

Research Use Only (RUO) Growth Factors & Cytokines for Cell-Based Applications

Sartorius provides a broad range of high quality RUO growth factors and cytokines which are produced using recombinant DNA technology and do not contain any animal-derived components or contaminants. These products are produced and distributed according to a Quality Management System that is certified for compliance with ISO 9001.

Features

- High quality and extensively validated reagents ensure reproducibility and lot-to-lot consistency
- Carrier protein-free
- Highest levels of purity and bioactivity
- Animal-free
- Every lot tested to confirm activity
- Low endotoxin levels

RUO Growth Factor and Cytokine Products

SKU	Description	Unit of Measure (µg)
CYK-0100-0001	Recombinant human EGF protein	100
CYK-0500-0001	Recombinant human EGF protein	500
CYK-0050-0002	Recombinant FGF2-G3 protein	50
CYK-0100-0002	Recombinant FGF2-G3 protein	100
CYK-0500-0002	Recombinant FGF2-G3 protein	500
CYK-0010-0003	Recombinant human KGF (FGF-7) protein	10
CYK-0025-0003	Recombinant human KGF (FGF-7) protein	25
CYK-0050-0003	Recombinant human KGF (FGF-7) protein	50
CYK-0500-0003	Recombinant human KGF (FGF-7) protein	500
CYK-0100-0004	Recombinant human IGF-1 LR3 protein	100
CYK-0500-0004	Recombinant human IGF-1 LR3 protein	500
CYK-0010-0005	Recombinant human VEGF 165 protein	10
CYK-0025-0005	Recombinant human VEGF 165 protein	25
CYK-0100-0005	Recombinant human VEGF 165 protein	100
CYK-0500-0005	Recombinant human VEGF 165 protein	500
CYK-0025-0006	Recombinant human BMP-4 protein	25
CYK-0100-0006	Recombinant human BMP-4 protein	100
CYK-0025-0007	Recombinant human Activin A protein	25
CYK-0100-0007	Recombinant human Activin A protein	100



RUO Growth Factor and Cytokine Products (Continued)

SKU	Description	Unit of Measure (µg)
CYK-0500-0007	Recombinant human Activin A protein	500
CYK-0010-0008	Recombinant human BDNF protein	10
CYK-0025-0008	Recombinant human BDNF protein	25
CYK-0100-0008	Recombinant human BDNF protein	100
CYK-0500-0008	Recombinant human BDNF protein	500
CYK-0025-0009	Recombinant human GDNF protein	25
CYK-0100-0009	Recombinant human GDNF protein	100
CYK-0500-0009	Recombinant human GDNF protein	500
CYK-0025-0010	Recombinant human TGF-β1 PLUS protein	25
CYK-0050-0010	Recombinant human TGF-β1 PLUS protein	50
CYK-0100-0010	Recombinant human TGF-β1 PLUS protein	100
CYK-0025-0011	Recombinant human PDGF-AA protein	25
CYK-0100-0011	Recombinant human PDGF-AA protein	100
CYK-0025-0012	Recombinant human LIF protein	25
CYK-0100-0012	Recombinant human LIF protein	100
CYK-0025-0013	Recombinant human BMP-2 protein	25
CYK-0100-0013	Recombinant human BMP-2 protein	100
CYK-0025-0014	Recombinant human NRG-1 protein	25
CYK-0100-0014	Recombinant human NRG-1 protein	100
CYK-0500-0014	Recombinant human NRG-1 protein	500
CYK-0025-0015	Recombinant human TGF-β3 PLUS pr	25
CYK-0050-0015	Recombinant human TGF-β3 PLUS protein	50
CYK-0100-0015	Recombinant human TGF-β3 PLUS protein	100
CYK-0025-0016	Recombinant human Noggin protein	25
CYK-0100-0016	Recombinant human Noggin protein	100
CYK-0500-0016	Recombinant human Noggin protein	500
CYK-0025-0017	Recombinant human FGF-10 protein	25
CYK-0100-0017	Recombinant human FGF-10 protein	100
CYK-0500-0017	Recombinant human FGF-10 protein	500
CYK-0025-0018	Recombinant human FGF-9 protein	25
CYK-0100-0018	Recombinant human FGF-9 protein	100
CYK-0025-0019	Recombinant human R-spondin 1 protein	25
CYK-0100-0019	Recombinant human R-spondin 1 protein	100
CYK-0500-0019	Recombinant human R-spondin 1 protein	500
CYK-0025-0020	Recombinant human NT-3 protein	25
CYK-0050-0020	Recombinant human NT-3 protein	50
CYK-0100-0020	Recombinant human NT-3 protein	100
CYK-0025-0021	Recombinant human HGF (NK1) protein	25
CYK-0100-0021	Recombinant human HGF (NK1) protein	100
CYK-0025-0022	Recombinant human OSM protein	25
CYK-0100-0022	Recombinant human OSM protein	100
CYK-0100-0023	Recombinant human FGF-2 protein 154 aa	100

