

ChemACX Vendor List

A B C D E F G H I J K L M N O P Q R S T U V W X Y Z
-#-

1K SCIENTIFIC

3A PHARMATECH

3B SCIENTIFIC CORP.

Updated May 2020

3RCHEM

8S LABORATORIES LLC

-A-

**A&A LIFE SCIENCE INC. DBA
AVACHEM SCIENTIFICS**

A&C PHARMTECH, INC

A&J PHARMATECH

A.G. SCIENTIFIC

A1 BIOCHEM LABS

Updated Dec. 2020

A2B CHEM LLC

Updated Apr. 2021

AA BLOCKS LLC

Updated Jul. 2020

AB CHEM TECHNOLOGIES LLC

ABCR GMBH & Co. KG

ABSCO LTD

**ABX ADVANCED BIOCHEMICAL
COMPOUNDS**

AEICHEM

AK SCIENTIFIC, INC.

AKL RESEARCH LLP

AOBICHEM USA

AOKBIO Co., LTD

AOKCHEM Co., LTD.

APAC PHARMACEUTICAL

AQCHEM LLC

ASW MEDCHEM, INC.

AURUM PHARMATECH LLC

AARON CHEMICALS LLC

New Feb. 2021

AARON CHEMISTRY

ABAMACHEM LTD

ABBEY COLOR

**ABBOTT LABORATORIES CHEMICAL
DIVISION**

ABBY PHARMATECH LLC

ABERJONA LABORATORIES

ABOVICHEM LLC

ABSOLUTE CHIRAL

ACADECHEM COMPANY LIMITED

ACADEMIQUE PHYTOCANADA

ACCELA CHEMBIO Co., LTD.

Updated Aug. 2020

ACCELEDEV CHEMICAL LLC

ACCUSTANDARD, INC.

ACE SYNTHESIS LLC

ACENTEX SCIENTIFIC, INC.

ACESCHEM INC.

ACESYS PHARMATECH

ACETO CORPORATION

ACHEMICA

ACINOPEPTIDE Co. LTD

ACME BIOSCIENCE, INC.

ACROS ORGANICS

ACROS ORGANICS – USA

ACTIVATE SCIENTIFIC GMBH

ADESIS INC.

ADVAMACS

ADVANCE RESEARCH CHEMICALS, INC.

ADVANCED AROMATICS, L.P.

ADVANCED ASYMMETRICS, INC.

ADVANCED CHEMBLOCKS INC.

ADVANCED CHEMTECH

**ADVANCED MOLECULAR
TECHNOLOGIES PTY LTD**

**ADVANCED SCIENTIFIC
INTERNATIONAL, LLC**

**ADVANCED SYNTHESIS
TECHNOLOGIES, S.A.**

**ADVANCED TECHNOLOGY &
INDUSTRIAL Co., LTD.**

AIMS FINE CHEMICALS INC.

AIR PRODUCTS & CHEMICALS, INC.

**AJINOMOTO - AMINO ACID
DEPARTMENT**

ALBANY MOLECULAR RESEARCH, INC.

ALBEMARLE CORPORATION

ALBRIGHT & WILSON AMERICAS

ALCHEM INTERNATIONAL LIMITED

ALCHEM PHARMTECH, INC.

ALDLAB CHEMICALS, LLC

ALDRICH

ALFA AESAR

ALFA CHEMISTRY

Updated Jul. 2020

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

ChemACX Vendor List

ALFA CHEMISTRY (MATERIALS)

New Aug. 2020

ALFA PYRIDINES

ALICHEM INC.

Updated Feb. 2021

ALLIEDSIGNAL INC. SPECIALTY CHEMICALS

AMBEED, INC.

Updated Feb. 2021

AMERICAN ADVANCED ORGANICS (DIV OF ALBANY MOLECULAR)

AMERICAN ADVANCED SCIENTIFIC INC.

AMERICAN BIOANALYTICAL

AMERICAN BIOCHEMICALS

AMERICAN CUSTOM CHEMICAL CORP

AMERICAN PEPTIDE COMPANY

AMERICAN RADIOLABELED CHEMICALS, INC.

AMERICHEM, INC.

AMRESKO INC.

ANGEL PHARMATECH, LTD.

ANGELINI FINE CHEMICALS

ANGENE INTERNATIONAL LIMITED

ANGUS CHEMICAL COMPANY

ANNKER ORGANICS

ANTIBIOTICOS SPA

ANWARD CHEMICALS

APEIRON SYNTHESIS

APEX SCIENTIFIC INC.

APEXMOL TECHNOLOGY

APITOXIN

APOLLO SCIENTIFIC LTD. (BIOCHEMICALS)

APOLLO SCIENTIFIC LTD. (FLUORO CHEMICALS)

APOLLO SCIENTIFIC LTD. (INTERMEDIATES)

APOLLO SCIENTIFIC LTD. (SUPPORT RESEARCH LIBRARY)

APPLICHEM

ARBOŃOVA OY AB

ARCTOM CHEMICALS

ARK PHARM, INC.

