

**A NOTE FOR CUSTOMERS PLANNING TO PERFORM DGGE, CDGE, TGTGE OR OTHER APPLICATIONS REQUIRING USE OF THE HEATER:**

**This cooling tank is not recommended for heating applications. To ensure best results, please order the standard DCode electrophoresis tank.**

<b>Catalog number</b>	<b>Product description</b>
<b>170-9222</b>	<b>DCode electrophoresis tank</b>