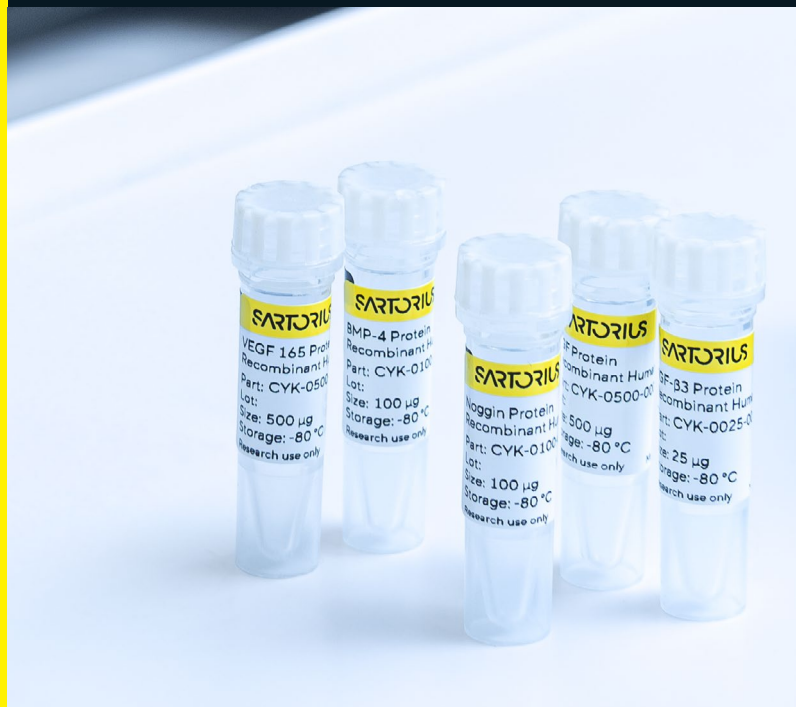


RUO Growth Factors & Cytokines



Reconstitution Instructions

1. Refer to the table below for specific reconstitution solution recommendations for your product.
2. Briefly centrifuge to make sure all the powder in the vial is at the bottom of the vial before opening.
3. Add the correct reconstitution solution to give a concentration of 0.5 – 1 mg/mL .
4. For larger sample sizes, reconstitute in ~75% of the total volume and then leave for 1 minute before transferring to a fresh tube, rinsing the original tube with the remaining solution before pooling the samples together.
5. Allow to stand for a minute at room temperature before gently agitating the mixture while being careful not to create foam – **do not** vortex.
6. Your stock solution is ready and can now be made into single-use or stock aliquots. At this point forward, all further dilutions should be made with water or PBS. It is advised that if you are storing at a lower concentration than <math>< 50 \mu\text{g/mL}</math>, if compatible with your work to add a carrier protein (all our products are carrier protein-free).
7. To avoid freeze-thaw cycles which can reduce the bioactivity of the protein and damage it, it is advisable to store in single-use aliquots and freeze between $-20\text{ }^{\circ}\text{C}$ and $-80\text{ }^{\circ}\text{C}$.