ARKÉ ORGANICS SRL

AROMA CIRCLE

ARONIS

ARRAN CHEMICAL CO.

ARYLCHEM CO.

ARZA BIOSCIENCE LIMITED

Updated Sep. 2020

ASAHI GLASS COMPANY - FLUORO CHEMICALS DIVISION

ASCENT SCIENTIFIC

ASHWOOD CHEMICALS LTD.

ASINEX LTD.

ASINEX LTD. PLATINUM COLLECTION

ASISCHEM INC.

ASTATECH, INC.

ASTAR PHARMACEUTICALS, LLC

ATOMOLE SCIENTIFIC Co., LTD.

AUGUSTUS OILS LTD

AURORA FINE CHEMICALS LTD

Updated Feb. 2021

AVANTI (A SIGMA-ALDRICH BRAND)

AVIABOR FINE BORON COMPOUNDS

AVONCHEM/CHROMOS EXPRESS LTD

AXISPHARM

New Aug. 2020

AXITEC Co., LTD

AZEPINE LTD

-B-

BACHEM AG

BASF CHEMICAL INTERMEDIATES

BE-ASSOCIATES, LLC

BLD PHARMATECH LTD

Updated Jan. 2021

BOC SCIENCES

BSC CHEMICALS

BAIHUA BIO-PHARMACEUTICAL CO. LTD.

BALAJI AMINES LTD

BAYER CORPORATION

BAYER INORGANICS

BEDOUKIAN RESEARCH, INC.

BEIJING CENTURY RICHAP CHEMISTRY CO. LTD

BEIJING GOLDEN OLIVE COMPANY

BEPHARM LTD.

BESTIPHARMA

BETA PHARMA, INC.

BEWE LEEPHARMA INC.

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

ChemACX Vendor List

BIDDLE SAWYER CORPORATION

BIDE PHARMATECH

BIO FARMA LTD.

BIO-RESEARCH PRODUCTS, INC.

BIOBLOCKS, INC.

BIOQUADRANT INC.

BIOVISION, INC.

BIOCHEMIE GMBH

BIOCHEMPEG

BIOCORE PHARMTECH LIMITED

BIOFINE INTERNATIONAL INC.

Updated Feb. 2021

BIOMOL

BIOSYNTH AG

BIOTIUM, INC.

BIOTREND CHEMIKALIEN GMBH
(NEUROCHEMICALS)

BIOTREND CHEMIKALIEN GMBH
(RADIOCHEMICALS)

BOERCHEM PHARMTECH CO., LTD

BOROCHEM

BORON MOLECULAR

BOSCHE SCIENTIFIC LLC

BOULDER SCIENTIFIC COMPANY

BRIDGE ORGANICS CO.

BRIGHT EVERGREEN

BROADPHARM

BURDICK & JACKSON, INC.

BUSH BOAKE ALLEN INC.

BUTT PARK LTD

-C-

C/D/N ISOTOPES INC.

CF PLUS CHEMICALS S.R.O.

CGENETECH, INC.

CMS CHEMICALS LTD --
CARBOHYDRATES

CMS CHEMICALS LTD -- ORGANICS

CNH TECHNOLOGIES, INC.

CU CHEMIE UETIKON

CACHESYN INC.

CALAIRE CHIMIE S.A.

CALLERY CHEMICAL COMPANY

CAMBREX CORPORATION

CAMBRIDGE ISOTOPE LABS

CAMBRIDGE MAJOR LABORATORIES,
INC.

CANCERPHARM GROUP

CAPOT CHEMICAL COMPANY LIMITED

CARBOMER INC.

CARBOSYNTH LTD.

CAYMAN CHEMICAL COMPANY

CELANESE, LTD.

CENTEC INTERNATIONAL

CENTRAL GLASS CO. LTD.- FINE
CHEMICAL DEPARTMENT

CERILLIANT CORP.

CHANGZHOU SHUOCHENG
BIOLOGICAL TECHNOLOGY CO. LTD

CHANGZHOU WELTON CHEMICAL
CO.,LTD.

CHARKIT CHEMICAL CORPORATION
CHATTEM CHEMICALS, INC.

CHEM-IMPEX INTERNATIONAL, INC.

CHEMALONG LABS LLC

CHEMBASICS PTY LTD

CHEMBRIDGE CORPORATION
(BUILDING BLOCKS)

Updated Feb. 2021

CHEMBRIDGE CORPORATION
(SCREENING COMPOUNDS)

Updated Mar. 2021

CHEMCONTRACT RESEARCH INC.

