technical	4L	Amber Glass	B395-4
n-Butyl Bromide	Certified	500mL	Amber Glass	B400-500
n-Butyl Chloride	Certified	1L	Amber Glass	B416-1
n-Butyl Chloride	Certified	4L	Amber Glass	B416-4
n-Butyl Chloride	HPLC	4L	Amber Glass/Safe-Cote	B429SK-4
n-Butyl Phosphate	Laboratory	500mL	Amber Glass	B404-500
n-Butyl Phosphate	Laboratory	4L	Amber Glass	B404-4
n-Butyl Phthalate	Reagent	500mL	Amber Glass	D30-500
n-Butyl Phthalate	Reagent	4L	Amber Glass	D30-4
n-Butylamine	Certified	500mL	Amber Glass	B415-500
Carbon Disulfide	Certified ACS	2.5L	Glass	C184-212
Carbon Disulfide	Laboratory	2.5L	Glass	C183-212
Carbon Disulfide (Low Benzene)		500mL	Amber Glass	C185-500
Chlorobenzene	Certified	500mL	Amber Glass	B255-500
Chlorobenzene	Certified	1L	Amber Glass	B255-1
Chlorobenzene	Laboratory	4L	Amber Glass	B254-4
Chlorobenzene	Laboratory	20L	Steel Pail	B254-20
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	500mL	Amber Glass	C298-500
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	1L	Amber Glass	C298-1
Chloroform (Approx. 0.75% Ethanol as Preservative)	HPLC	1L	Amber Glass	C606-1
Chloroform (Approx. 0.75% Ethanol as Preservative)	HPLC	1L	Amber Glass/Safe-Cote	C606SK-1
Chloroform (Approx. 0.75% Ethanol as Preservative)	Spectranalyzed*	1L	Amber Glass	C574-1
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	4L	Amber Glass	C298-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	4L	SafeTin	C298S-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	4L	Amber Glass/Safe-Cote	C298SK-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	HPLC	4L	Amber Glass	C606-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	HPLC	4L	Amber Glass/Safe-Cote	C606SK-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Spectranalyzed*	4L	Amber Glass	C574-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Spectranalyzed*	4L	Amber Glass/Safe-Cote	C574SK-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Technical	4L	Amber Glass	C295-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Technical	4L	SafeTin	C295S-4
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	20L	Steel Pail	C298-20
Chloroform (Approx. 0.75% Ethanol as Preservative)	Technical	20L	Steel Pail	C295-20
Chloroform (Approx. 0.75% Ethanol as Preservative)	Certified ACS	200L	Steel Drum	C298-200
Chloroform (Approx. 50ppm Amylene as Preservative)	Optima	4L	Amber Glass	C297-4
Chloroform (Approx. 50ppm Pentene as Preservative)	HPLC	1L	Amber Glass	C607-1
Chloroform (Approx. 50ppm Pentene as Preservative)	HPLC	4L	Amber Glass	C607-4
Chloroform (Approx. 50ppm Pentene as Preservative)	HPLC	4L	Amber Glass/Safe-Cote	C607SK-4
Chloroform (With Amylene Preservative)	Pesticide	4L	Amber Glass	C603-4
2-Chloro-2-methylpropane	Certified	500mL	Amber Glass	O1875-500
Cyclohexane	Certified ACS	500mL	Amber Glass	C556-500
Cyclohexane	Certified ACS	1L	Amber Glass	C556-1
Cyclohexane	HPLC	1L	Amber Glass	C620-1
Cyclohexane	HPLC	1L	Amber Glass/Safe-Cote	C620SK-1

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Cyclohexane	Spectranalyzed	1L	Amber Glass	C555-1
Cyclohexane	Certified ACS	4L	Amber Glass	C556-4
Cyclohexane	HPLC	4L	Amber Glass	C620-4
Cyclohexane	HPLC	4L	Amber Glass/Safe-Cote	C620SK-4
Cyclohexane	Pesticide	4L	Amber Glass	C553-4
Cyclohexane	Reagent	4L	Amber Glass	O2093-4
Cyclohexane	Spectranalyzed	4L	Amber Glass	C555-4
Cyclohexane	Reagent	20L	Steel Pail	O2093-20
Cyclohexanol	Reagent	500mL	Amber Glass	C558-500
Cyclohexanone	Certified	1L	Amber Glass	O2109-1
Cyclohexanone	Certified	4L	Amber Glass	O2109-4
Cyclohexanone	Laboratory	4L	Amber Glass	C550-4
Cyclohexanone	Certified	200L	Stainless Steel Drum	O2109SS200
Cyclohexene (Contains 0.01% tert-Butylcresol as an Inhibitor)	Reagent	1L	Amber Glass	O2111-1
Decane	Certified	500mL	Amber Glass	O2128-500
Diacetone Alcohol	Technical	4L	Amber Glass	D17-4
o-Dichlorobenzene	Certified	1L	Amber Glass	O2231-1
1,2-Dichloroethane	Certified ACS	500mL	Amber Glass	E175-500
1,2-Dichloroethane	Certified ACS	4L	Amber Glass	E175-4
1,2-Dichloroethane	Spectranalyzed	4L	Amber Glass	E190-4
1,2-Dichloroethane	Certified ACS	20L	Steel Pail	E175-20
Diethylenetriamine	Technical	500mL	Amber Glass	D126-500
N,N-Dimethylacetamide	Certified	500mL	Amber Glass	D135-500
Dimethyl Sulfoxide	Certified ACS	500mL	Amber Glass	D128-500
Dimethyl Sulfoxide	Certified ACS	1L	Amber Glass	D128-1
Dimethyl Sulfoxide	Spectranalyzed	1L	Amber Glass	D136-1
Dimethyl Sulfoxide	Certified ACS	4L	Amber Glass	D128-4
Dimethyl Sulfoxide	HPLC	4L	Amber Glass	D159-4
Diethylene Glycol	Reagent	1L	Amber Glass	D49-1
Diisopropylamine	Certified	4L	Amber Glass	O2412-4
1,2-Dimethoxyethane	Certified	1L	Amber Glass	O2430-1
1,2-Dimethoxyethane	Certified	4L	Amber Glass	O2430-4
1,4-Dioxane	Certified ACS	500mL	Amber Glass	D111-500
1,4-Dioxane	Certified ACS	4L	Amber Glass	D111-4
1,4-Dioxane	Histological	4L	SafeTin	D56S-4
1,4-Dioxane	Scintanalyzed	4L	Amber Glass	D116-4
1,4-Dioxane	Scintanalyzed	200L	Steel Drum	D116-200
Ethanol, Anhydrous	Histological	4L	Poly Bottle	A405P-4
Ethanol, Anhydrous	Histological	20L	Steel Pail	A405-20
Ethanol, CDA 19	Histological	1 gal.	Poly Bottle	A406F1GAL
Ethanol, CDA 19	Histological	4L	Poly Bottle	A406P-4
Ethanol, CDA 19	Histological	20L	Steel Pail	A406-20
Ethanolamine	Laboratory	1L	Amber Glass	M251-1
Ethanolamine	Laboratory	4L	Amber Glass	M251-4
Ethyl Acetate	Certified ACS	1L	Amber Glass	E145-1
Ethyl Acetate	HPLC	1L	Amber Glass	E195-1
Ethyl Acetate	Certified ACS	4L	Amber Glass	E145-4
Ethyl Acetate	Certified ACS	4L	SafeTin	E145S-4
Ethyl Acetate	Certified ACS	4L	Amber Glass/Safe-Cote	E145SK-4
Ethyl Acetate	HPLC	4L	Amber Glass	E195-4
Ethyl Acetate	HPLC	4L	Amber Glass/Safe-Cote	E195SK-4
Ethyl Acetate	NF	4L	Amber Glass	E124-4

