

SOURCE LIST

Company	Phone Number	International Phone	Website
Ambion (Applied Biosystems)	800-888-8804	See website	www.ambion.com OR www.appliedbiosystems.com
Amersham Biosciences (GE Healthcare Life Sciences)	800-526-3593	See website	www.amershambiosciences.com
Barnstead International	800-553-0039	See website	www.barnstead.com
BioTek Instruments, Inc.	888-451-5171	See website	www.biotek.com
ColeParmer	800-323-4340	See website	www.coleparmer.com
Corning (Costar)	607-974-9000	1-607-974-9000	www.corning.com/lifesciences
Dexter Magnetic Technologies	See website	See website	www.lifeseq.com
Eppendorf	800-645-3050	See website	www.eppendorfn.com
Fisher Scientific	800-766-7000	See website	www.fishersci.com
Invitrogen	800-955-6288	See website	www.invitrogen.com
Millipore	800-645-5476	See website	www.millipore.com
MJ Research (Bio-Rad)	200-424-6723	See website	www.mjr.com
Molecular Probes (Invitrogen)	800-955-6288	See website	probes.invitrogen.com
MOSS Inc.	800-932-6677	See website	www.mosssubstrates.com
New England Biolabs	800-632-5227	See website	www.neb.com
Pall Life Sciences	800-521-1520	See website	www.pall.com
PerkinElmer	800-762-4000	See website	www.perkinelmer.com
Pierce	800-874-3723	See website	www.piercenet.com
Prozyme	800-457-9444	See website	www.prozyme.com
Rainin	800-472-4646	See website	www.rainin.com
Roche Applied Science	800-428-5433, Ext. 14659	See website	www.roche-applied-science.com (see Industrial Products)
Samco Scientific	800-522-3359	1-818-838-2400	www.samcosci.com
Sigma	800-325-3010	See website	www.sigmaaldrich.com
Qiagen	800-426-8157	See website	www.qiagen.com
USA Scientific	800-522-8477	1-352-237-6288	www.usascientific.com
VWR	800-932-5000	See website	www.vwr.com