



Berthold Technologies USA Representative Directory

United States

<i>Coverage Area</i>	<i>Company Name</i>	<i>Phone/Webite</i>
Steel Measurement Systems	Radiametrics, Inc.	440 245 9977 www.radiametrics.com Kathy Huerner kh@radiametrics.com
Alabama	Hatfield & Company	205-620-4000 www.hatfieldandcompany.com Cameron Poe cpoe@hatfieldandcompany.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.iprocess.com David DiDomizio david@iprocess.com
Colorado	Techstar	832 599 3794 www.techstar.com Phil Deaton pdeaton@techstar.com
Connecticut	Allied Instrumentation	207 378 1474 Alliedinst.com Bob Yale bob.yale@alliedinst.com
Delaware	North East Technical Sales	215 513 1000 www.netechsales.com Doug Reichert dreichert@netechsales.com
Florida Zip Code 324 & 325	Hatfield & Company	205-620-4000 www.hatfieldandcompany.com Cameron Poe cpoe@hatfieldandcompany.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 Didmomio 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	zip codes 461, 462, 470-472, 474-478	Henry M Wood	513 755 4060 www.henrymwood.com Eric Yaeger eric@henrymwood.com
	Zip Codes: 460, 463 - 469, 473, 479	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	All except zip codes 411, 412, 415, 416	Henry M Wood	513 755 4060 www.henrymwood.com Eric Yaeger eric@henrymwood.com
	Zip codes 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		Allied Instrumentation	207 378 1474 Alliedinst.com Bob Yale bob.yale@alliedinst.com
Maryland	zip codes 215 & 217	Hareza Technical Sales	(724) 942-4966 www.hareza.com Walt Lemunyon office@hareza.com
	all zip codes except 215 & 217	North East Technical Sales	215 513 1000 www.netechsales.com Doug Reichert dreichert@netechsales.com

Massachusetts	Allied Instrumentation	207 378 1474 Alliedinst.com Bob Yale bob.yale@alliedinst.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	Regional Sales Manager:	Mark Klee 865 274 9742 mark.klee@berthold-us.com
Zip Codes: 894; 895; 897	Trailblazer Controls Corporation	801 955 3900 www.trailblazercontrols.com Jason Snedaker jason@trailblazercontrols.com
Zip Codes: 893; 898	Innovative Process Controls	702 759 0290 www.iprocessc.com David DiDomizio david@iprocessc.com
Zip Codes: 889-891		
New Hampshire	Allied Instrumentation	207 378 1474 Alliedinst.com Bob Yale bob.yale@alliedinst.com
New Jersey	North East Technical Sales	215 513 1000 www.netechsales.com Doug Reichert dreichert@netechsales.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; 106-119	North East Technical Sales	215 513 1000 www.netechsales.com Doug Reichert dreichert@netechsales.com
North Carolina		System Service Corporation	704 226 1869 www.systemservicecorp.com Tim Good tgood@systemservicecorp.com
North Dakota		Allied Instrumentation	844 398 9724 alliedinst.com Dawn Whittenburg dawnw@alliedinst.com
Ohio	430-433, 435 Defiance Co., 437 except Belmont, 438, 446 Wayne and Holmes Co,	Henry M Wood	513 755 4060 www.henrymwood.com Brad Bossong brad@henrymwood.com
	Zip codes 434-436, except Defiance Counties, 448-449	Allied Instrumentation	219 924 4545 Alliedinst.com Lorenzo Nirchi lorenzo.nirchi@alliedinst.com
	437 Belmont, 439-445, 446 Carroll & Tuscarawas, 447, 457	Hareza Technical Sales	(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	(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		Allied Instrumentation	207 378 1474 Alliedinst.com Bob Yale bob.yale@alliedinst.com
South Carolina		System Service Corporation	704 226 1869 www.systemservicecorp.com Tim Good tgood@systemservicecorp.com
South Dakota		Allied Instrumentation	844 398 9724 alliedinst.com Dawn Whittenburg dawnw@alliedinst.com

Tennessee

Zip codes 370-374, 376-379, 384-385

Process Supplies (PSA)

865-271-1093

www.processsupplies.com/

John Higdon

jhigdon@processsupplies.com

380-383

Regional Sales Manager:

Becky Olliges

becky.olliges@berthold-us.com

865 323 0876

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

Allied Instrumentation

207 378 1474

Alliedinst.com

Bob Yale

bob.yale@alliedinst.com

Virginia

PTLF Process, LLC

434 981 2515

www.ptlfprocess.com

John Douglas

jdouglas@ptlfprocess.com

Washington

Engineered Process Controls

425 686 4450

www.engineeredprocess.com

Adam DiDomizio

adam@engineeredprocess.com

West Virginia

Hareza Technical Sales

(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; 824-828

Techstar

832 599 3794

www.techstar.com

Phil Deaton

pdeaton@techstar.com

Zip Codes: 823; 829-831

Trailblazer Controls Corporation

801 955 3900

www.trailblazercontrols.com

Jason Snedaker

jason@trailblazercontrols.com