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Ethyl Acetate	Optima	4L	Amber Glass	E196-4
Ethyl Acetate	HPLC	19L	NOWPak I	E195N1-19
Ethyl Acetate	Optima	4L	Safe-Cote	E196SK-4
Ethyl Acetate	Pesticide	4L	Amber Glass	E191-4
Ethyl Acetate	Spectranalyzed	4L	Amber Glass	E189-4
Ethyl Acetate	Certified ACS	20L	Steel Pail	E145-20
Ethyl Acetate	NF	20L	Steel Pail	E124-20
Ethyl Acetate	Certified ACS	200L	Steel Drum	E145-200
Ethyl Acetoacetate	Reagent	500mL	Amber Glass	E146-500
Ethyl Alcohol, Denatured (Proprietary Solvent)		500mL	Amber Glass	A407-500
Ethyl Alcohol, Denatured (Proprietary Solvent)		1L	Amber Glass	A407-1
Ethyl Alcohol, Denatured (Proprietary Solvent)		4L	Amber Glass	A407-4
Ethyl Alcohol, Denatured (Proprietary Solvent)		4L	Poly Bottle	A407P-4
Ethyl Alcohol, Denatured (Proprietary Solvent)		4L	SafeTin	A407S-4
Ethyl Alcohol, Denatured (Proprietary Solvent)		4L	Amber Glass/Safe-Cote	A407SK-4
Ethyl Alcohol, Denatured (Proprietary Solvent)		20L	Steel Pail	A407-20
Ethyl Alcohol, Denatured (Proprietary Solvent)		200L	Steel Drum	A407-200
Ethylene Glycol	Certified	500mL	Amber Glass	E178-500
Ethylene Glycol	Certified	1L	Amber Glass	E178-1
Ethylene Glycol	Certified	4L	Amber Glass	E178-4
Ethylene Glycol	Laboratory	4L	Amber Glass	E177-4
Ethylene Glycol	Laboratory	20L	Steel Pail	E177-20
Ethylene Glycol	Certified	200L	Steel Drum	E178-200
Ethylene Glycol Monomethyl Ether	Certified	500mL	Amber Glass	E182-500
Ethylene Glycol Monomethyl Ether	Certified	4L	Amber Glass	E182-4
Ethylene Glycol Monomethyl Ether	Certified	20L	Steel Pail	E182-20
Ethylene Glycol Monoethyl Ether	Laboratory	1L	Amber Glass	E180-1
Ethylene Glycol Monoethyl Ether	Laboratory	4L	Amber Glass	E180-4
Ethylene Glycol Monoethyl Ether	Laboratory	20L	Steel Pail	E180-20
Ethylene Glycol Monoethyl Ether Acetate	Laboratory	4L	Amber Glass	E181-4
Ethyl Ether	Pesticide	4L	Amber Glass	E199-4
Ethyl Ether	Spectranalyzed	4L	Tin Can	E197-4
Ethyl Ether	Laboratory	20L	Steel Pail	E134-20
Ethyl Ether, Fat Extraction	Laboratory	4L	Tin Can	E492-4
Ethyl Ether, Fat Extraction	Laboratory	20L	Steel Pail	E492-20
Ethyl Ether, Anhydrous (BHT Stabilized)	Certified ACS	500mL	Aluminum Can	E138500
Ethyl Ether, Anhydrous (BHT Stabilized)	Certified ACS	1L	Aluminum Can	E138-1
Ethyl Ether, Anhydrous (BHT Stabilized)	Certified ACS	4L	Aluminum Can	E138-4
Ethyl Ether, Anhydrous (BHT Stabilized)	Certified ACS	20L	Steel Pail	E138-20
Ethyl Ether, Anhydrous (BHT Stabilized)	HPLC	4L	Tin Can	E198-4
Ethylbenzene	Certified	1L	Amber Glass	O2751-1
Formaldehyde, 40% by volume	USP	4L	Poly Bottle	F77P-4
Formaldehyde, 40% by volume	USP	20L	PolyPac	F77P-20
Formaldehyde, 40% by volume	USP	200L	Steel Drum/Poly Liner	F77-200
Formaldehyde, 37% by weight	Histological	1 gal.	Poly Bottle	F75F1GAL
Formaldehyde, 37% by weight	Histological	1 gal.	Poly Bottle	F75P1GAL
Formaldehyde, 37% by weight	Histological	20L	PolyPac	F75P-20
Formaldehyde, 37% by weight (with preservative)	Certified ACS	500mL	Amber Glass	F79-500
Formaldehyde, 37% by weight (with preservative)	Certified ACS	1L	Amber Glass	F79-1
Formaldehyde, 37% by weight (with preservative)	Certified ACS	4L	Amber Glass	F79-4
Formaldehyde, 37% by weight (with preservative)	Certified ACS	4L	Poly Bottle	F79P-4

