



- Rep - RSM**
- Abbey Mec-tric - SA
 - Advanced Technical Sales- AH
 - Advanced Technology Associates- BO
 - Allied Instrumentation
 - Central Pacific Controls- AH
 - Empire- BO
 - Engineered Process Controls - AH
 - Henry M Wood- BO
 - Hile Controls - SA
 - Innovative Process Controls- AH
 - North East Technical Sales- BO
 - Open- AH
 - Hareza Technical Sales - BO
 - Open - SA
 - PTLF Process- BO
 - Pond- BO
 - System Services- SA
 - Techstar- SA
 - Trailblazer- AH
 - Techstar - AH
 - Open - BO