

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.

3RCHEM

8S LABORATORIES LLC

-A-

**A&A LIFE SCIENCE INC. DBA
AVA CHEM SCIENTIFICS**

A&C PHARMTECH, INC

A&J PHARMATECH

A.G. SCIENTIFIC

A1 BIOCHEM LABS
Updated Sep. 2018

AA BLOCKS LLC
New Sep. 2018

AB CHEM TECHNOLOGIES LLC

ABCR GMBH & Co. KG

ABSCO LTD

**ABX ADVANCED BIOCHEMICAL
COMPOUNDS**

AE CHEM

AK SCIENTIFIC, INC.
Updated Oct. 2018

AKL RESEARCH LLP

AOBCHEM USA

AOKBIO Co., LTD
Updated Apr. 2018

AOKCHEM Co., LTD.

APAC PHARMACEUTICAL

AQCHEM LLC

ASW MEDCHEM, INC.

AURUM PHARMATECH LLC

AARON CHEMISTRY

ABAMACHEM LTD

ABBEY COLOR

**ABBOTT LABORATORIES CHEMICAL
DIVISION**

ABBY PHARMATECH LLC

ABERJONA LABORATORIES

ABOVCHEM LLC

ABSOLUTE CHIRAL
Updated July 2018

ACADECHEM COMPANY LIMITED

ACADEMIQUE PHYTOCANADA

ACCELA CHEMBIO Co., LTD.
Updated June 2018

ACCELEDEV CHEMICAL LLC

ACCUSTANDARD, INC.

ACE SYNTHESIS LLC

ACENTEX SCIENTIFIC, 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.
Updated Aug. 2018

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
New June 2018

ALFA PYRIDINES

All specifications subject to change

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

EU 00 800 875 UK +44 1223

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

FAX 1 617 588-

JP 0120 73 800

EMAIL

informatics.info@perkinelmer.com

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2018 PerkinElmer Inc. All Rights

ChemACX Vendor List

ALLIEDSIGNAL INC. SPECIALTY CHEMICALS

AMBEED, INC.

New Dec. 2018

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.

ANGELINI FINE CHEMICALS

ANGENE INTERNATIONAL LIMITED

Updated Aug. 2018

ANGUS CHEMICAL COMPANY

ANNKER ORGANICS

ANTIBIOTICOS SPA

ANWARD CHEMICALS

APEIRON SYNTHESIS

APEX SCIENTIFIC INC.

APEXMOL TECHNOLOGY

APITOXIN

APOLLO SCIENTIFIC LTD. (BIOCHEMICALS)

Updated Sep. 2018

APOLLO SCIENTIFIC LTD. (FLUROCHEMICALS)

Updated Oct. 2018

APOLLO SCIENTIFIC LTD. (INTERMEDIATES)

Updated Nov. 2018

APOLLO SCIENTIFIC LTD. (SUPPORT RESEARCH LIBRARY)

Updated Dec. 2018

APPLICHEM

ARBONOVA OY AB

ARCTOM CHEMICALS

New Apr. 2018

ARKÉ ORGANICS SRL

AROMA CIRCLE

ARONIS

New July 2018

ARRAN CHEMICAL CO.

ARYLCHEM CO.

ARZA BIOSCIENCE LIMITED

Updated Sep. 2018

ASAHI GLASS COMPANY - FLUROCHEMICALS 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 Aug. 2018

AVIABOR FINE BORON COMPOUNDS

AVONCHEM/CHROMOS EXPRESS LTD

AXITEC CO., LTD

AZEPINE LTD

-B-

BACHEM AG

BASF CHEMICAL INTERMEDIATES

BE-ASSOCIATES, LLC

BLD PHARMATECH LTD

Updated Jan. 2019

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.

BIDDLE SAWYER CORPORATION

BIDE PHARMATECH

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 1998-2018 PerkinElmer Inc. All Rights

ChemACX Vendor List

BIO FARMA LTD.

BIO-RESEARCH PRODUCTS, INC.

BIOBLOCKS, INC.

BIOQUADRANT INC.

BIOVISION, INC.

BIOCHEMIE GMBH

BIOCHEMPEG

BIOCORE PHARMTECH LIMITED

BIOFINE INTERNATIONAL INC.