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Formaldehyde, 37% by weight (with preservative)	Certified ACS	20L	Rigid Poly Bottle/ Corrugated Box	F79-20
Formaldehyde, 37% by weight (with preservative)	Certified ACS	20L	PolyPac	F79P-20
Formaldehyde, 37% by weight (with preservative)	Certified ACS	200L	Poly Drum	F79-200
Formalin, Buffered, 10% (Acetate Buffer)	Certified	4L	Poly Bottle	SF99-4
Formalin, Buffered, 10% (Acetate Buffer)	Certified	20L	PolyPac	SF99-20
Formalin Solution, 10%	Histological	4L	Poly Bottle	SF98-4
Formalin Solution, 10%	Histological	20L	PolyPac	SF98-20
Formamide	Certified ACS	1L	Amber Glass	F84-1
Formic Acid	OPTIMA LC/MS	50mL	Poly Bottle	A117-50
Formic Acid	OPTIMA LC/MS	10 x 1mL	Ampule	A117- 10X1AMP
Formic Acid	OPTIMA LC/MS	1mL	Ampule	A117-1AMP
Formic Acid	OPTIMA LC/MS	0.5mL	Ampule	A117-05AMP
Formic Acid	OPTIMA LC/MS	2mL	Ampule	A117-2AMP
Glycerol	Certified ACS	500mL	Amber Glass	G33-500
Glycerol	Certified ACS	1L	Amber Glass	G33-1
Glycerol	Spectranalyzed	1L	Amber Glass	G153-1
Glycerol	Certified ACS	4L	Amber Glass	G33-4
Glycerol	Spectranalyzed	4L	Amber Glass	G153-4
Glycerol	Certified ACS	20L	Rigid Poly Bottle/ Corrugated Box	G33-20
Glycerol	Certified ACS	200L	Steel Drum	G33-200
Glycerol	USP/FCC	500mL	Amber Glass	G31-500
Glycerol	USP/FCC	1L	Amber Glass	G31-1
Glycerol	USP/FCC	4L	Amber Glass	G31-4
Glycerol	USP/FCC/EP/BP	4L	Amber Glass	G30-4
Glycerol	USP/FCC	20L	Rigid Poly Bottle	G31-20
Glycerol	USP/FCC/EP/BP	20L	Rigid Poly Bottle/ Corrugated Box	G30-20
Glycerol	USP/FCC	200L	Poly Drum	G31-200
Glycerol	USP/FCC/EP/BP	200L	Steel Drum	G30200
Heptane	Certified	1L	Amber Glass	O3008-1
Heptane	HPLC	1L	Amber Glass	H350-1
Heptane	HPLC	1L	Amber Glass/Safe-Cote	H350SK-1
Heptane	Certified	4L	Amber Glass	O3008-4
Heptane	HPLC	4L	Amber Glass	H350-4
Heptane	HPLC	4L	Amber Glass/Safe-Cote	H350SK-4
Heptane	Spectranalyzed	4L	Amber Glass	H340-4
Heptane	Technical	20L	Steel Pail	H20-20
Heptane	Technical	200L	Steel Drum	H20-200
n-Hexane	GC <i>Resolv</i>	4L	Amber Glass	H307-4
n-Hexane, 95%	Optima	4L	Amber Glass	H306-4
n-Hexane, 95%	Optima	4L	Amber Glass/Safe-Cote	H306SK-4
Hexanes	Certified	500mL	Amber Glass	H291-500
Hexanes	Certified ACS	500mL	Amber Glass	H292-500
Hexanes	Certified ACS	1L	Amber Glass	H292-1
Hexanes	HPLC	1L	Amber Glass	H302-1
Hexanes	HPLC	1L	Amber Glass/Safe-Cote	H302SK-1
Hexanes	Optima	1L	Amber Glass	H303-1
Hexanes	Spectranalyzed	1L	Amber Glass	H334-1
Hexanes	Certified	4L	Amber Glass	H291-4
Hexanes	Certified	4L	SafeTin	H291S-4

