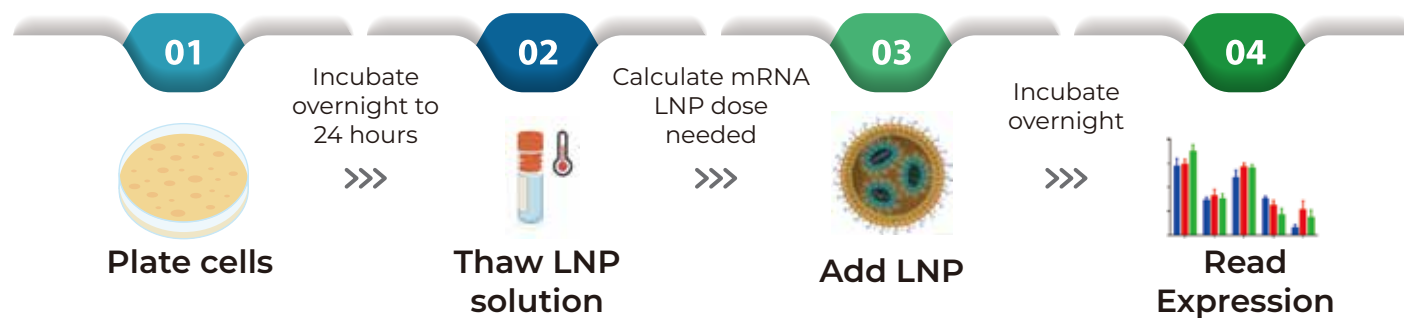


mRNA LNP *in vitro* Cell Transfection

1. LNP can be added directly to cells incubated in complete media
2. Cells grown and transfected in complete media (DMEM +10% FBS) has higher expression than reduced serum medium
3. For suspension cell lines, we recommend using low fluorescent background media such as Flourobrite® with 10% FBS, 1X GlutaMax® and 1X sodium pyruvate
4. For adherent cell line, we recommend replace complete media with Opti-mem right before reading the cell eGFP expression
5. For higher expression, we recommend incubate the LNP with FBS at 1:2 volume ratio (LNP: FBS) for 30 minutes at 37C before transfecting



[Learn more - www.genscript.com/LNP](http://www.genscript.com/LNP)