CHEMDESIGN CORP. (A BAYER
COMPANY)

CHEMDIV (BUILDING BLOCKS)

Updated Apr. 2021

CHEMDIV (DISCOVERY CHEMICALS)

New Jan. 2021

CHEMFOCUS, LLC

CHEMPACIFIC CORPORATION

CHEMSAMPCo, LLC

CHEMSERVICE, INC.

CHEMADA

CHEMATEK S.P.A.

CHEMBO PHARMA Co LTD

CHEMCIA SCIENTIFIC, LLC

CHEMFUTURE PHARMATECH

CHEMGENX LLC

CHEMI S.P.A.

CHEMICAL TECHNOLOGIES &
INVESTIGATIONS, LTD.

CHEMICREA INC.

CHEMIELIVA PHARMACEUTICAL CO.,
LTD

Updated Sep. 2020

CHEMINISCHE LABORTORIEN
DR. SOENKE PETERSEN

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

ChemACX Vendor List

CHEMISPHERE LIMITED

CHEMIVATE LIMITED

CHEMIZON INC.

CHEMODEX LTD (BUILDING BLOCKS)

CHEMODEX LTD (SCREENING COMPOUNDS)

CHEMSCENE LLC (BUILDING BLOCKS)

CHEMSCENE LLC (REFERENCE COMPOUNDS)

CHEMTEC LEUNA - SPECIALTY CHEMICALS AND FINE CHEMICALS

CHEMTOS, LLC

CHEMTURA CORPORATION

CHEMZON SCIENTIFIC

CHENGDU GT BIOCHEM Co.,LTD.

CHESS GMBH

CHEVRON PHILLIPS CHEMICAL COMPANY LLC

CHIREX LTD.

CHINA NATIONAL CHEMICAL CONSTRUCTION CORPORATION

CHIRACON GMBH

CHIRAGENE CORPORATION (FORMERLY CELGENE)

CHIRALSTAR INC

CHIROBLOCK GMBH

CHIROTECH TECHNOLOGY LIMITED

CHIRON AS

Updated Oct. 2020

CHONGQING JIULONGPO YUHUA CHEMICALS Co., LTD.

CHUGAI KASEI Co., LTD.

CILAG AG

CITY CHEMICAL LLC

CIVENTI CHEM

CLARIANT CORPORATION

CLASSIC FLAVORS & FRAGRANCES

CLEARSYNTH INC.

COALITE CHEMICALS

COMBI-BLOCKS, INC.

COMBIPHOS CATALYSTS, INC.

COOL PHARM, LTD.

CORVINUS CHEMICALS

CREANOVA, INC. (FORMERLY HÜLS AMERICA, INC.)

CUSCHEM INC.

CYMTEK LABORATORIES, LIMITED

CYTH CHEMICAL TRADE Co., LTD.

-D-

D&O CHEMICALS, INC.

D.F. GOLDSMITH CHEMICALS & METALS Co.

DC CHEMICALS

New Aug. 2020

DSM FINE CHEMICALS INC.

DAICEL CHEMICAL INDUSTRIES, LTD.

DAIKIN INDUSTRIES LTD

DAISO Co., LTD.

DALTON PHARMA SERVICES INC.

DANNIER CHEMICAL, INC.

DANYANG HENGAN CHEMICAL Co., LTD.

DAVOS CHEMICAL CORP.

DAYBIOCHEM TECHNOLOGY Co., LTD
DE VOIT ASSOCIATES

DEBYE SCIENTIFIC

Updated Aug. 2020

DEEPAK CHEMICALS GROUP

DEEPWATER CHEMICALS

DEGUSSA CORP.

DHRUVTARA CHEMICALS PVT LTD

DIAGNOSTIC CHEMICALS LIMITED

DIGITAL SPECIALTY CHEMICALS, INC.

DINAMITE DIPHARMA SPA

DIOSYNTH B.V.

DISCOVERY MOLECULES

DISHMAN USA

DIVERCHIM

DIXIE CHEMICAL COMPANY, INC.

DOTTIKON EXCLUSIVE SYNTHESIS AG

DYSTAR LP

DYNAMIT NOBEL GMBH

-E-

ECA INTERNATIONAL CORPORATION

EMD CHEMICALS INC.

EMKA-CHEMIE

EASTAR CHEMICAL CORPORATION

EASTMAN CHEMICAL COMPANY

ECHELON BIOSCIENCES INC.

ELAN CHEMICAL COMPANY, INC.

ELECTRON MICROSCOPY SERVICES

ELEMENT STORE

ELKIMIA INC.

ELPAR INTERNATIONAL INC.

ENAMINE LTD.