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Hexanes	Certified ACS	4L	Amber Glass	H292-4
Hexanes	Certified ACS	4L	Amber Glass/Safe-Cote	H292SK-4
Hexanes	HPLC	4L	Amber Glass	H302-4
Hexanes	HPLC	4L	Amber Glass/Safe-Cote	H302SK-4
Hexanes	Optima	4L	Amber Glass	H303-4
Hexanes	Optima	4L	Amber Glass/Safe-Cote	H303SK-4
Hexanes	Pesticide	4L	Amber Glass	H300-4
Hexanes	Spectranalyzed	4L	Amber Glass	H334-4
Hexanes	Technical	4L	SafeTin	N3S-4
Hexanes	Certified	20L	Steel Pail	H291-20
Hexanes	Certified ACS	20L	Steel Pail	H292-20
Hexanes	Technical	20L	Steel Pail	N3-20
Hexanes	Certified	200L	Steel Drum	H291-200
Hexanes	Certified ACS	200L	Steel Drum	H292-200
Hexanes	Technical	200L	Steel Drum	N3-200
IPA (Isopropyl Alcohol), 70% (v/v)		500mL	Poly Bottle	A459-500
IPA (Isopropyl Alcohol), 70% (v/v)		1L	Poly Bottle	A459-1
IPA (Isopropyl Alcohol), 70% (v/v)		4L	Poly Bottle	A459-4
IPA (Isopropyl Alcohol), 70% (v/v)		20L	PolyPac	A459-20
Isoamyl Alcohol	Certified	500mL	Amber Glass	A393-500
Isoamyl Alcohol	Certified	4L	Amber Glass	A393-4
Isobutyl Alcohol	Certified ACS	1L	Amber Glass	A397-1
Isobutyl Alcohol	Certified ACS	4L	Amber Glass	A397-4
Isooctane	Certified ACS	1L	Amber Glass	O299-1
Isooctane	HPLC	1L	Amber Glass	O296-1
Isooctane	HPLC	1L	Amber Glass/Safe-Cote	O296SK-1
Isooctane	Spectranalyzed	1L	Amber Glass	O300-1
Isooctane	Certified ACS	4L	Amber Glass	O299-4
Isooctane	HPLC	4L	Amber Glass	O296-4
Isooctane	HPLC	4L	Amber Glass/Safe-Cote	O296SK-4
Isooctane	Optima	4L	Amber Glass	O301-4
Isooctane	Pesticide	4L	Amber Glass	O297-4
Isooctane	Spectranalyzed	4L	Amber Glass	O300-4
Isopropanol, 10% (v/v)	Certified	1 gal.	Poly Bottle	A460-1GAL
Isopropyl Acetate	Certified	1L	Poly Bottle	O61111
Isopropyl Acetate	Certified	2.5L	Poly Bottle	O6111-212
Isopropyl Ether	Certified	4L	Amber Glass	E141-4
Kerosene (Odorless)		4L	Amber Glass	K10-4
Kerosene (Odorless)		200L	Steel Drum	K10-200
Ligroine, 100% Heptane	Laboratory	4L	Amber Glass	O3387-4
Ligroine, Hexane	Laboratory	20L	Steel Pail	O3386-20
Methanol	Certified ACS	500mL	Amber Glass	A412-500
Methanol	NF	500mL	Amber Glass	A413500
Methanol	Optima LC/MS	500mL	Amber Glass	A456-500
Methanol	Certified ACS	1L	Amber Glass	A412-1
Methanol	HPLC	1L	Amber Glass	A452-1
Methanol	HPLC	1L	Amber Glass/Safe-Cote	A452SK-1
Methanol	Optima	1L	Amber Glass	A454-1
Methanol	Optima LC/MS	1L	Amber Glass	A456-1
Methanol	Spectranalyzed	1L	Amber Glass	A408-1
Methanol	Histological	1 gal.	Poly Bottle	A433F1GAL
Methanol	Certified ACS	4L	Amber Glass	A412-4
Methanol	Certified ACS	4L	Poly Bottle	A412P-4

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Methanol	Certified ACS	4L	Amber Glass/Safe-Cote	A412SK-4
Methanol	Electronic	4L	Poly Bottle	A947-4
Methanol	GC <i>Resolv</i>	4L	Amber Glass	A457-4
Methanol	Histological	4L	Poly Bottle	A433P-4
Methanol	Histological	4L	SafeTin	A433S-4
Methanol	HPLC	4L	Amber Glass	A452-4
Methanol	HPLC	4L	Amber Glass/Safe-Cote	A452SK-4
Methanol	Laboratory	4L	Amber Glass	A411-4
Methanol	NF	4L	Amber Glass	A4134
Methanol	Optima	4L	Amber Glass	A454-4
Methanol	Optima	4L	Amber Glass/Safe-Cote	A454SK-4
Methanol	Optima LC/MS	4L	Amber Glass	A456-4
Methanol	Pesticide	4L	Amber Glass	A450-4
Methanol	Spectranalyzed	4L	Amber Glass	A408-4
Methanol	Spectranalyzed	4L	Amber Glass/Safe-Cote	A408SK-4
Methanol	HPLC	19L	NOWPak I	A452N1-19
Methanol	Certified ACS	20L	Unlined Steel Pail	A412-20
Methanol	Certified ACS	20L	Unichrome Steel Pail	A434-20
Methanol	Histological	20L	Steel Pail	A433S-20
Methanol	Laboratory	20L	Steel Pail	A411-20
Methanol	NF	20L	Steel Pail	A41320
Methanol	Certified ACS	200L	Steel Drum	A412-200
Methanol	Histological	200L	Steel Drum	A433S-200
Methanol	NF	200L	Steel Drum	A413-200
Methanol (Low Water)	Low Water	4L	Amber Glass	A935-4
Methanol		500mL	Amber Glass	A453-500
(Purge and Trap Suitable for Volatile Organic Residue Analysis)				
Methanol		1L	Amber Glass	A453-1
(Purge and Trap Suitable for Volatile Organic Residue Analysis)				
Methanol	HPLC	19L	NOWPak I	A452N1-19
Methylal	Certified	500mL	Amber Glass	M222-500
Methylene Chloride	Optima	1L	Amber Glass	D151-1
Methylene Chloride	GC <i>Resolv</i>	4L	Amber Glass	D154-4
Methylene Chloride	Optima	4L	Amber Glass	D151-4
Methylene Chloride	Pesticide	4L	Amber Glass	D142-4
Methylene Chloride (Not Stabilized)	HPLC	1L	Amber Glass	D150-1
Methylene Chloride (Not Stabilized)	HPLC	1L	Amber Glass/Safe-Cote	D150SK-1
Methylene Chloride (Not Stabilized)	HPLC	4L	Amber Glass	D150-4
Methylene Chloride (Not Stabilized)	HPLC	4L	Amber Glass/Safe-Cote	D150SK-4
Methylene Chloride (Stabilized)	Certified ACS	500mL	Amber Glass	D37-500
Methylene Chloride (Stabilized)	Certified ACS	1L	Amber Glass	D37-1
Methylene Chloride (Stabilized)	Certified ACS	4L	Amber Glass	D37-4
Methylene Chloride (Stabilized)	Certified ACS	4L	Amber Glass/Safe-Cote	D37SK-4
Methylene Chloride (Stabilized)	Certified ACS	20L	Steel Pail	D37-20
Methylene Chloride (Stabilized)	Certified ACS	200L	Steel Drum	D37-200
Methylene Chloride (Stabilized)	HPLC	1L	Amber Glass	D143-1
Methylene Chloride (Stabilized)	HPLC	1L	Amber Glass/Safe-Cote	D143SK-1
Methylene Chloride (Stabilized)	HPLC	4L	Amber Glass	D143-4
Methylene Chloride (Stabilized)	HPLC	4L	Amber Glass/Safe-Cote	D143SK-4
Methylene Chloride (Stabilized)	Spectranalyzed	1L	Amber Glass	D35-1
Methylene Chloride (Stabilized)	Spectranalyzed	4L	Amber Glass	D35-4
Methylene Chloride (with Cyclohexene Preservative)	HPLC	1L	Amber Glass	D138-1
Methylene Chloride (with Cyclohexene Preservative)	HPLC	4L	Amber Glass	D138-4

