



Customized solutions for cell cultures.

Protocol

Production of **liquid medium** from TCX10D powder

Please note, this document may be periodically updated in order to ensure the most current practices are in place. It is the user's responsibility to ensure the latest release of this protocol is applied. Valid versions are made available via Xell's webshop.

Production of liquid medium from TCX10D powder

Guideline:

We recommend preparing the whole powder container in a single batch. For that, please adjust the amounts/volumes per liter given in this protocol according to your batch size.

Material:

- TCX10D powder (23.76 g/L; Cat.No. 1150-XXXXDPM)
- H₂O (WFI or equivalent quality)
- 2.40 g/L NaHCO₃ Ph. Eur.

We recommend wearing a dust mask during preparation.



Visual control:

Check:

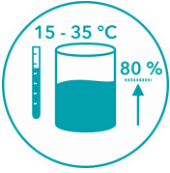


- A. Container **Sealed and without any damage.**
- B. Appearance **Free flowing powder** (record color).



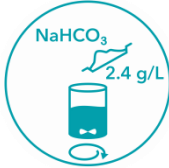






Color: _____

Procedure:

Check:

1.		Fill 0.8 L per 1 L final medium solution 15-35°C water (WFI or equivalent quality) into the stirred tank/blending vessel. Note: <i>Deviating temperature may alter dissolution rate. An adaption of time for solubilization might be necessary.</i>	<input type="checkbox"/>
2.		Start the stirrer of the system. Due to foam formation during medium production, the vortex should not reach the stirrer.	<input type="checkbox"/>
3.		Add 23.76 g/L of TCX10D powder slowly to the stirred water. Avoid clumping. Note: <i>We recommend preparing the whole powder container at once.</i>	<input type="checkbox"/>

4.		<p>Rinse the emptied powder container with a suitable amount of water (WFI or equivalent quality) and pour liquid into the stirred tank.</p>	<input type="radio"/>
5.		<p>Stir for 30 minutes with lid closed.</p> <p>Note: <i>The powder should be completely dissolved and the solution should be clear.</i></p>	<input type="radio"/>
6.		<p>Add 2.4 g/L NaHCO₃ Ph. Eur. to the stirred tank.</p>	<input type="radio"/>
7.		<p>Add an appropriate volume of water (WFI or equivalent quality) to reach the final volume.</p> <p>Note: <i>Final volume depends on batch/container size.</i></p>	<input type="radio"/>
8.		<p>Stir for 10 minutes with lid closed.</p> <p>Note: <i>The solution should be clear, without precipitates. If not, stepwise increase mixing time by 10 min.</i></p>	<input type="radio"/>
9.		<p>Measure pH (7.1 - 7.6) and osmolality (300 ± 15 mOsmol/kg) of the medium.</p>	<input type="radio"/>
10.		<p>The medium can now be sterile filtered (0.45 µm + 0.2 µm or 0.45 µm + 0.1 µm) and bottled.</p>	<input type="radio"/>

For further information or assistance
contact us.

www.xell.de
info@xell.de

Xell AG
Alte Verler Str. 1
33689 Bielefeld

Fon: +49 (0)521 96989-200
Fax: +49 (0)521 96989-201