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

BURDICK & JACKSON, INC.

BUSH BOAKE ALLEN INC.

BUTT PARK LTD

-C-

C/D/N ISOTOPES INC.

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.

Updated Nov. 2018

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 Dec. 2018

CHEMBRIDGE CORPORATION
SCREENING COMPOUNDS

Updated Nov. 2018

CHEMCONTRACT RESEARCH INC.

CHEMDESIGN CORP. (A BAYER
COMPANY)

CHEMFOCUS, LLC

CHEMPACIFIC CORPORATION

CHEMSAMP Co, LLC

CHEMSERVICE, INC.

CHEMADA

CHEMATEK S.P.A.

CHEMBO PHARMA Co LTD

Updated May 2018

CHEMCIA SCIENTIFIC, LLC

CHEMFUTURE PHARMATECH

CHEMGENX LLC

CHEMI S.P.A.

CHEMICAL TECHNOLOGIES &
INVESTIGATIONS, LTD.

CHEMICREA INC.

CHEMIELIVA PHARMACEUTICAL Co.,
LTD

Updated June 2018

CHEMINISCHE LABORTORIEN
DR. SOENKE PETERSEN

CHEMISPHERE LIMITED

CHEMIVATE LIMITED

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 1998-2018 PerkinElmer Inc. All Rights

ChemACX Vendor List

CHEMIZON INC.

CHEMSCENE LLC (BUILDING BLOCKS)

New Nov. 2018

CHEMSCENE LLC (REFERENCE COMPOUNDS)

New Nov. 2018

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

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.

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

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.

ENDOTHERM GMBH

ENNOVA MEDCHEM GROUP, INC.

ERNESTO VENTOS S.A.

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 1998-2018 PerkinElmer Inc. All Rights

ChemACX Vendor List

ERREGIERRE INDUSTRIA CHIMICA
S.P.A

ESPRIT CHEMICAL COMPANY

ESSTECH, INC.

EUROLABS LIMITED

EUROPEAN TECHNOLOGIES INC.

EVOBLOCKS

EXCHEM ORGANICS

EXCLUSIVE CHEMISTRY LTD

EXPANSIA S.A.

-F-

F2 CHEMICALS LTD.

FCH GROUP
Updated Aug. 2018

FTECH, INC.

FABBRICA ITALIANA SINTETICI

FAR RESEARCH

FARCHEMIA DIVISION OF
TESSENDERLO CHEMIE

FARMAK, A.S.

FINE CHEMICAL CO., LTD. LANZHOU
BELFAST

FINE CHEMICALS CORPORATION
(PTY) LTD.

FINE ORGANICS LTD., SEAL SANDS

FINE TECH INDUSTRY LTD

FINORNIC CHEMICALS (I) PVT. LTD.

FIRMENICH SA

FIRST CHEMICAL CORP.

FISHER SCIENTIFIC (U.S.)

FISHER SCIENTIFIC UK

FLEXCHEM BV

FLUKA CHEMICAL CORP.

FLUOROCHEM GROUP -- SCALE UP
SOLUTIONS

FLUOROCHEM UK
Updated June 2018

FLUOROUS TECHNOLOGIES, INC.

FOCUS SYNTHESIS LLC

FRAMOCHEM (CHAYSECHEM, NA
DISTRIBUTOR)

FRINTON LABORATORIES
Updated Sep. 2018

FRONTIER SCIENTIFIC SERVICES INC.
Updated Dec. 2018

FRONTIER SCIENTIFIC, INC.
Updated May 2018

FRUTAROM (LABORATORY CHEMICALS)

FRUTAROM LTD (AROMA CHEMICALS)

FUJI HUNT SPECIALTY PRODUCTS

-G-

G&J RESEARCH CHEMICALS LTD.

G.J. CHEMICAL COMPANY, INC.

GFS CHEMICALS, INC.

GNI CHEMICALS CORPORATION

GANES CHEMICALS INC.

GAYLORD CHEMICAL CORPORATION

GELEST INC.

GENERAL INTERMEDIATES OF
CANADA

GENERICHEM CORPORATION

GENOSYS (A SIGMA ALDRICH BRAND)

GENZYME PHARMACEUTICALS

GEORGE UHE COMPANY, INC.