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Methylene Chloride (with Cyclohexene Preservative)	HPLC	4L	Amber Glass/Safe-Cote	D138SK-4
Methyl Acetate	Certified	500mL	Amber Glass	M203-500
Methyl Acetate	Certified	4L	Amber Glass	M203-4
Methyl Ethyl Ketone	Certified ACS	1L	Amber Glass	M209-1
Methyl Ethyl Ketone	Technical	1L	Amber Glass	M208-1
Methyl Ethyl Ketone	Certified ACS	4L	Amber Glass	M209-4
Methyl Ethyl Ketone	Certified ACS	4L	SafeTin	M209S-4
Methyl Ethyl Ketone	Technical	4L	Amber Glass	M208-4
Methyl Ethyl Ketone	Certified ACS	20L	Steel Pail	M209-20
Methyl Ethyl Ketone	Technical	20L	Steel Pail	M208-20
Methyl Ethyl Ketone	Certified ACS	200L	Steel Drum	M209-200
2-Methylbutane	Certified	4L	Amber Glass	O3551-4
Methyl iso-Butyl Ketone	Certified ACS	1L	Amber Glass	M213-1
Methyl iso-Butyl Ketone	Certified ACS	4L	Amber Glass	M213-4
Methyl iso-Butyl Ketone	Certified ACS	20L	Steel Pail	M213-20
Methyl iso-Butyl Ketone	Certified ACS	200L	Steel Drum	M213-200
Methyl tert-Butyl Ether	HPLC	4L	Amber Glass	E127-4
Morpholine	Certified ACS	1L	Amber Glass	M263-1
N,N-Dimethylformamide	Certified ACS	500mL	Amber Glass	D119-500
N,N-Dimethylformamide	Certified ACS	1L	Amber Glass	D119-1
N,N-Dimethylformamide	Spectranalyzed	1L	Amber Glass	D131-1
N,N-Dimethylformamide	Certified ACS	4L	Amber Glass	D119-4
N,N-Dimethylformamide	Certified ACS	4L	SafeTin	D119S-4
N,N-Dimethylformamide	Spectranalyzed	4L	Amber Glass	D131-4
N,N-Dimethylformamide	Certified ACS	20L	Poly Pail	D119-20
Nitrobenzene	Certified ACS	500mL	Amber Glass/Poison Pack	N911-500
Nitrobenzene	Certified ACS	4L	Amber Glass/Poison Pack	N911-4
Nitromethane	Certified ACS	500mL	Amber Glass	N98-500
Octane	Reagent	1L	Amber Glass	O3980-1
1-Octanol	Certified	500mL	Amber Glass	A402-500
1-Octanol	Certified	4L	Amber Glass	A402-4
2-Octanol	Laboratory	500mL	Amber Glass	O269-500
Pentane	HPLC	1L	Amber Glass	P399-1
Pentane	HPLC	1L	Amber Glass/Safe-Cote	P399SK-1
Pentane	Spectranalyzed	1L	Amber Glass	P393-1
Pentane	Certified	4L	Amber Glass	O4062-4
Pentane	HPLC	4L	Amber Glass	P399-4
Pentane	HPLC	4L	Amber Glass/Safe-Cote	P399SK-4
Pentane	Pesticide	4L	Amber Glass	P400-4
Pentane	HPLC	19L	Steel Drum	P399RS19
Pentane	Certified	20L	Steel Pail	O4062-20
Petroleum Ether	Certified ACS	500mL	Amber Glass	E139-500
Petroleum Ether	Certified ACS	1L	Amber Glass	E139-1
Petroleum Ether	Certified ACS	4L	Amber Glass	E139-4
Petroleum Ether	Certified ACS	4L	SafeTin	E139S-4
Petroleum Ether	Certified ACS	4L	Amber Glass/Safe-Cote	E139SK-4
Petroleum Ether	Optima	4L	Amber Glass	E120-4
Petroleum Ether	Optima	4L	Amber Glass/Safe-Cote	E120SK-4
Petroleum Ether	Pesticide	4L	Amber Glass	P480-4
Petroleum Ether	Certified ACS	20L	Steel Pail	E139-20