RUO Growth Factor and Cytokine Products (Continued)

SKU	Description	Unit of Measure (µg)
CYK-0500-0023	Recombinant human FGF-2 protein 154 aa	500
CYK-0025-0024	Recombinant human SHH protein	25
CYK-0100-0024	Recombinant human SHH protein	100
CYK-0025-0025	Recombinant human FGF-4 protein	25
CYK-0100-0025	Recombinant human FGF-4 protein	100
CYK-0025-0026	Recombinant human R-spondin 3 protein	25
CYK-0100-0026	Recombinant human R-spondin 3 protein	100
CYK-0025-0027	Recombinant human FGF-8a protein	25
CYK-0100-0027	Recombinant human FGF-8a protein	100
CYK-0025-0028	Recombinant human FGF-8b protein	25
CYK-0100-0028	Recombinant human FGF-8b protein	100
CYK-0025-0030	Recombinant human G-CSF protein	25
CYK-0100-0030	Recombinant human G-CSF protein	100
CYK-0025-0031	Recombinant human M-CSF protein	25
CYK-0100-0031	Recombinant human M-CSF protein	100
CYK-0025-0032	Recombinant human PDGF-BB protein	25
CYK-0100-0032	Recombinant human PDGF-BB protein	100
CYK-0025-1001	Recombinant human Interleukin-2	25
CYK-0100-1001	Recombinant human Interleukin-2	100
CYK-0025-1002	Recombinant human Stem Cell Factor	25
CYK-0100-1002	Recombinant human Stem Cell Factor	100
CYK-0025-1003	Recombinant human Interleukin-15	25
CYK-0100-1003	Recombinant human Interleukin-15	100
CYK-0025-1004	Recombinant human Tumour Necrosis Factor Alpha	25
CYK-0100-1004	Recombinant human Tumour Necrosis Factor Alpha	100
CYK-0025-1005	Recombinant human Fms-like tyrosine kinase 3 ligand	25
CYK-0100-1005	Recombinant human Fms-like tyrosine kinase 3 ligand	100
CYK-0025-1006	Recombinant human Interleukin-4	25
CYK-0100-1006	Recombinant human Interleukin-4	100
CYK-0025-1007	Recombinant human Interleukin-10	25
CYK-0100-1007	Recombinant human Interleukin-10	100
CYK-0025-1008	Recombinant human Interleukin-7	25
CYK-0100-1008	Recombinant human Interleukin-7	100
CYK-0025-1009	Recombinant human Interleukin-12	25
CYK-0100-1009	Recombinant human Interleukin-12	100
CYK-0100-1010	Recombinant human Vitronectin	100
CYK-0025-1011	Recombinant human Interleukin-3	25
CYK-0100-1011	Recombinant human Interleukin-3	100
CYK-0025-1012	Recombinant human Thrombopoietin	25
CYK-0100-1012	Recombinant human Thrombopoietin	100
CYK-0025-1013	Recombinant human Erythropoietin	25
CYK-0100-1013	Recombinant human Erythropoietin	100

RUO Growth Factor and Cytokine Products (Continued)

SKU	Description	Unit of Measure (µg)
CYK-0025-1014	Recombinant human Interleukin-6	25
CYK-0100-1014	Recombinant human Interleukin-6	100
CYK-0025-1015	Recombinant human Interleukin-18	25
CYK-0100-1015	Recombinant human Interleukin-18	100
CYK-0025-1016	Recombinant human Interleukin-1 Beta	25
CYK-0100-1016	Recombinant human Interleukin-1 Beta	100
CYK-0025-1017	Recombinant human Interferon gamma	25
CYK-0100-1017	Recombinant human Interferon gamma	100
CYK-0025-1018	Recombinant human Interleukin-21	25
CYK-0100-1018	Recombinant human Interleukin-21	100
CYK-0025-1019	Recombinant human Granulocyte-macrophage colony-stimulating factor	25
CYK-0100-1019	Recombinant human Granulocyte-macrophage colony-stimulating factor	100

Germany:

Sartorius Lab Instruments
Phone +49 551 308 0

USA:

Sartorius Corporation
Phone +1 631 254 4249
Toll-free +1 800 635 2906



Visit our online store: www.sartorius.com/shop



For questions, email: