

## Product Information

<b>Description:</b>	This kit contains all the components required for Human iPSC Derived Cardiomyocytes (BPS Bioscience #78529) thaw.
<b>Formulation:</b>	<ul style="list-style-type: none"><li>• <b>Component A:</b> Maintenance Medium C17 (BPS Bioscience #78509)</li><li>• <b>Component B:</b> Serum Replacement (BPS Bioscience #82245)</li><li>• <b>Component C:</b> 10 mM Thiazovivin (BPS Bioscience #82246)</li></ul>
<b>Quality:</b>	Mycoplasma-free. All components are low endotoxin.
<b>Storage:</b>	Upon receipt, Thaw Medium Kit C17 should be stored in the dark at -20°C.
<b>Stability:</b>	When stored in the dark at -20°C, the product is stable for six months. Component A can be stored at 4°C for a month after thaw. Thaw components at 4°C or room temperature (do not use a water bath at 37°C to thaw). Avoid repeated freeze/thaw cycles.
<b>Preparation:</b>	Thaw Medium C17 (prepare immediately before use): add 5 ml of Component B and 50 µl of 10 mM of Component C to 45 ml of Component A. Prepare under aseptic conditions. Do not filter.
<b>Culture Protocol:</b>	For detailed thawing and cell culture protocol, see Human iPSC Derived Cardiomyocytes (BPS Bioscience #78529) datasheet.