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Petroleum Ether	Certified ACS	200L	Steel Drum	E139-200
1-Propanol	Certified	500mL	Amber Glass	A414-500
1-Propanol	Certified	1L	Amber Glass	A414-1
1-Propanol	Certified	4L	Amber Glass	A414-4
1-Propanol	Certified	4L	SafeTin	A414S-4
1-Propanol	Certified	20L	Steel Pail	A414-20
2-Propanol	Certified ACS Plus	500mL	Amber Glass	A416-500
2-Propanol	Optima LC/MS	500mL	Amber Glass	A461-500
2-Propanol	USP	500mL	Amber Glass	A516500
2-Propanol	Certified	1L	Amber Glass	A417-1
2-Propanol	HPLC	1L	Amber Glass	A451-1
2-Propanol	Spectranalyzed	1L	Amber Glass	A419-1
2-Propanol	Histological	1 gal.	Poly Bottle	A426F1GAL
2-Propanol	Certified	4L	Amber Glass	A417-4
2-Propanol	Certified ACS Plus	4L	Amber Glass	A416-4
2-Propanol	Certified ACS Plus	4L	Poly Bottle	A416P-4
2-Propanol	Certified ACS Plus	4L	SafeTin	A416S-4
2-Propanol	Certified ACS Plus	4L	Amber Glass/Safe-Cote	A416SK-4
2-Propanol	Histological	4L	Poly Bottle	A426P-4
2-Propanol	Histological	4L	SafeTin	A426S-4
2-Propanol	HPLC	4L	Amber Glass	A451-4
2-Propanol	HPLC	4L	Amber Glass/Safe-Cote	A451SK-4
2-Propanol	Laboratory	4L	Amber Glass	A415-4
2-Propanol	Optima	4L	Amber Glass	A464-4
2-Propanol	Optima	4L	Amber Glass/Safe-Cote	A464SK-4
2-Propanol	Pesticide	4L	Amber Glass	A519-4
2-Propanol	Spectranalyzed	4L	Amber Glass	A419-4
2-Propanol	USP	4L	Amber Glass	A516-4
2-Propanol	Certified ACS Plus	20L	Steel Pail	A416-20
2-Propanol	Histological	20L	Steel Pail	A426S-20
2-Propanol	Laboratory	20L	Steel Pail	A415-20
2-Propanol	USP	20L	Steel Pail	A516-20
2-Propanol	Certified ACS Plus	200L	Steel Drum	A416-200
2-Propanol	Histological	200L	Steel Drum	A426S-200
2-Propanol	USP	200L	Steel Drum	A516-200
2-Propanol (Low Water)	Low Water	4L	Amber Glass	A520-4
Propylene Glycol	USP/FCC	1L	Amber Glass	P355-1
Propylene Glycol	USP/FCC	4L	Amber Glass	P355-4
Propylene Glycol	USP/FCC	20L	Steel Pail	P355-20
Propylene Glycol	USP/FCC	200L	Steel Drum	P355-200
Propylene Oxide	Reagent	1L	Amber Glass	O4332-1
p-Xylene	Certified	4L	Amber Glass	O5082-4
Pyridine	Certified ACS	500mL	Amber Glass	P368-500
Pyridine	Certified ACS	1L	Amber Glass	P368-1
Pyridine	Certified ACS	4L	Amber Glass	P368-4
Pyridine	Certified ACS	4L	SafeTin	P368S-4
Pyridine	Technical	4L	Amber Glass	P369-4
sec-Butanol	Certified	4L	Amber Glass	O1664-4
Stoddard Solvent		4L	Amber Glass	S457-4
tert-Butanol	Certified	1L	Poly Bottle	A401-1
1,1,2,2-Tetrabromoethane	Certified	100mL	Amber Glass	A29-100
1,1,2,2-Tetrabromoethane	Laboratory	4L	Amber Glass	A33-4

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Tetrachloroethylene	Technical	4L	Amber Glass	C182-4
Tetrachloroethylene	Technical	20L	Steel Pail	C182-20
Tetrachloroethylene (Stabilized)	Reagent	4L	Amber Glass	O4586-4
Tetrahydrofuran (BHT as preservative)	Certified	500mL	Amber Glass	T397-500
Tetrahydrofuran (BHT as preservative)	Histological	500mL	Amber Glass	T400-500
Tetrahydrofuran (BHT as preservative)	Certified	1L	Amber Glass	T397-1
Tetrahydrofuran (no preservative)	HPLC	1L	Amber Glass	T425-1
Tetrahydrofuran (no preservative)	HPLC	1L	Amber Glass/Safe-Cote	T425SK-1
Tetrahydrofuran (no preservative)	Optima	1L	Amber Glass	T427-1
Tetrahydrofuran (BHT as preservative)	Certified	4L	Amber Glass	T397-4
Tetrahydrofuran (BHT as preservative)	Certified	4L	Amber Glass/Safe-Cote	T397SK-4
Tetrahydrofuran (BHT as preservative)	Histological	4L	Amber Glass	T400-4
Tetrahydrofuran (no preservative)	HPLC	4L	Amber Glass	T425-4
Tetrahydrofuran (no preservative)	HPLC	4L	Amber Glass/Safe-Cote	T425SK-4
Tetrahydrofuran (no preservative)	Optima	4L	Amber Glass	T427-4
Tetrahydrofuran (no preservative)	Optima	4L	Amber Glass/Safe-Cote	T427SK-4
Tetrahydrofuran (no preservative)	Spectranalyzed	4L	Amber Glass	T424-4
Tetrahydrofuran (BHT as preservative)	Certified	20L	Steel Pail	T397-20
Tetrahydrofuran (BHT as preservative)	Histological	20L	Steel Pail	T400-20
Tetrahydrofuran (BHT as preservative)	Certified	200L	Steel Drum	T397-200
Toluene	Certified ACS	500mL	Amber Glass	T324-500
Toluene	Certified ACS	1L	Amber Glass	T324-1
Toluene	HPLC	1L	Amber Glass	T290-1
Toluene	HPLC	1L	Amber Glass/Safe-Cote	T290SK-1
Toluene	Histological	1 gal.	Poly Bottle	T326F1GAL
Toluene	Certified ACS	4L	Amber Glass	T324-4
Toluene	Certified ACS	4L	SafeTin	T324S-4
Toluene	Certified ACS	4L	Amber Glass/Safe-Cote	T324SK-4
Toluene	Histological	4L	Poly Bottle	T326P-4
Toluene	HPLC	4L	Amber Glass/Safe-Cote	T290SK-4
Toluene	Laboratory	4L	Amber Glass	T323-4
Toluene	Optima	4L	Amber Glass	T291-4
Toluene	Scintanalyzed	4L	Amber Glass	T313-4
Toluene	Scintanalyzed	4L	Amber Glass/Safe-Cote	T313SK-4
Toluene	Spectranalyzed	4L	Amber Glass	T330-4
Toluene	Certified ACS	20L	Steel Pail	T324-20
Toluene	Histological	20L	Steel Pail	T326S20
Toluene	Laboratory	20L	Steel Pail	T323-20
Toluene	HPLC	4L	Amber Glass	T290-4
1,2,4-Trichlorobenzene	HPLC	4L	Amber Glass	O4846-4
Trichloroethylene	Electronic	4L	Amber Glass	T403-4
Trichloroethylene	Technical	4L	Amber Glass	T340-4
Trichloroethylene (Stabilized)	Certified ACS	500mL	Amber Glass	T341-500
Trichloroethylene (Stabilized)	Certified ACS	20L	Steel Pail	T341-20
Triethanolamine	Certified	500mL	Amber Glass	T407-500
Triethanolamine	NF	500mL	Amber Glass	T350-500
Triethanolamine	Certified	1L	Amber Glass	T407-1
Triethanolamine	Certified	4L	Amber Glass	T407-4
Triethanolamine	NF	4L	Amber Glass	T350-4
Triethylamine	HPLC	100mL	Amber Glass	O4884-100
Triethylamine	Reagent	1L	Amber Glass	O4885-1