GREATFIELD MINERAL AND
CHEMICAL INC.

GREENCHEM USA

GUANGZHOU TOPWORK
INTERNATIONAL CO. LTD.

GWALIOR CHEMICAL INDUSTRIES
LTD.

-H-

H&S CHEMICAL CO. INC.

HAISO PHARMCHEM

HALOCARBON PRODUCTS CORP.

HANDE SCIENCES US

HANGZHOU APICHEM TECHNOLOGY
CO., LTD.
Updated July 2018

HANGZHOU PHARMA & CHEM. CO. LTD

HANGZHOU TRYLEAD CHEMICAL
TECHNOLOGY CO., LTD

HANGZHOU YUHAO CHEMICAL
TECHNOLOGY CO., LTD

HANSHAN AGILEBIOCHEM

HETCAT

HETEROCYCLICS INC.

HEXAGON LABS INC.

HIGH PERFORMANCE PRODUCT
ENGINEERING

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 1998-2018 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)

Updated Aug. 2018

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 Oct. 2018

J.P. PHARMA-CHEM INDUSTRIES

J.H. CALO COMPANY

J.T. BAKER

JPM2 PHARMACEUTICALS, INC.

JRD FLUOROCEMICALS LTD

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

KEY ORGANICS BIONET BUILDING BLOCKS

KEY ORGANICS BIONET COMBINED-SCREENING-FRAGMENTS

KEY ORGANICS BIONET FRAGMENT LIBRARY

KEYINTERMEDIATE

KINENTIA BIOSCIENCES
New Sep. 2018

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.

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 1998-2018 PerkinElmer Inc. All Rights

ChemACX Vendor List

-L-

LC LABORATORIES

LGM PHARMA

LKT LABORATORIES, INC.

LABNETWORK

LABORATORI MAG SPA

LANCASTER SYNTHESIS INC.

LANSDOWNE CHEMICALS PLC

LANXESS AG

LANZHOU TECH BIOCHEMICAL CO.
LTD.

LAPORTE PERFORMANCE CHEMICALS
(FORMERLY INSPEC GROUP)

LATOXAN

LEADGEN LABS, LLC
Updated Sep. 2018

LIFE CHEMICALS INC.

LIVCHEM LOGISTICS GMBH

LOBA FEINCHEMIE AG

LONZA LTD.

LUKNOVA INC

LUMAC*LSC B.V.

LYONDELL CHEMICAL COMPANY
(FORMERLY ARCO CHEMICAL
COMPANY)

-M-

MP BIOMEDICALS (FORMERLY
ICN BIOMEDICALS, INC.)

MV LABORATORIES, INC.

MACFARLAN SMITH

MALLINCKRODT LABORATORY
CHEMICALS

All specifications subject to change

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

EU 00 800 875 UK +44 1223

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MANCHESTER ORGANICS LTD
Updated Oct. 2018

MARKER GENE TECHNOLOGIES, INC

MARSHALLTON RESEARCH
LABORATORIES, INC.

MATERIA INC.

MATRIX SCIENTIFIC
Updated Jan. 2019

MAYBRIDGE CHEMICAL COMPANY,
LTD.

MCULE
Updated Aug. 2018

MEDCHEMEXPRESS LLC (SCREENING
LIBRARIES)
New Nov. 2018

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.

-N-

NSC TECHNOLOGIES

NACALAI TESQUE INC

NAMIKI-SHOJI
Updated Nov. 2018

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.

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.

NOVEL CHEMICAL SOLUTIONS

NOVELKORE

NOVOCHEMY LTD.

-O-

OAKWOOD PRODUCTS, INC.

OCCIDENTAL PETROLEUM
CORPORATION

OMEGACHEM, INC.
Updated Apr. 2018

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
Updated July 2018

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
Updated Dec. 2018

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)
Updated Oct. 2018

PRINCETON BIOMOLECULAR
RESEARCH, INC. (SCREENING
COMPOUNDS)
Updated Dec. 2018

PROCHEM, INC.

PRODOTTI CHIMICI E ALIMENTARI
SPA

PURE CHEMISTRY SCIENTIFIC INC.

