

Product Information

Description:	NK Cell Basal Medium is optimized to expand human natural killer (NK) cells, using BPS Bioscience Growth-Arrested NK Feeder Cells (#78912).
Formulation:	Contains Heat Inactivated Fetal Bovine Serum (HiFBS) and other supplements. Growth factors required for the optimal expansion of NK cells have not been added to NK Cell Basal Medium.
Quality:	Sterile filtered (0.2 μ m) and Mycoplasma-free. All components are low endotoxin. FBS is ISIA-certified and sourced from the United States of America.
Storage:	Upon receipt, the medium should be stored in the dark at 4°C.
Stability:	When stored in the dark at 4°C, the product is stable for six months after the first thaw. Avoid repeated freeze/thaw cycles.

Directions for Use:

1. Add desired cytokines, growth factors, or other components to NK Cell Basal Medium. Mix thoroughly. Set up cultures as desired.

Note: The use of Growth-Arrested NK Feeder Cells (#78912) is highly recommended.