Solvents (Continued)

Solvent	Purity Grade	Quantity	Packaging	Cat. No.
Trifluoroacetic Acid	Optima LC/MS	50mL	Poly Bottle	A116-50
Trifluoroacetic Acid	Optima LC/MS	10 x 1mL	Ampule	A116-10X1AMP
Trifluoroacetic Acid	Optima LC/MS	1mL	Ampule	A116-1AMP
Trifluoroacetic Acid	Optima LC/MS	0.5mL	Ampule	A116-05AMP
Trifluoroacetic Acid	Optima LC/MS	2mL	Ampule	A116-2AMP
Triethylamine	Reagent	4L	Amber Glass	O4885-4
Triethylamine	Reagent	20L	Steel Pail	O488520
Triethylene Glycol	Laboratory	4L	Amber Glass	T346-4
Triethylenetetramine	Technical	500mL	Amber Glass	T410-500
Triethylenetetramine	Technical	1L	Amber Glass	T410-1
Water	Optima LC/MS	500mL	Amber Glass	W6-500
Water	Plasma	500mL	LDPE Bottle	W9-500
Water	HPLC	1L	Amber Glass	W5-1
Water	HPLC	1L	Amber Glass/Safe-Cote	W5SK-1
Water	Environment	1L	Amber Glass	W11-1
Water	Optima LC/MS	1L	Amber Glass	W6-1
Water	Plasma	1L	LDPE Bottle	W9-1
Water	Plasma	2L	LDPE Bottle	W9-2
Water	HPLC	4L	Amber Glass	W5-4
Water	HPLC	4L	Amber Glass/Safe-Cote	W5SK-4
Water	Environment	4L	Amber Glass	W11-4
Water	Optima	4L	Amber Glass	W7-4
Water	Optima LC/MS	4L	Amber Glass	W6-4
Water	HPLC	19L	NOWPak I	W5-N119
Water, DIUF		4L	Poly Bottle	W2-4
Water, DIUF		20L	PolyPac	W2-20
0.1% FA in Water	OPTIMA LC/MS	500mL	Amber Glass	LS118-500
0.1% FA in Water	OPTIMA LC/MS	1L	Amber Glass	LS118-1
0.1% FA in Water	OPTIMA LC/MS	2.5L	Amber Glass	LS118-212
0.1% FA in Water	OPTIMA LC/MS	4L	Amber Glass	LS118-4
0.1% TFA in Water	OPTIMA LC/MS	55mL	Amber Glass	LS119-500
0.1% TFA in Water	OPTIMA LC/MS	1L	Amber Glass	LS119-1
0.1% TFA in Water	OPTIMA LC/MS	2.5L	Amber Glass	LS119-212
0.1% TFA in Water	OPTIMA LC/MS	4L	Amber Glass	LS119-4
Xylene	Histological	1 gal.	Poly Bottle	X3P-1GAL
Xylene	Certified ACS	4L	Amber Glass	X5-4
Xylene	Certified ACS	4L	SafeTin	X5S-4
Xylene	Certified ACS	4L	Amber Glass/Safe-Cote	X5SK-4
Xylene	Histological	4L	SafeTin	X3S-4
Xylene	Laboratory	4L	Amber Glass	X4-4
Xylene	Certified ACS	20L	Steel Pail	X5-20
Xylene	Histological	20L	Steel Pail	X3S-20
Xylene	Laboratory	20L	Steel Pail	X4-20
Xylene	Certified ACS	200L	Steel Drum	X5-200
Xylene	Histological	200L	Steel Drum	X3S-200
o-Xylene	Certified	4L	Amber Glass	O5081-4
p-Xylene	Certified	500mL	Amber Glass	O5082-500

AMERICAS**Canada**

Fisher Scientific Canada
112 Colonnade Road
Ottawa, Ontario
Post Code: K2E 7L6
Toll-Free Number: 800-234-7437
Fax: 800-463-2996
www.fishersci.ca

Latin America

Fisher Scientific Global Export, Latin America
3970 Johns Creek Court
Suite 500
Suwanee, GA
Post Code: 30024
Toll-Free Number: 770-871-4725
Fax: 770-871-4726
www.fishersci.com

United States

Fisher HealthCare
9999 Veterans Memorial Drive
Houston, TX
Post Code: 77038
Toll-Free Number: 800-640-0640
Fax: 800-290-0290
www.fisherhealthcare.com

Fisher Scientific
2000 Park Lane Drive
Pittsburgh, PA
Post Code: 15275
Toll-Free Number: 800-766-7000
Fax: 800-926-1166
www.fishersci.com

ASIA**China**

Fisher Scientific China
Toll-Free Number: 400 881 5117
sales.china@thermofisher.com
www.fishersci.com.cn

Shanghai Corporate Office
6/F Long Life Mansion
No. 1566 Yan An West Rd.
Shanghai, China
Post Code: 200052
Tel: (8621) 5258 1100
Fax: (8621) 5258 0119

Beijing Office
Units 702-715, 7th Floor
Tower West, Yonghe Plaza
No. 28 Andingmen East Street
Beijing, China
Post Code: 100007
Tel: (8610) 8419 3588
Fax: (8610) 8419 3580

Guangzhou Office
Room 2405-2406, JianLiBao Mansion
No. 410-412 Middle DongFeng Rd.
Guangzhou, China
Post Code: 510030
Tel: (8620) 8314 5288
Fax: (8620) 3877 1941

India

Fisher Scientific India
403-404, B-Wing, Delphi
Hiranandani Business Park
Powai, Mumbai - 400 076
Tel: 1 800 209 7001
Fax: +91 22 6680 3001/3002
qfc.customer@thermofisher.com
http://www.fishersci.in

Japan

Fisher Scientific Japan
Thermo Fisher Scientific K.K.
C-2F, 3-9 Moriya-cho
Kanawaga-ku, Yokohama
221-0022 Japan
Tel: 81 45 450 6310
Fax: 81 45 450 6316
support@fishersci.co.jp
www.fishersci.co.jp