Updated Jan. 2021

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

ChemACX Vendor List

ENDOTHERM GMBH	FINETECH INDUSTRY LTD	GENERAL INTERMEDIATES OF CANADA
ENNOVA MEDCHEM GROUP, INC.	FINORNIC CHEMICALS (I) PVT. LTD.	GENERICHEM CORPORATION
E NOVATION CHEMICALS LLC	FIRMENICH SA	GENOSYS (A SIGMA ALDRICH BRAND)
ERNESTO VENTOS S.A.	FIRST CHEMICAL CORP.	GENZYME PHARMACEUTICALS
ERREGIERRE INDUSTRIA CHIMICA S.P.A	FISHER SCIENTIFIC (U.S.)	GEORGE UHE COMPANY, INC.
ESPRIT CHEMICAL COMPANY	FISHER SCIENTIFIC UK	GREATFIELD MINERAL AND CHEMICAL INC.
ESSTECH, INC.	FLEXCHEM BV	GREENCHEM USA
EUROLABS LIMITED	FLUKA CHEMICAL CORP.	GUANGZHOU TOPWORK INTERNATIONAL CO. LTD.
EUROPEAN TECHNOLOGIES INC.	FLUOROCEM GROUP -- SCALE UP SOLUTIONS	GWALIOR CHEMICAL INDUSTRIES LTD.
EVOBLOCKS	FLUOROCEM UK	
EXCENEN PHARMATECH (BIOCHEMICALS) Updated Jan 2021	FLUOROUS TECHNOLOGIES, INC.	-H-
EXCHEM ORGANICS	FOCUS SYNTHESIS LLC	H&S CHEMICAL CO. INC.
EXIMED LABORATORY	FRAMOCEM (CHAYSECEM, NA DISTRIBUTOR)	HAISO PHARMCEM
EXCLUSIVE CHEMISTRY LTD	FRINTON LABORATORIES Updated Jun. 2020	HALOCARBON PRODUCTS CORP.
EXPANSIA S.A.	FRONTIER SCIENTIFIC SERVICES INC. Updated Apr. 2021	HANDE SCIENCES US
-F-	FRONTIER SCIENTIFIC, INC.	HANGZHOU APICHEM TECHNOLOGY CO., LTD. Updated Aug. 2020
F2 CHEMICALS LTD.	FRUTAROM (LABORATORY CHEMICALS)	HANGZHOU PHARMA & CHEM. CO. LTD
FCH GROUP	FRUTAROM LTD (AROMA CHEMICALS)	HANGZHOU TRYLEAD CHEMICAL TECHNOLOGY CO., LTD
FTECH, INC.	FUJI HUNT SPECIALTY PRODUCTS	HANGZHOU YUHAO CHEMICAL TECHNOLOGY CO., LTD
FABBRICA ITALIANA SINTETICI	-G-	HANSHAN AGILEBIOCHEM
FAR RESEARCH	G&J RESEARCH CHEMICALS LTD.	HETCAT
FARCHEMIA DIVISION OF TESSENDERLO CHEMIE	G.J. CHEMICAL COMPANY, INC.	HETEROCYCLICS INC.
FARMAK, A.S.	GFS CHEMICALS, INC.	HEXAGON LABS INC.
FINE CHEMICAL Co., LTD. LANZHOU BELFAST	GNI CHEMICALS CORPORATION	HIGH PERFORMANCE PRODUCT ENGINEERING
FINE CHEMICALS CORPORATION (PTY) LTD.	GANES CHEMICALS INC.	
FINE ORGANICS LTD., SEAL SANDS	GAYLORD CHEMICAL CORPORATION	
	GELEST INC.	

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588- FAX 1 617 588- EMAIL
EU 00 800 875 UK +44 1223 JP 0120 73 80C informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

ChemACX Vendor List

HOKKO CHEMICAL INDUSTRY CO., LTD.
- FINE CHEMICALS DIVISION

HOVIONE GROUP

HUHUGENE INC.

-I-

IAN HETEROCYCLE CO., LTD.

ISP FINE CHEMICALS (FORMERLY
GAF)

IVD (A SIGMA-ALDRICH BRAND)

ICON SERVICES

IHARA CHEMICAL INDUSTRY CO., LTD.

IHARA NIKKEI CHEMICAL INDUSTRY
CO., LTD.

INFARMATIK, INC.

INABATA AMERICA CORP.

INDOFINE CHEMICAL COMPANY,
INC.

INDUSTRIA CHIMICA MILANESE S.P.A.

INNOCEHEM CHEMICAL CO., LTD.

INSIGHT CHEMICAL SOLUTIONS

INTERBIOSCREEN LTD.
(BIOLOGICALS)

INTERBIOSCREEN LTD. (BUILDING
BLOCKS)

INTERBIOSCREEN LTD (SCREENING
COMPOUNDS)

INTERCHEM CORP.

INTERNATIONAL FLAVORS &
FRAGRANCES INC.

IOLiTEC GMBH & Co KG

IRIS BIOTECH GMBH

ISOCHEM, (CHAYSECHEM, NA
DISTRIBUTOR)

ISOTEC INC.