Technical Specifications

Product Name	SKU Number(s)	From	Recommended Reconstitution Buffer
Recombinant human EGF protein	CYK-0100-0001 CYK-0500-0001	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant FGF2-G3 protein	CYK-0050-0002 CYK-0100-0002 CYK-0500-0002	Lyophilized from Tris/NaCl/CyS/mannitol	Water
Recombinant human KGF (FGF-7) protein	CYK-0010-0003 CYK-0025-0003 CYK-0050-0003 CYK-0500-0003	Lyophilized from PBS	Water
Recombinant human IGF-1 LR3 protein	CYK-0100-0004 CYK-0500-0004	Lyophilized from acetonitrile/TFA	Water
Recombinant human VEGF 165 protein	CYK-0010-0005 CYK-0025-0005 CYK-0100-0005 CYK-0500-0005	Lyophilized from acetonitrile/TFA	Water
Recombinant human BMP-4 protein	CYK-0025-0006 CYK-0100-0006	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human activin A protein	CYK-0025-0006 CYK-0100-0006	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human BDNF protein	CYK-0010-0008 CYK-0025-0008 CYK-0100-0008 CYK-0500-0008	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human GDNF protein	CYK-0025-0009 CYK-0100-0009 CYK-0500-0009	Lyophilized from acetonitrile/TFA	Water
Recombinant human TGF- β 1 PLUS protein	CYK-0025-0010 CYK-0050-0010 CYK-0100-0010	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human PDGF-AA protein	CYK-0025-0011 CYK-0100-0011	Lyophilized from acetonitrile/TFA	Water
Recombinant human LIF protein	CYK-0025-0012 CYK-0100-0012	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human BMP-2 protein	CYK-0025-0013 CYK-0100-0013	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human NRG-1 protein	CYK-0025-0014 CYK-0100-0014 CYK-0500-0014	Lyophilized from acetonitrile/TFA	Water
Recombinant human TGF- β 3 protein	CYK-0025-0015 CYK-0050-0015 CYK-0100-0015	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human Noggin protein	CYK-0025-0016 CYK-0100-0016 CYK-0500-0016	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human FGF-10 protein	CYK-0025-0017 CYK-0100-0017 CYK-0500-0017	Lyophilized from PBS	Water
Recombinant human FGF-9 protein	CYK-0025-0018 CYK-0100-0018	Lyophilized from PBS	Water
Recombinant human R-spondin 1 protein	CYK-0025-0019 CYK-0100-0019 CYK-0500-0019	Lyophilized from acetonitrile/TFA	10 mM HCl

Product Name	SKU Number(s)	From	Recommended Reconstitution Buffer
Recombinant human NT-3 protein	CYK-0025-0020 CYK-0050-0020 CYK-0100-0020	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human HGF protein	CYK-0025-0021 CYK-0100-0021	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human OSM protein	CYK-0025-0022 CYK-0100-0022	Lyophilized from acetonitrile/TFA	Water
Recombinant human FGF-2 protein (154 aa)	CYK-0100-0023 CYK-0500-0023	Lyophilized from Tris/NaCl/cysteine/mannitol	NA
Recombinant human Shh protein	CYK-0025-0024 CYK-0100-0024	Lyophilized from PBS	Water
Recombinant human FGF-4 protein	CYK-0025-0025 CYK-0100-0025	Lyophilized from PBS	Water
Recombinant human R-spondin 3 protein	CYK-0025-0026 CYK-0100-0026	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human FGF-8a protein	CYK-0025-0027 CYK-0100-0027	Lyophilized from PBS	Water
Recombinant human FGF-8b protein	CYK-0025-0028 CYK-0100-0028	Lyophilized from HEPES/NaCl/mannitol	Water
Recombinant human G-CSF protein	CYK-0025-0030 CYK-0100-0030	Lyophilized from Tris/mannitol	Water
Recombinant human M-CSF protein	CYK-0025-0031 CYK-0100-0031	Lyophilized from acetonitrile/TFA	10 mM HCl
Recombinant human PDGF-BB protein	CYK-0025-0032 CYK-0100-0032	Lyophilized from acetonitrile/TFA	Water
Recombinant Human Interleukin-2	CYK-0025-1001 CYK-0100-1001	Lyophilized from a 0.2 µm-filtered solution containing sodium acetate and sodium chloride, pH 5.0	Acetic acid 0,1%
Recombinant Human Stem Cell Factor	CYK-0025-1002 CYK-0100-1002	Lyophilized from a 0.2 µm-filtered solution containing sodium phosphate, sodium chloride and EDTA, pH 7.4	Water
Recombinant Human Interleukin-15	CYK-0025-1003 CYK-0100-1003	Lyophilized from a 0.2 µm-filtered solution containing potassium phosphate, sodium chloride and polysorbate 20, pH 7.5	Water
Recombinant human Tumour Necrosis Factor Alpha	CYK-0025-1004 CYK-0100-1004	Lyophilized from a 0.2 µm-filtered solution containing potassium phosphate, sodium phosphate, potassium chloride and sodium chloride, pH 7.5	Water
Recombinant Human Fms-like tyrosine kinase 3 ligand	CYK-0025-1005 CYK-0025-1005	Lyophilized from a 0.2 µm-filtered solution containing potassium phosphate, sodium phosphate, potassium chloride and sodium chloride, pH 7.5	Water
Recombinant Human Interleukin-4	CYK-0025-1006 CYK-0100-1006	Lyophilized from a 0.2 µm-filtered solution containing potassium phosphate, sodium phosphate, potassium chloride and sodium chloride, pH 7.5	Water
Recombinant Human Interleukin-10	CYK-0025-1007 CYK-0100-1007	Lyophilized from a 0.2 µm-filtered solution containing potassium phosphate, sodium phosphate, potassium chloride and sodium chloride, pH 7.5	Water
Recombinant Human Interleukin-7	CYK-0025-1008 CYK-0100-1008	Lyophilized from a 0.2 µm-filtered solution containing sodium phosphate, sodium chloride and polysorbate 20, pH 7.5	Water


