



Suggested Tips for Running 41580 with Cell Lysates

Sample Preparation:

For measuring hIFN γ R1 in cell culture:

All cell culture samples should be quick frozen and stored below -20°C until being tested.

Measuring hIFN γ R1 of adherent cells:

- Remove cell culture supernatant and wash cells twice with cold PBS.
- Add Sample Diluent to adhered cells at 10 ml per T75 flask or at desired volume. Use a cell scraper to dislodge the cells.
- For accurate quantitation of total cellular hIFN γ R1, check under a microscope to confirm that all cells are dislodged from the flask.
- Incubate the cell suspension on ice for 20 minutes. If necessary, break up cellular clusters by gently pipetting up and down.
- Centrifuge the solution at 500 X g for 15 minutes and collect the supernatant (cell lysate solution).
- Add 100 μ l of the lysate to designated wells on the ELISA plate. Cell lysate may be diluted with Sample Diluent as desired to within the assay range of the ELISA.

Measuring hIFN γ R1 of suspension cells:

- Centrifuge cell suspension at 500 X g for 5 minutes. Wash the cells twice with cold PBS.
- Suspend cells in Sample Diluent at 10 ml per 30,000,000 cells or at desired volume, and incubate on ice for 20 minutes.
- Centrifuge the solution at 500 X g for 15 minutes and collect the supernatant (cell lysate solution).
- Add 100 μ l of the lysate to designated wells on the ELISA plate. Cell lysate may be diluted with Sample Diluent as desired to within the assay range of the ELISA.