-J-

J&K SCIENTIFIC LTD

J&W PHARMLAB, LLC
Updated Jul. 2020

J.P. PHARMA-CHEM INDUSTRIES

J.H. CALO COMPANY

J.T. BAKER

JPM2 PHARMACEUTICALS, INC.

JRD FLUOROCEMICALS LTD
Updated Sep. 2020

JIANGSU INSTITUTE OF ECOMONES

JIANGSU ZHONGDAN CHEMICAL
IMP. & EXP. CO., LTD.

JOST CHEMICAL

JUNSEI CHEMICAL CO., LTD

JUPITER BIOSCIENCES LTD

JUPITER CHEMICALS, INC.

-K-

K-I CHEMICAL INDUSTRY CO.,
LTD.

KADEN BIOCHEMICALS GMBH

KANEKA CORPORATION - FINE
CHEMICALS DIVISION

KANTO CHEMICAL CO., INC.

KATWIJK CHEMIE BV

KAIRONKEM

KEMIRA FINE CHEMICALS

KEY MOLECULAR CORPORATION

KEY ORGANICS BIONET
(BIOCHEMICALS)
Updated May. 2020

KEY ORGANICS BIONET (BUILDING
BLOCKS)

KEY ORGANICS BIONET (COMBINED-
SCREENING-FRAGMENTS)

KEY ORGANICS BIONET (FRAGMENT
LIBRARY)
Updated Oct. 2020

KEYINTERMEDIATE

KINENTIA BIOSCIENCES

KINGCHEM INC.

KINGSCIENTIFIC

KINGSH CHEMICALS INC.

KISHIDA BUILDING BLOCKS

KISHIDA CHEMICAL CO., LTD

KLAUS F. MEYER GMBH

KOEI CHEMICAL COMPANY, LTD

KYOWA HAKKO KOGYO Co., LTD.

-L-

LC LABORATORIES

LGM PHARMA

LKT LABORATORIES, INC.

LABNETWORK

LABORATORI MAG SPA

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

ChemACX Vendor List

LAN PHARMATECH LTD.

Updated Apr. 2021

LANCASTER SYNTHESIS INC.

LANSDOWNE CHEMICALS PLC

LANXESS AG

**LANZHOU TECH BIOCHEMICAL CO.
LTD.**

**LAPORTE PERFORMANCE CHEMICALS
(FORMERLY INSPEC GROUP)**

LATOXAN

LEADGEN LABS, LLC

**LIFE CHEMICALS INC. (BUILDING
BLOCKS)**

**LIFE CHEMICALS INC. (SCREENING
COMPOUNDS)**

Updated May 2020

LIVCHEM LOGISTICS GMBH

LOBA FEINCHEMIE AG

LONZA LTD.

LUKNOVA INC

LUMAC*LSC B.V.

LYONDELL CHEMICAL COMPANY

-M-

MM (A SIGMA-ALDRICH BRAND)

MP BIOMEDICALS

MV LABORATORIES, INC.

MACFARLAN SMITH

**MALLINCKRODT LABORATORY
CHEMICALS**

MANCHESTER ORGANICS LTD

Updated Mar. 2021

MARKER GENE TECHNOLOGIES, INC

**MARSHALLTON RESEARCH
LABORATORIES, INC.**

MATERIA INC.

MATRIX SCIENTIFIC

**MAYBRIDGE CHEMICAL COMPANY,
LTD.**

MCULE

Updated Nov. 2020

MCULE ULTIMATE

Updated Apr. 2021

**MEDCHEMEXPRESS LLC (SCREENING
LIBRARIES)**

MEDKOO BIOSCIENCES

MEDICAL ISOTOPES, INC.

MEGAZYME INTERNATIONAL

MELFORD LABORATORIES LTD.

MENAI ORGANICS

MERCK SCHUCHARDT OHG

**METALL-CHEMIE GOERRIG GMBH &
CO. KG**

MICROCOMBICHEM E.K.

MILESTONE PHARMTECH LLC

**MILLIONLINK (TIANJIN)
INTERNATIONAL TRADE CO., LTD.**

MINOL ACIDS & CHEMICALS PVT. LTD.

MITCHELL COTTS CHEMICALS

MITENI SPA

MITSUBISHI RAYON CO., LTD.

**MITSUI CHEMICALS, INC.
MOLBRIDGE**

MOLE CHEMICAL CORPORATION

MOLECULAR BIOSCIENCES

MOLECULAR PROBES, INC.

MOLEKULA

MOLEPEDIA Co., LTD.

MORAVEK BIOCHEMICALS

MORFLEX, INC.

MORRE-TEC INDUSTRIES, INC.

MORTON INTERNATIONAL, INC.

