



Berthold Technologies USA Representative Directory United States

<i>Coverage Area</i>	<i>Company Name</i>	<i>Phone/Website</i>
Steel Measurement Systems	Radiometrics, Inc.	440 245 9977 www.radiometrics.com Kathy Huerner kh@radiometrics.com
Alabama	Hile Controls Inc.	205-620-4000 www.hilecontrolsinc.com Grady Andrews grady.andrews@hilecontrolsinc.com
Alaska	Engineered Process Controls	425 844 8901 www.engineeredprocess.com Greg LeBrun greg@engineeredprocess.com
Arizona	Techstar	832 599 3794 www.techstar.com Phil Deaton pdeaton@techstar.com
Arkansas	TechStar	832 599 3794 www.techstar.com Phil Deaton pdeaton@techstar.com
California	Innovative Process Controls	310 414 3700 www.iprocessc.com David DiDomizio david@iprocessc.com
Colorado	Techstar	832 599 3794 www.techstar.com Phil Deaton pdeaton@techstar.com
Connecticut	Pond Technical Sales, Inc.	860 316 2570 www.pondtechnicalsales.com Bob Osnoe rosnoe@pondtechnicalsales.com
Delaware	North East Technical Sales	215 513 1000 www.netechsales.com Doug Reichert dreichert@netechsales.com
Florida Zip Code 324 & 325	Hile Controls Inc.	205-620-4000 www.hilecontrolsinc.com Grady Andrews grady.andrews@hilecontrolsinc.com
Zip Codes Except 324 & 325	Abbey Mec-Tric	770-840-9191 https://www.abbeymec-tric.com/ Will Thurman sales@abbeymec-tric.com

Georgia		Abbey Mec-Tric	770-840-9191 https://www.abbeymec-tric.com/ Will Thurman sales@abbeymec-tric.com
Hawaii		Central Pacific Controls	808 450 2483 www.cpacificc.com David Didmomizio david@cpacificc.com
Idaho	Zip Codes 832-834; 836-837	Trailblazer Controls Corporation	801 955 3900 www.trailblazercontrols.com Jason Snedaker jason@trailblazercontrols.com
	Zip Codes: 835; 838	Engineered Process Controls	425 686 4450 www.engineeredprocess.com Adam DiDomizio adam@engineeredprocess.com
Illinois		Allied Instrumentation	219 924 4545 Alliedinst.com Lorenzo Nirchi lorenzo.nirchi@alliedinst.com
Indiana	all zip codes except below	Henry M Wood	513 755 4060 www.henrymwood.com Eric Yaeger eric@henrymwood.com
	Zip Codes: 463, 465, 467, & 468	Allied Instrumentation	219 924 4545 Alliedinst.com Lorenzo Nirchi lorenzo.nirchi@alliedinst.com
Iowa		Allied Instrumentation	844 398 9724 alliedinst.com Dawn Whittenburg dawnw@alliedinst.com
Kansas		Techstar	832 599 3794 www.techstar.com Phil Deaton pdeaton@techstar.com
Kentucky	except zip codes 411, 412, 415, 416	Henry M Wood	513 755 4060 www.henrymwood.com Eric Yaeger eric@henrymwood.com
	zip code 411, 412, 415, 416	Regional Sales Manager:	Becky Olliges becky_olliges@berthold-us.com 865 323 0876
Louisiana		Techstar	832 599 3794 www.techstar.com Phil Deaton pdeaton@techstar.com
Maine		Pond Technical Sales, Inc. d/b/a NECI	507 851 3560 neci.com Alic Albright aalbright@neci.com

Maryland	zip codes 215 & 217	Hareza Technical Sales I	(724) 942-4966 www.hareza.com Walt Lemunyon office@hareza.com
	all other areas	North East Technical Sales	215 513 1000 www.netechsales.com Doug Reichert dreichert@netechsales.com
Massachusetts		Pond Technical Sales, Inc.	860 316 2570 www.pondtechnicalsales.com Bob Osnoe rosnoe@pondtechnicalsales.com
Michigan		Allied Instrumentation	(248) 478-8892 Alliedinst.com Lorenzo Nirchi lorenzo.nirchi@alliedinst.com
Minnesota		Allied Instrumentation	844 398 9724 alliedinst.com Dawn Whittenburg dawnw@alliedinst.com
Mississippi		Techstar	832 599 3794 www.techstar.com Phil Deaton pdeaton@techstar.com
Missouri		Allied Instrumentation	(636) 343-0808 alliedinst.com Charlie Greise charlie.griese@alliedinst.com
Montana		Advanced Technical Sales, Inc.	406 861 2156 www.advancedtechmt.com Scott Nielsen scottn@advancedtechmt.com
Nebraska		Allied Instrumentation	844 398 9724 alliedinst.com Dawn Whittenburg dawnw@alliedinst.com
Nevada	Zip Codes: 894; 895; 897	Berthold Regional Sales Manager: Andrea Hand	865 483 1488 andrea.hand@berthold-us.com
	Zip Codes: 893; 898	Trailblazer Controls Corporation	801 955 3900 www.trailblazercontrols.com Jason Snedaker jason@trailblazercontrols.com
	Zip Codes: 889-891	Innovative Process Controls	702 759 0290 www.iprocessc.com David DiDomizio david@iprocessc.com
New Hampshire		Pond Technical Sales, Inc.	860 316 2570 www.pondtechnicalsales.com

		Bob Osnoe rosnoe@pondtechnicalsales.com
New Jersey	North East Technical Sales	215 513 1000 www.netechsals.com Doug Reichert dreichert@netechsals.com
New Mexico	Techstar	832 599 3794 www.techstar.com Phil Deaton pdeaton@techstar.com
New York	Empire Instruments LLC	716 304 6843 www.empireinst.com Bruno Albano brunoa@empireinst.com
	Zip Codes: 105; 120-149	
	Zip Codes: 100-104; 106- North East Technical Sales 119	215 513 1000 www.netechsals.com Doug Reichert dreichert@netechsals.com
North Carolina	System Service Corporation	704 226 1869 www.systemservicecorp.com Tim Good tgood@systemsservicecorp.com
North Dakota	Allied Instrumentation	844 398 9724 alliedinst.com Dawn Whittenburg dawnw@alliedinst.com
Ohio		
	Zip codes 430-432, 450- Henry M Wood 456	513 755 4060 www.henrymwood.com Brad Bossong brad@henrymwood.com
	Zip codes 433-438, 440- Regional Sales Manager: 449, 458, Harrison & Jefferson counties	Becky Olliges 865-323-0876 becky.olliges@berthold-us.com
	Zip code 457, Gallia, Lawrence, Monroe & Belmont counties	Hareza Technical Sales I (724) 942-4966 www.hareza.com Walt Lemunyon office@hareza.com
Oklahoma	Techstar	832 599 3794 www.techstar.com Phil Deaton pdeaton@techstar.com
Oregon	Engineered Process Controls	509 521 0971 www.engineeredprocess.com Greg LeBrun lebrun@engineeredprocess.com
Pennsylvania		
	zip codes 150-168	Hareza Technical Sales I (724) 942-4966 www.hareza.com

		Walt Lemunyon office@hareza.com
zip codes 169-196	North East Technical Sales	215 513 1000 www.netechsales.com Doug Reichert dreichert@netechsales.com
Rhode Island	Pond Technical Sales, Inc.	860 316 2570 www.pondtechnicalsales.com Bob Osnoe rosnoe@pondtechnicalsales.com
South Carolina	System Service Corporation	704 226 1869 www.systemservicecorp.com Tim Good tgood@systems-servicecorp.com
South Dakota	Allied Instrumentation	844 398 9724 alliedinst.com Dawn Whittenburg dawnw@alliedinst.com
Tennessee	Berthold Regional Sales Manager: Sam Arwood	865 274 9742 sam.arwood@berthold-us.com
Texas	Techstar	832 599 3794 www.techstar.com Phil Deaton pdeaton@techstar.com
Utah	Trailblazer Controls Corporation	801 955 3900 www.trailblazercontrols.com Jason Snedaker jason@trailblazercontrols.com
Vermont	Pond Technical Sales, Inc.	860 316 2570 www.pondtechnicalsales.com Bob Osnoe rosnoe@pondtechnicalsales.com
Virginia	PTLF Process, LLC	434 981 2515 www.ptlprocess.com John Douglas jdouglas@ptlprocess.com
Washington	Engineered Process Controls	425 686 4450 www.engineeredprocess.com Adam DiDomizio adam@engineeredprocess.com
West Virginia	Hareza Technical Sales I	(724) 942-4966 www.hareza.com Walt Lemunyon office@hareza.com
Wisconsin	Allied Instrumentation	219 924 4545 Alliedinst.com Lorenzo Nirchi lorenzo.nirchi@alliedinst.com

Wyoming

Zip Codes: 820-822; 824- Techstar
828

832 599 3794
www.techstar.com
Phil Deaton
pdeaton@techstar.com

Trailblazer Controls
Zip Codes: 823; 829-831 Corporation

801 955 3900
www.trailblazercontrols.com
Jason Snedaker
jason@trailblazercontrols.com

Canada

Coverage Area

Company Name

Phone/Webite

Canada

Vanko Analytics

780-433-0281
www.vanko.net
Tim Jones
timj@vanko.net

Steel Measurement Systems

Radiametrics, Inc.

440 245 9977
www.radiametrics.com
Kathy Huerner
kh@radiametrics.com

Mexico, Central and South America*

*All phone numbers should be preceded with 011 if calling from the US

Coverage Area

Company Name

Phone/Webite

Argentina, Bolivia, and Uruguay

Dastec SRL

54 11 5352 2500
www.dastecsrl.com.ar
Ignacio Calzadilla
icalzadilla@dastecsrl.com.ar

Brazil-Steel-Siderurgy-Metallurgy Industry Only

Vika Controls

5511 3159 3674
www.vikacontrols.com.br
Vinicius Kabakian
vk@vikacontrols.com.br

Brazil-All Industries except Steel-Siderurgy-Metallurgy Industry Only

Lince

21 2573-2344
www.instrumentos-lince.com.br
Marcel de Roure
marcel.roure@instrumentos-lince.com.br

Chile

ASPA

56 9 9442 5425

Boris Paéz
boris.paez@apsaspa.cl

Colombia, Dominican Republic, Ecuador, and Panama

Continental Process Instruments

57 1 467 3233
www.cpinstruments.com
Miguel Rodriquez
miguel.rodriquez@cpinstruments.com

Guatemala, El Salvador, Nicaragua, Honduras, Costa Rica, and Belize

Reset S.A

502 2315 0000
www.reseta.com
Julio Edgar Castillo Rojas
juliocastillo@reseta.com

Mexico

All Industries except Sugar

Medidores Industriales y Medicos S.A

52 55 5658 0766
www.medidores.com
Roberto Marquez Lopez

rmarquez@medidores.com

Sugar Industry

Biser

52 1 271 749 1030

biserindustrial.com

Sergio Martínez Hernández

s.martinez@biserindustrial.com

Peru

Marpatech Representaciones
S.A.C

511 224 9779

www.marvitech.com

Marco Paretto

mparetto@marpatech.com

**Venezuela, Trinidad, Tobago,
Caribbean (excluding the
Dominican Republic and Cuba),
and US Virgin Islands**

Renain

58 286 962 0628

www.renain.net

Reinaldo Nino

reinaldo.nino@renain.net

Pacific

Coverage Area

Company Name

Phone/Webite

**Guam, The Commonwealth of the
Northern Marianas Island, and the
Federated States of Micronesia**

Central Pacific Controls

808 450 2483

www.cpacifcc.com

David Didomizio

david@cpscifcc.com