

Reconstituting Your Vial

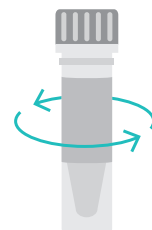
1. Receiving Delivery of Your Growth Factor | Cytokine

Once you receive delivery of your cytokine place it immediately in the freezer at -20°C or -80°C where it will remain stable for 2 years.



2. Centrifuging Your Vial

Briefly centrifuge to make sure all the powder in the vial is at the bottom of the vial before opening and be aware that the cytokine material may not always be visible.



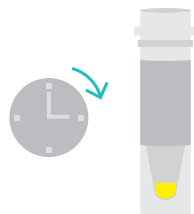
3. Adding the Buffer

Look at the table (RUO-Growth-Factors-And-Cytokines-Datasheet-en-Sartorius.pdf) and add the correct reconstitution solution to give a concentration of $0.5 - 1 \text{ mg/mL}$. For larger sample sizes, reconstitute in $\sim 75\%$ 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.



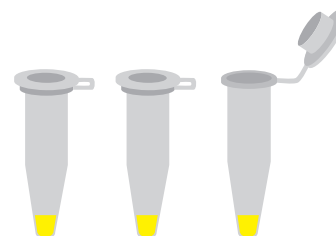
4. Allow Time to Reconstitute and Then Mix

Allow to stand for a minute at room temperature before gently agitating the mixture while being careful not to create foam – do not vortex.



5. Aliquoting for Storage

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 $< 50 \mu\text{g/mL}$, if compatible with your work to add a carrier protein (all our products are carrier protein-free).



6. Storage

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°C and -80°C .



Germany

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

For further information, visit
www.sartorius.com

Specifications subject to change without notice.
Copyright Sartorius Lab Instruments GmbH & Co. KG.
Printed in the EU or US on paper bleached without chlorine.
Version: 09 | 2023