MUSECHEM

Updated Jan. 2021

-N-

NSC TECHNOLOGIES

NACALAI TESQUE INC

NAMIKI-SHOJI Co., LTD.

Updated Sep. 2020

**NANJING DAXIAN CHEMICAL
INSTITUTE LTD.**

**NANJING PHARMACEUTICAL FACTORY
Co., LTD.**

NANJING RAYMON BIOTECH Co., LTD

NANJING REDLINE LTD.

NEOMPS, INC.

NEOSYN LABORATORIES INC.

NETQEM

**NEW ZEALAND PHARMACEUTICALS
LIMITED**

NICHIMEN AMERICA INC.

NIGU CHEMIE GMBH

NIHON NOHYAKU Co., LTD.

NIPA HARDWICKE INC.

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

ChemACX Vendor List

NIPPOH CHEMICALS CO., LTD.

NIPPON SHOKUBAI CO. LTD.

NORCHIM S.A.

NORQUAY TECHNOLOGY INC. USA

NORTHEAST PHARMACEUTICAL
GROUP CO. LTD

NOVABIOCHEM (EMD BIOSCIENCES,
INC.)

NOVASYN ORGANICS PVT. LTD.

NOVELKORE

NOVOCHEMY LTD.

-O-

OAKWOOD PRODUCTS, INC.

OCCIDENTAL PETROLEUM
CORPORATION

OMEGACHEM, INC.

ONTARIO PHARMA RESEARCH
CHEMICALS

ONYX SCIENTIFIC LIMITED

OPTIMA CHEMICAL GROUP LLC

ORGANICA FEINCHEMIE GMBH
WOLFEN

ORGANIX, INC.

ORGANOCEM LTD

ORGAPHARM-AXYNTIS GROUP

ORGASYNTH

ORIBASE PHARMA

ORION CORPORATION FERMION

OSAKA ORGANIC CHEMICAL INDUSTRY,
LTD.

OTAVA

OXCHEM CORPORATION

-P-

PFL FINE CHEMICALS GMBH

PK CHEM

PNK UNITED, INC.

PPG-SIPSY

PALINDENT ORGANIC
INTERMEDIATES

PANCHIM

PARADIGM SCIENTIFIC

PARISH CHEMICAL COMPANY

PARKWAY SCIENTIFIC LLC

PEAKDALE FINE CHEMICALS LIMITED

PENG TENG SHANGHAI FINE CHEMICAL
CO.

PENN SPECIALTY CHEMICALS, INC.

PENTAGON CHEMICALS

PEPTeCH CORPORATION

PERKINELMER CHEMICALS

PFALTZ & BAUER, INC.

PFANSTIEHL LABORATORIES, INC.

PHARMABLOCK R&D Co. LTD.

PHARMACORE, INC.

PHARMABRIDGE

PHARMACEUTICAL WORKS POLPHARMA
S.A.

PHARMACIA & UPJOHN COMPANY

PHARMCO PRODUCTS INC.

PHARMTEN CHEMICAL Co.,
LIMITED

PHOENIX CHEMICALS

PHYTOLAB GMBH & Co. KG

PLATTE VALLEY SCIENTIFIC, LLC

POLI INDUSTRIA CHIMICA S.P.A.

POLIUM TECHNOLOGIES, INC.

POLYCARBON INDS., INC.

POLYMER LABORATORIES LTD

POLYORGANIX, INC.

POLYSCIENCES, INC.

PORSE FINE CHEMICAL Co., LTD.

PRATAP ORGANICS

PRECISION BIOCHEMICALS, INC.

PRECURSOR CHEMICALS, INC.

PRINCETON BIOMOLECULAR
RESEARCH, INC. (BUILDING BLOCKS)

PRINCETON BIOMOLECULAR
RESEARCH, INC. (SCREENING
COMPOUNDS)

Updated Oct. 2020

PROCHEM, INC.

PRODOTTI CHIMICI E ALIMENTARI
SPA

PURE CHEMISTRY SCIENTIFIC INC.

-Q-

QUANTA BIODESIGN, LTD.

QUEEN'S UNIVERSITY OF BELFAST

-R-

R&S CHEMICALS, INC.

R.I. CHEMICAL, INC.

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

ChemACX Vendor List

RCA KG (REUTER CHEMISCHE
APPARATEBAU KG)

RSP AMINO ACID ANALOGUES, INC.

RARE CHEMICALS GMBH
Updated Oct. 2020

RARE CHEMICALS GMBH (SCREENING
LIBRARY)

RASCHIG AG

REAXON GBR

REILLY INDUSTRIES, INC.

RELAX LTD.

RELIABLE BIOPHARMACEUTICAL
CORPORATION

RESEARCH ORGANICS

RHODIA FINE ORGANICS

RICHMAN CHEMICAL INC.

