








ExpiCHO and Expi293 System shaking speed recommendations

Maximize your protein yield by following these guidelines

Small scale			Large scale
125 mL–2 L flasks Recommended shaking speeds: 125 ± 5 rpm (19 mm orbital diameter) 120 ± 5 rpm (25 mm orbital diameter) 95 ± 5 rpm (50 mm orbital diameter)			2.8–3 L flasks Recommended shaking speeds: 75 ± 5 rpm (19 mm orbital diameter) 80 ± 5 rpm (25 mm orbital diameter) 80 ± 5 rpm (50 mm orbital diameter)
 125 mL	 250 mL		
<p>Culture volume to transfect 100 mL</p> <p>Ratio of medium to flask volume 1:3.6</p> <p>Final culture volume ~140 mL</p>	<p>Culture volume to transfect 200 mL</p> <p>Ratio of medium to flask volume 1:3.6</p> <p>Final culture volume ~280 mL</p>	<p>Culture volume to transfect 400 mL</p> <p>Ratio of medium to flask volume 1:3.6</p> <p>Final culture volume ~560 mL</p>	<p>Culture volume to transfect 600–750 mL</p> <p>Ratio of medium to flask volume 1:3.0–3.5</p> <p>Final culture volume ~850 mL–1 L</p>
 500 mL	 1 L	 2 L	 2.8–3 L
 Vented, nonbaffled for all volumes			

Find out more at thermofisher.com/proteinexpression

ThermoFisher
SCIENTIFIC