Product Name	SKU Number(s)	From	Recommended Reconstitution Buffer
Recombinant Human Interleukin-12	CYK-0025-1009 CYK-0100-1009	Lyophilized from a 0.2 µm-filtered solution containing sodium phosphate, mannitol, pH 6.5	Water
Recombinant Human Vitronectin	CYK-0100-1010	Lyophilized from a 0.2 µm-filtered solution containing sodium phosphate, sodium chloride, pH 7.5	Water
Recombinant Human Interleukin-3	CYK-0025-1011 CYK-0100-1011	Lyophilized from a 0.2 µm-filtered solution containing sodium phosphate and sodium chloride, pH 7.2	Water
Recombinant Human Thrombopoietin	CYK-0025-1012 CYK-0100-1012	Lyophilized from a 0.2 µm-filtered aqueous solution	Water
Recombinant Human Erythropoietin	CYK-0025-1013 CYK-0100-1013	Lyophilized from a 0.2 µm-filtered solution containing sodium phosphate, sodium chloride, mannitol and Tween-20, pH 7.0	Water
Recombinant Human Interleukin-6	CYK-0025-1014 CYK-0100-1014	Lyophilized from a 0.2 µm-filtered solution containing potassium phosphate, sodium phosphate, potassium chloride and sodium chloride, pH 7.5	Water
Recombinant Human Interleukin-18	CYK-0025-1015 CYK-0100-1015	Lyophilized from a 0.2 µm-filtered solution containing sodium phosphate, sodium chloride, mannitol and Tween-20, pH 7.2	Water
Recombinant Human Interleukin-1 Beta	CYK-0025-1016 CYK-0100-1016	Lyophilized from a 0.2 µm-filtered solution containing potassium phosphate, sodium phosphate, potassium chloride and sodium chloride, pH 7.5	Water
Recombinant Human Interferon gamma	CYK-0025-1017 CYK-0100-1017	Lyophilized from a 0.2 µm-filtered solution containing succinate, pH 5.0	Water
Recombinant Human Interleukin-21	CYK-0025-1018 CYK-0100-1018	Lyophilized from a 0.2 µm-filtered solution containing sodium acetate and sodium chloride, pH 5.5	Water
Recombinant Human Granulocyte-macrophage colony-stimulating factor	CYK-0025-1019 CYK-0100-1019	Lyophilized from a 0.2 µm-filtered solution containing potassium phosphate, sodium phosphate, potassium chloride and sodium chloride, pH 7.5	Water

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further contacts, visit
www.sartorius.com