

Preparation of Lysates from Cultured Cells

Step 1. Prepare adherent cultured cells

1. Culture 1×10^6 to 1×10^7 cells.
2. Trypsinize the cells and remove the cells using a cell scraper.
3. Transfer cells to a 15 mL conical tube and spin at $500 \times g$ at 4°C for 5 mins.
4. Carefully aspirate the supernatant and resuspend the cells in 5 mL of ice-cold PBS and spin again at $500 \times g$ at 4°C . Repeat this step once with 5 mL of ice-cold PBS.
5. Remove as much supernatant as possible without disturbing the cell pellet. Proceed to Step 3 Lyse Cells.

Step 2. Prepare suspension cells

1. Culture 1×10^6 to 1×10^7 cells.
2. Transfer cells to a 15 mL conical tube and spin at $500 \times g$ at 4°C for 5 mins.
3. Carefully aspirate the supernatant and resuspend the cells in 5 mL of ice-cold PBS and spin again at $500 \times g$ at 4°C .
4. Remove as much supernatant as possible without disturbing the cell pellet. Proceed to Step 3 Lyse Cells.

Step 3. Lyse cells

1. To the prepared cultured cells from Step 1 or Step 2, add 200 μL of ice-cold ProcartaPlex™ Cell Lysis Buffer (EPX-99999-000) to the cell pellet.
2. Pipette up and down 8-10 times and incubate on ice for 5 mins.
3. Transfer the entire contents to a 1.5 mL microcentrifuge tube.
4. Centrifuge at $14,000 \times g$ (benchtop centrifuge) for 10 mins at 4°C .
5. Transfer the supernatant to a new tube. Assay the samples immediately or store them at -80°C .

Optional




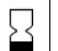


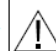
Perform a protein assay using a small aliquot of the lysate sample by following the manufacturer's protocol of a protein assay kit. We recommend Lowry Protein Assay for measuring protein concentrations of cell lysates. We recommend adjusting all samples to equal protein concentration at a range of 4-8 mg protein/mL using ice-cold ProcartaPlex™ Cell Lysis Buffer.

To proceed with the ProcartaPlex Protocol, add 25 μL of Universal Assay Buffer (EPX-11111-000) to 25 μL of the prepared sample to each sample well.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

Product label explanation of symbols and warnings

 REF	Catalog Number	 LOT	Batch code		Temperature limitation		Use by		Manufacturer		Consult instructions for use		Caution, consult accompanying documents
--	----------------	---	------------	---	------------------------	---	--------	---	--------------	---	------------------------------	---	---

Manufacturer's address: Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. s and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html.

Visit thermofisher.com/procartaplex for additional technical protocols.

Research Use Only. Not for use in diagnostic procedures.

MAN0017835 Rev. A.0 21 April 2018