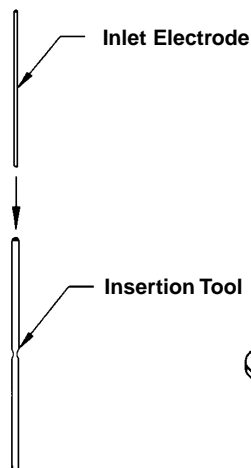
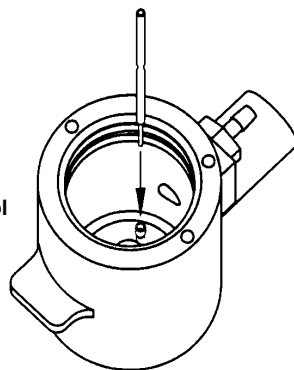


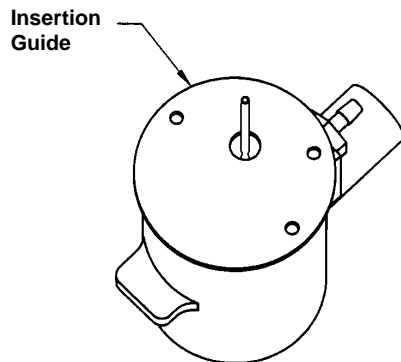
## Instructions, Inlet Electrode Installation



1. Push the Inlet Electrode all the way into the Inlet Electrode Insertion Tool.



2. Push the Inlet Electrode all the way into the Socket.



3. Use the Inlet Electrode Insertion Guide to position the Electrode. Remove the Inlet Electrode Insertion Tool before removing the Inlet Electrode Insertion Guide.