-Q-

QUANTA BIODESIGN, LTD.

QUEEN'S UNIVERSITY OF BELFAST

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 1998-2018 PerkinElmer Inc. All Rights

ChemACX Vendor List

-R-

R&S CHEMICALS, INC.

R.I. CHEMICAL, INC.

RCA KG (REUTER CHEMISCHE
APPARATEBAU KG)

RSP AMINO ACID ANALOGUES, INC.

RARE CHEMICALS GMBH

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.

ROBINSON BROTHERS LIMITED

ROCHE
Updated Nov. 2018

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.

SEQUOIA RESEARCH PRODUCTS LTD

SHANDONG YOUNG BIOCHEMICAL
TECHNOLOGY CO., LTD.

SHANGHAI ACT CHEMICAL LIMITED
New Apr. 2018

SHANGHAI AOBO BIO-PHARMACEUTICAL
TECHNOLOGY CO., LTD.

SHANGHAI CHEMHERE CO., LTD

SHANGHAI ELITES ORGANICS CO., LTD.

SHANGHAI EXCELLENT CHEMICAL CO., LTD.

SHANGHAI FWD CHEMICALS CO., LTD.

SHANGHAI HAUYUAN 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
Updated Oct. 2018

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

SIGMA ALDRICH LIBRARY OF RARE
CHEMICALS

SIGMA CHEMICAL COMPANY

SIGMA-ALDRICH BRAND CHEMICALS
Updated Dec. 2018

SIGMA-ALDRICH JAPAN K.K.

SILAR LABORATORIES - A DIVISION OF
WRIGHT CORPORATION

SIMAFEX

SINOCHEM HEBEI TIANJIN IMPORT -
EXPORT CORPORATION

SINON CORPORATION

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 1998-2018 PerkinElmer Inc. All Rights

ChemACX Vendor List

SMALL MOLECULES

SOHVE CORPORATION

SOLTEC VENTURES, INC.

SPECS BUILDING BLOCKS

SPECS NATURAL PRODUCTS

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
Updated Mar. 2018

SUZHOU YIELD PHARMACEUTICAL R&D Co., LTD

SYNAFFIX BV

SYNCHEM, INC.

SYNFINE RESEARCH

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

SYNTEZIA 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
Updated Aug. 2018

TOKUYAMA CORPORATION

TOKYO CHEMICAL INDUSTRY Co., LTD.

TOMAH PRODUCTS, INC.

TORAY INDUSTRIES, INC.

TORONTO RESEARCH CHEMICALS, INC.
Updated July 2018

TOSOH CORPORATION

TRACTUS COMPANY LIMITED

TRANS WORLD CHEMICALS INC.

TRATAMIENTOS GUADALQUIVIR S.A. (TRAGUSA)

TRIFARMA SPA

TRINAPCO, 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 73 800 informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2018 PerkinElmer Inc. All Rights

ChemACX Vendor List

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 PHARMACOPEIA (A
SIGMA ALDRICH BRAND)

-V-

VWR INTERNATIONAL GMBH

VWR INTERNATIONAL, LLC

VANDEMARK GROUP

VARSAL INSTRUMENTS, INC.

VESINO INDUSTRIAL CO.

VETEC QUIMICA FINA LTDA
Updated Oct. 2018

VIJAYA PHARMACEUTICALS

VIS FARMACEUTICI SPA

VITAS-M LABORATORY, LTD.

VIVA CORPORATION

VIWIT PHARMATECH Co., LTD.

-W-

WACKER CHEMICALS (USA)

WAKO PURE CHEMICAL INDUSTRIES,
LTD.

WATANABE CHEMICAL INDUSTRIES,
LTD.

WATEC LABORATORIES, INC.
Updated Mar. 2018

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

XINCHEM COMPANY

-Y-

YAMAKAWA CHEMICAL

YUANHONG SINO CHEMICAL

YUHAN CORPORATION

-Z-

ZAMBON GROUP

ZENECA SPECIALTIES

ZEON CORP

ZERENEX BIOACTIVES

ZERENEX MOLECULAR LIMITED

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 1998-2018 PerkinElmer Inc. All Rights