



Human Cell Expressed (HCE) Cytokines and Growth Factors

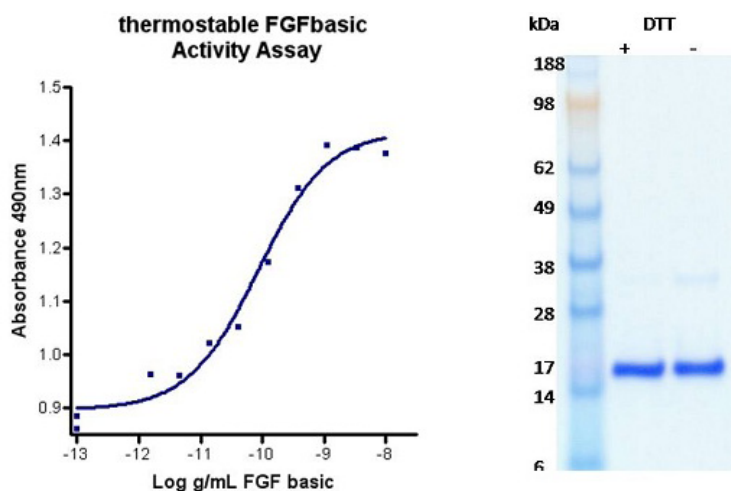
As ***Human*** As You Can Get

- Authentic human glycosylation & native human post-translational modifications
- Animal-component-free, protein-free production
- High bioactivity and stability; lyophilized
- Heterodimeric cytokines are formed like endogenous versions
- Highly pure, low endotoxin proteins

PBL Assay Science offers authentic recombinant, human cell expressed (HCE) cytokines and growth factors which closely mimic native human proteins. These affordable, tag-free, animal-component-free proteins are produced in stable, serum-free, chemically defined cell culture helping ensure lot-to-lot consistency and reproducibility.

These HCE cytokines retain native properties and possess authentic human N- and/or O-linked glycosylation and other post-translational modifications preserving appropriate biological activities and epitopes.

Figures 1 & 2. FGF Basic-TS (thermostable) Activity Assay & SDS-PAGE



Product Information for Human Cell Expressed (HCE) Cytokines and Growth Factors

Fibroblast Growth Factors (FGFs)

11631	Human FGF Basic-TS (thermostable)	10, 50, 1000 µg
11632	Human FGF-4	10, 100, 1000 µg
11633	Human FGF-7 (Keratinocyte Growth Factor, KGF)	10, 100, 1000 µg
11634	Human FGF-8b	10, 100, 1000 µg

Hematopoietin and IL Family

11600	Human Serum Albumin (HSA)	100, 1000 µg
11601	Human IL-1 beta	10, 100, 1000 µg
11602	Human IL-2	10, 100, 1000 µg
11603	Human IL-3	10, 100, 1000 µg
11604	Human IL-4	10, 100, 1000 µg
11605	Human IL-6	10, 100, 1000 µg
11606	Human IL-7	10, 100, 1000 µg
11607	Human IL-9	10, 100, 1000 µg
11608	Human IL-10	10, 100, 1000 µg
11609	Human IL-12	10, 100, 1000 µg
11610	Human IL-17A	10, 100, 1000 µg
11611	Human IL-17F	10, 100, 1000 µg
11612	Human IL-23	10, 100, 1000 µg
11613	Human IL-27	10, 100, 1000 µg
11614	Human Erythropoietin (EPO)	10, 100, 1000 µg
11615	Human Thrombopoietin (TPO)	10, 100, 1000 µg
11616	Human G-CSF	10, 100, 1000 µg
11617	Human GM-CSF	10, 100, 1000 µg
11618	Human Hepatocyte Growth Factor (HGF)	10, 100, 1000 µg
11619	Human Growth Hormone (HGH)	10, 100, 1000 µg

TGF Beta Superfamily

11621	Human Activin A	2, 10, 100, 1000 µg
11622	Human BMP-2	10, 100, 1000 µg
11623	Human BMP-4	10, 100, 1000 µg
11624	Human BMP-7	10, 100, 1000 µg
11625	Human LEFTY-1	10, 100, 1000 µg
11627	Human TGF β1	5, 100, 1000 µg
11628	Human TGF β2	5, 100, 1000 µg
11629	Human TGF β3	5, 100, 1000 µg

Tyrosine Kinase Receptor Family

11640	Human FLT3 Ligand	10, 100, 1000 µg
11641	Human M-CSF	10, 100, 1000 µg
11642	Human Stem Cell Factor (SCF)	10, 100, 1000 µg
11643	Human VEGF121	10, 100, 1000 µg
11644	Human VEGF165	10, 100, 1000 µg
11645	Human PDGF-AA	10, 100, 1000 µg
11646	Human Leukemia Inhibitory Factor (LIF)	10, 100, 1000 µg

Other HCE Proteins

11650	TNF α	10, 100, 1000 µg
11660	Human Cystatin C	10, 100, 1000 µg
11661	Human Oncostatin M (OSM)	10, 100, 1000 µg
11662	Human Pro-IGF-II	10, 100, 1000 µg
11663	Human Pleiotrophin (PTN)	10, 100, 1000 µg
11670	Human β NGF	10, 100, 1000 µg
11671	Human Noggin	10, 100, 1000 µg