Korea

Fisher Scientific Korea
Sambu Bldg. 13
676 Yeoksam-dong,
Kangnam-Gu
Seoul 135-979, Korea
Dir (02) 527-0300
Customer Service Center: (02) 527-0300
Fax: (02) 527-0311
sales@fishersci.co.kr
www.acros.co.kr

Malaysia

Fisher Scientific Malaysia Sdn Bhd
No.3 Jalan Sepadu 25/123
Taman Perindustrian Axis Seksyen 25
40400 Shah Alam
Selangor Oarul Ehsan Malaysia
Technical Service Hotline: 1-300-88-7868
Tel: (603) 51218888
Fax: (603) 51218899
enquiry.my@thermofisher.com
www.fishersci.com.my

Singapore

Fisher Scientific Singapore
Fisher Scientific Pte Ltd.
8 Pandan Crescent
LL4, #05-05 UE Tech Park
Singapore 128464
Tel: (65) 6873 6006
Fax: (65) 6873 5005
enquiry.sg@thermofisher.com
www.fishersci.com.sg

EUROPE**Austria**

Fisher Scientific (Austria) GmbH
Rudolf von Alt-Platz 1
A-1030 Wien
Phone: 0800 20 88 40
Fax: 0800 20 66 90
info.austria@thermofisher.com
www.at.fishersci.com

Belgium

Fisher Scientific
BP 567
B-7500 Tournai 1
Tel: 056 260 260
Fax: 056 260 270
be.fisher@thermofisher.com
www.be.fishersci.com

Czech Republic

Fisher Scientific, spol. s r.o.
Kosmonautu 324
Pardubice
CZ-530 09
Tel: 466 798 230
Fax: 466 435 008
info.cz@thermofisher.com
www.thermofisher.cz

Denmark

Fisher Scientific Biotech Line A/S
Industrivej 3
Postboks 60
OK-3550 Slangerup
Tel: +45 70 27 99 20
Fax: +45 70 27 99 29
kundeservice@thermofisher.com
www.fishersci.dk

Finland

Fisher Scientific Oy
Ratastie 2
FI-01620 Vantaa
Tel: +358 802 76 237
Fax: +358 802 76 235
fisher.fi@thermofisher.com
www.fishersci.fi

France

Fisher Scientific
Parc d'Innovation BP 50111
67403 Illkirch Cedex
Tel: 03 88 67 53 20
Fax: 03 88 67 11 68
fr.commande@thermofisher.com
www.fr.fishersci.com

Germany

Fisher Scientific GmbH
Im Heiligen Feld 17
D-58239 Schwerte
Phone: 0800 3 47 43 70
Fax: 0800 3 47 43 71
info.germany@thermofisher.com
www.de.fishersci.com

Ireland

Fisher Scientific Ireland
Suite 3 Plaza 212
Blanchardstown Corporate
Park 2
Ballycoolin
Dublin 15
Tel: +353 01 885 5854
Fax: +353 01 899 1855
fsie.sales@thermofisher.com
www.ie.fishersci.com

Italy

Fisher Scientific
Tel: 02 953 28 258
fax: 02 953 27 374
it.fisher@thermofisher.com
www.it.fishersci.com

The Netherlands

Fisher Scientific
Postbus 4
1120 AA Landsmeer
Tel: 020 487 70 00
Fax: 020 487 70 70
nl.info@thermofisher.com
www.nl.fishersci.com

Norway

Fisher Scientific
Frysjaveien 33E
0884 Oslo
Tel: +47 22 95 59 59
Fax: +47 22 95 59 40
fisher.no@thermofisher.com
www.fishersci.no

Portugal

Fisher Scientific
Rua Pedro Álvares Cabral, n°24, 3°D
Edifício Euro - Infatado
2670-391 Loures
Telephone: +351 21 425 33 50/4
Fax: +351 21 425 33 51
pt.fisher@thermofisher.com
www.pt.fishersci.com

Spain

Fisher Scientific
C/ Luis I, 9
28031 MADRID
Tfno: 91 380 67 10
Fax: 91 380 85 02
es.fisher@thermofisher.com
www.es.fishersci.com

Sweden

Fisher Scientific
Box 9193
40094 Göteborg
Tel: +46 31 - 68 94 30
Fax: +46 31 - 68 07 17
gtf.info@thermofisher.com
www.fishersci.se

Switzerland

Fisher Scientific
Wilstrasse 57 - Postfach 1006
5610 Wohlen
Tel: 056 618 41 11
Fax: 056 618 41 41
info.ch@thermofisher.com
www.ch.fishersci.com

United Kingdom

Fisher Scientific UK
Bishop Meadow Road
Loughborough
Leicestershire LE11 5RG
Tel: +44 (0)1509 231166
Fax: +44 (0)1509 231893
info@fisher.co.uk
www.fisher.co.uk

Rest of Europe

Thermo Fisher Scientific
Geel West Zone 2
Janssen Pharmaceuticaalaa 3a
2440 Geel - Belgium
Tel: +32 14 57 52 11
fax: +32 14 59 26 10
www.acros.com

MIDDLE EAST AND AFRICA**Fisher Scientific Global Export,****Latin America**

3970 Johns Creek Court
Suite 500
Suwanee, GA
Post Code: 30024
Toll-Free Number: 770-871-4725
Fax: 770-871-4726
www.fishersci.com

OCEANIA**Australia**

Thermo Fisher Scientific Australia Pty Ltd
5 Caribbean Drive
Scoresby, VIC 3179
Tel: +1300 735 292
Fax: +61 3 9763 1169
AUinfo@thermofisher.com
www.thermofisher.com.au

New Zealand

Thermo Fisher Scientific New Zealand Ltd 244
Bush Road,
Albany, Auckland 0632
Tel: 0800 933 966
Fax: +64 9 980 6788
infonz@thermofisher.com
www.thermofisher.co.nz

©2009 Thermo Fisher Scientific Inc.

All rights reserved. Litho in U.S.A.

All brands and trademarks are part of
Thermo Fisher Scientific Inc. and its
subsidiaries

10_3047 KE/LY 10M-IW-11/11
BN1221095