RIEDEL-DE HAËN

RIEKE METALS, INC.
Updated Apr. 2021

RIEKE METALS, INC. (NOVEL
CHEMICAL SOLUTIONS)

ROBINSON BROTHERS LIMITED

ROCHE

ROCKWOOD LITHIUM GMBH

ROHM & HAAS PROCESS CHEMICALS

ROHNER LTD.

ROMIL LTD

RUETGERS-NEASE CORPORATION

RYAN SCIENTIFIC, INC.

RYSCOR SCIENCE, INC.

-S-

S.I.M.S. S.R.L.

S.K. CHEMICAL INDUSTRIES

SEAC

SK ENERGY AND CHEMICAL, INC.

SKW CHEMICALS

SNPE CHEMICALS, GROUPE SNPE

SALFORD ULTRAFINE CHEMICALS &
RESEARCH LTD.

SAMCHULLY PHARMACY COMPANY,
LTD.

SANFUR USA, INC.

SANOCHEMIA PHARMAZEUTIKA AG

SANTAI LABS

SANTIAGO

SAUDI BASIC INDUSTRIES
CORPORATION

SCHERING AG

SELENA CHEMICALS

SELLECKCHEM

SENN CHEMICALS

SENSACHEM, INC.

SHAANXI BONO BIOTECH Co., LTD

SEQUOIA RESEARCH PRODUCTS LTD

SHANDONG YOUBANG BIOCHEMICAL
TECHNOLOGY Co., LTD.

SHANGHAI ACT CHEMICAL LIMITED

SHANGHAI AOBO BIO-PHARMACEUTICAL
TECHNOLOGY Co., LTD.

SHANGHAI CHEMHERE Co., LTD

SHANGHAI ELITTES ORGANICS Co., LTD.

SHANGHAI EXCELLENT CHEMICAL Co., LTD.

SHANGHAI FWD CHEMICALS Co., LTD.

SHANGHAI HAORYUAN CHEMEXPRESS
Co., LTD

SHANGHAI INTERNATIONAL
PHARMACEUTICAL Co., LTD

SHANGHAI PI CHEMICALS LTD.

SHANGHAI PHARMA TECH Co., LTD.
(PARLING)

SHANGHAI PU-JIE FRAGRANCE Co.,
LTD.

SHANGHAI RACE CHEMICAL Co., LTD

SHANGHAI REITO CHEMICAL
TECHNOLOGY Co., LTD.

SHANGHAI RONGKANG FINE
CHEMICAL Co., LTD

SHANGHAI SINOFLUORO SCIENTIFIC
CORPORATION LTD.

SHANGHAI SPECBIOCHEM CHEMISTRY
GROUP LIMITED

SHELL CHEMICAL LP

SICHUAN TONGSHENG AMINO ACID
Co., LTD

SIGALD (A SIGMA-ALDRICH BRAND)

SIGMA ALDRICH DISCOVERY
CHEMICALS

SIGMA CHEMICAL COMPANY

SIGMA-ALDRICH BRAND CHEMICALS

SIGMA-ALDRICH JAPAN K.K.

SILAR LABORATORIES - A DIVISION OF
WRIGHT CORPORATION

SIMAFEX

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

ChemACX Vendor List

SINOCHEM HEBEI TIANJIN IMPORT -
EXPORT CORPORATION

SINON CORPORATION

SMALL MOLECULES

SOHVE CORPORATION

SOLTEC VENTURES, INC.

SPECS (BUILDING BLOCKS)

SPECS (NATURAL PRODUCTS)

SPECS (SCREENING COMPOUNDS)
New May 2020

SPECTRUM CHEMICAL & LABORATORY
PRODUCTS

SPECTRUM INFO LTD.

SPHEROTECH, INC.

SPIROCHEM AG

STANFORD MATERIALS COMPANY

STREM CHEMICALS, INC.

SUMIKIN CHEMICAL Co. LTD.

SUMITOMO SEIKA CHEMICALS Co.,
LTD.

SUN BIOCHEMICALS, INC.

SUNOCO CHEMICALS

SUNSHINE CHEMICALS COMPANY

SUNTTO CHEMICAL

SUPELCO (A SIGMA ALDRICH BRAND)

SUPRIYA CHEMICALS

SUZHOU CHONTECH PHARMACHEM
TECHNOLOGY Co., LTD.

SUZHOU ROVATHIN PHARMA TECH Co.,
LTD.

SUZHOU RYAN PHARMACHEM
TECHNOLOGY Co., LTD.

SUZHOU SINOERA CHEM Co., LTD

SUZHOU YIELD PHARMACEUTICAL
R&D Co., LTD

SYNAFFIX BV

SYNCHEM, INC.

SYNFINE RESEARCH

SYNHET INC.
Updated Apr. 2021

SYNINNOVA

SYNQUEST LABORATORIES, INC.

