



# Human IFN Alpha **All Subtype** ELISA Kit, High Sensitivity

for Tissue Culture Media (TCM)

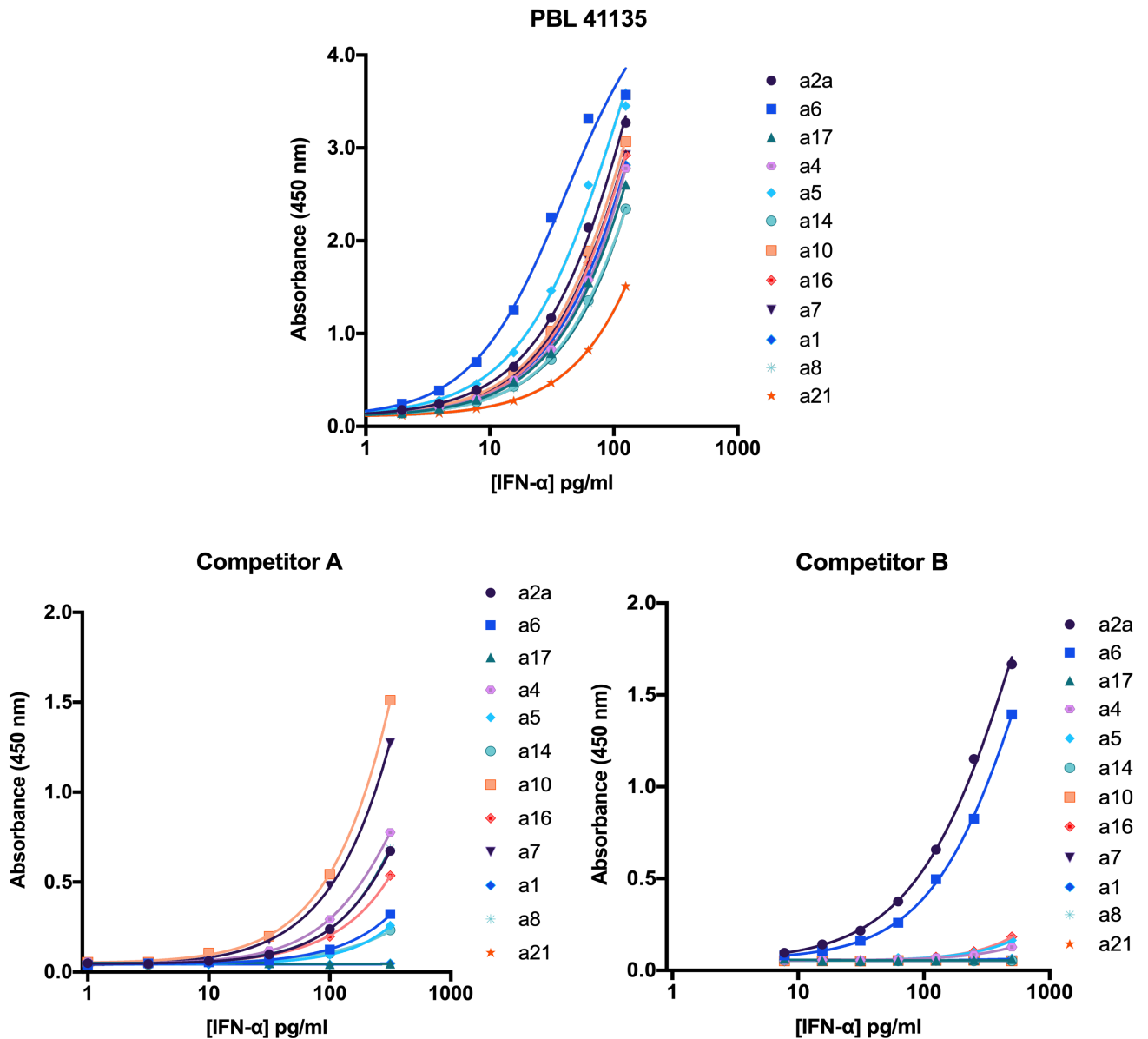
- Most sensitive ELISA on the market for global IFN- $\alpha$  detection in tissue culture media (TCM)
- Low pg/ml detection of **12 of 12 human IFN- $\alpha$  subtypes**
- Assay standard range of **1.95-125 pg/ml**
- Reproducible results with < 10% inter- and < 8% intra-assay CVs

PBL's **VeriKine-HS Human Interferon Alpha All Subtype ELISA (Catalog No. 41135)** features highly sensitive detection and quantification of all 12 human IFN- $\alpha$  subtypes and provides robust results in tissue culture media (TCM).

**Figure 1.** LLOQ and LLOD in Tissue Culture Media (TCM)

Subtype	LLOQ	LLOD
$\alpha 1$ ( $\alpha D$ )	1.17	0.36
$\alpha 2a$ ( $\alpha A$ )	1.22	0.37
$\alpha 4a$ ( $\alpha M1$ )	1.76	0.53
$\alpha 5$ ( $\alpha G$ )	1.14	0.36
$\alpha 6$ ( $\alpha K$ )	0.73	0.25
$\alpha 7$ ( $\alpha J1$ )	1.11	0.35
$\alpha 8$ ( $\alpha B2$ )	1.51	0.46
$\alpha 10$ ( $\alpha C$ )	1.89	0.61
$\alpha 14$ ( $\alpha H$ )	2.92	0.97
$\alpha 16$ ( $\alpha WA$ )	2.29	0.77
$\alpha 17$ ( $\alpha I$ )	1.97	0.62
$\alpha 21$ ( $\alpha F$ )	2.88	0.93

**Figure 2.** Comparison of IFN- $\alpha$  Subtypes Analyses (TCM)



- Both competitors' kits quantitate some, but **not all**, of the 12 IFN- $\alpha$  subtypes
- High sensitivity allows for **early detection** of low level IFN- $\alpha$  subtypes in your study

Visit our website for more information: [www.pbl assaysci.com](http://www.pbl assaysci.com)