



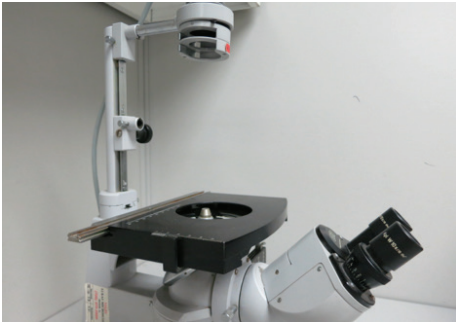
How do I visualize my cells in a Nunc Cell Factory system?

The Thermo Scientific™ Nunc™ Cell Factory™ systems are a proven solution for the cultivation of adherent cells. An inverted microscope is the perfect solution for monitoring the condition of your culture.

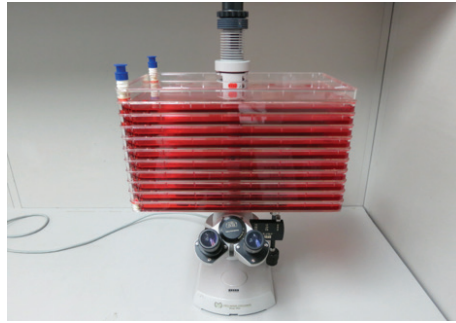
To help ensure success, the microscope must have a flat stage. When viewing larger formats such as a Cell Factory 10-layer system, a removable light source may be necessary.



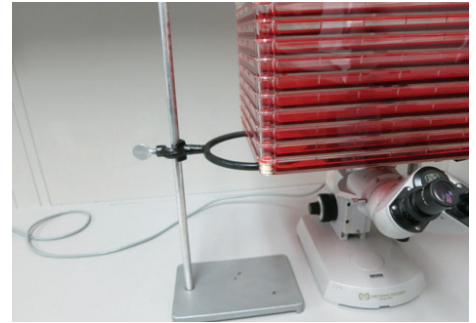
Process for visualizing and monitoring and monitor cells in the Nunc Cell Factory system



Begin by removing all stage accessories to provide a flat surface for the Cell Factory system to rest upon. If additional space is required, remove the light source.

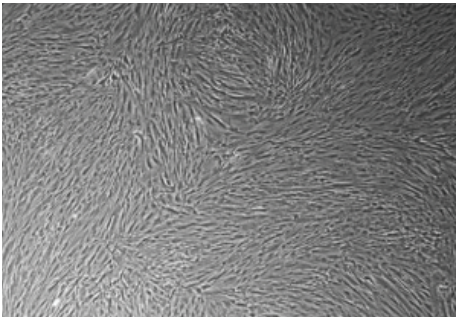


Place the Cell Factory system on the stage with the desired location to be visualized over the objective lens. If necessary, replace the light source or place an alternative light source above the Cell Factory system.

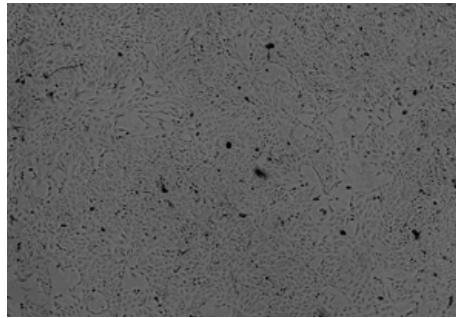


If the microscope stage is not capable of supporting the weight of the Cell Factory system, it is advisable to use an adjustable stand to support some of the weight of the Cell Factory system.

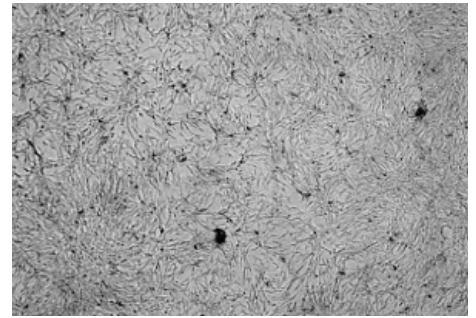
Photographs of cells cultured in a Cell Factory system viewed with an inverted microscope.



95% confluent MRC-5 cells
MRC-5 cells are a human diploid cell line commonly used in the manufacturing of vaccines for hepatitis A, rubella, varicella, and rabies.



80-90% confluent Vero cells
Vero cells are derived from the kidney of an African green monkey and are commonly used for the production of vaccines for rotavirus, smallpox, and inactivated poliovirus.



60-70% confluent chicken embryo fibroblast (CEF) cells
CEFs are primary cells used in the production of vaccines for measles, mumps, rabies, and influenza. They are also commonly used in the manufacturing of avian vaccines.

Find out more at thermofisher.com/cellfactory