SYNCHEM LABORGEMEINSCHAFT UG &
Co. KG

SYNCOZYMES (SHANGHAI) Co., LTD.

SYNERGY-SCIENTIFIC LTD.

SYNGANIX SCIENCE INC.

SYNOVEL LABORATORY Co., LTD.

SYNOVO GMBH

SYNTAI CHEMICALS &
PHARMACEUTICALS, LTD.

SYNTASTIC AB

SYNTECH SOLUTION

SYNTESIA CHEMIE GMBH

SYNTHE163

SYNTHETECH

SYNTHON CHEMICALS GMBH & Co. KG

SYNTHON CHIRAGENICS
CORPORATION

SYNTHONIX CORPORATION

SYNTIDE MEDI-CHEM TECHNOLOGY
Co., LTD.

SÄUREFABRIK SCHWEIZERHALL
(SF-CHEM)

-T-

TC CHEMICALS Co., LTD.

TOSLAB

TAIZHOU KANGDUO SEA BIOCHEMICAL &
PHARMACEUTICAL Co., LTD.

TAKASAGO INTERNATIONAL CORP.

TAOKA CHEMICAL Co., LTD.

TECOLAND CORPORATION

TENOVA PHARMACEUTICALS INC

TEVA GROUP API DIVISION

TEXAS BIOCHEMICALS INC.

THE NEW CENTURY CHEMICAL &
PHARMACEUTICAL ENTERPRISE

THE SHEPHERD CHEMICAL COMPANY

TOCRIS BIOSCIENCE

TOKUYAMA CORPORATION

TOKYO CHEMICAL INDUSTRY Co., LTD.

TOMAH PRODUCTS, INC.

TORAY INDUSTRIES, INC.

TORONTO RESEARCH CHEMICALS,
INC.

TOSOH CORPORATION

TRACTUS COMPANY LIMITED

TRANS WORLD CHEMICALS INC.

TRATAMIENTOS GUADALQUIVIR S.A.
(TRAGUSA)

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588- FAX 1 617 588- EMAIL
EU 00 800 875 UK +44 1223 JP 0120 73 800 informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

ChemACX Vendor List

TRIFARMA SPA

TRINAPCO, INC.

TRIOMPHE FINE CHEMISTRY CO.

TRIPLEBOND CORP.

TYGER SCIENTIFIC INC.

-U-

USDAC COMPANY

UBE INDUSTRIES, LTD. - FINE
CHEMICALS DIVISION

UBICHEM PLC

UNAVERA CHEMLAB GMBH

UNITED STATES PHARMA COPEIA (A
SIGMA ALDRICH BRAND)

-V-

VWR INTERNATIONAL GMBH

VWR INTERNATIONAL, LLC

VANDEMARK GROUP

VARSAL INSTRUMENTS, INC.

VESINO INDUSTRIAL CO.

VETEC QUIMICA FINA LTDA

VIJAYA PHARMACEUTICALS

VIS FARMACEUTICI SPA

VITAS-M LABORATORY, LTD.
Updated Aug. 2020

VITAS-M LABORATORY, LTD.
(SCREENING COMPOUNDS)
Updated Apr. 2021

VIVA CORPORATION

VIWIT PHARMATECH CO., LTD.

-W-

WACKER CHEMICALS (USA)

WAKO PURE CHEMICAL INDUSTRIES,
LTD.

WATANABE CHEMICAL INDUSTRIES,
LTD.

WATEC LABORATORIES, INC.

WATERSTONE TECHNOLOGY, LLC

WHYTE CHEMICALS LIMITED

WILLIAM BLYTHE LTD.

WILSHIRE CHEMICAL CO., INC.

WINNER GROUP INTERNATIONAL LTD.

WONDA SCIENCE INC.

WONDERCHEM CO., LIMITED

WUXI APPTec Co., LTD

WUHAN ZHONGFU BIO-TECH CO.,
LTD.

WUJIN LINCHUAN CHEMICAL
FACTORY

WUXI D-STONE INTERNATIONAL CO.,
LTD.

WYCHEM LIMITED

-X-

XPR INDUSTRIAL (CHINA) LIMITED

XIMO HUNGARY, LTD.

XINCHEM COMPANY

-Y-

YAMAKAWA CHEMICAL

YUANHONG SINO CHEMICAL

YUHAN CORPORATION

-Z-

ZAMBON GROUP

ZENECA SPECIALTIES

ZEON CORP

ZERENEX BIOACTIVES

ZERENEX MOLECULAR LIMITED
Updated Mar. 2021

ZHEJIANG KANGYU PHARMACEUTICAL
CO.,LTD

ZHEJIANG OUHUA CHEMICAL
IMP.&EXP.CO.,LTD

ZHENGTAI PHARMA & CHEM CO., LTD.

ZYLEXA PHARMA LTD

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 800

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights