

PREADYPORT™

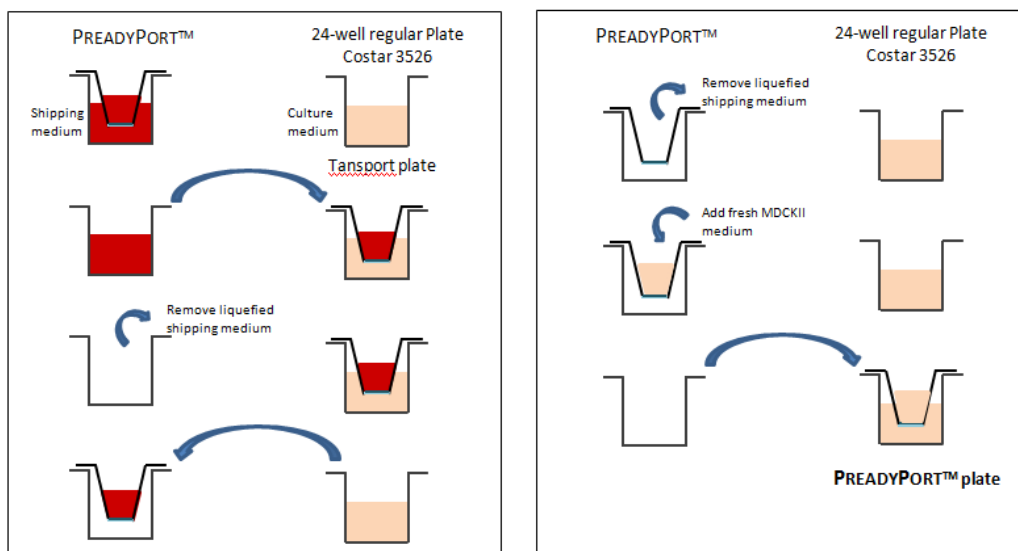
A CELL-BASED and READY-TO-USE CONCEPT FOR *IN VITRO* TRANSPORT EVALUATION

UPON RECEIPT OF PREADYPORT™ ASSAY KIT

- Remove the plastic bag containing PreadyPort™ plates from the outer and inner shipping boxes, and place the bag in a safe place protected from vibration, jarring and light.
- Store plates at 20-25°C

CHANGING THE SHIPPING MEDIUM

- To change the shipping medium, please use recommended basal plates (COSTAR ref. 3526 for 24 wells and Millipore ref PSHT004S5 for 96 wells). These plates are not provided.
- Conditions for liquefying the shipping medium of PreadyPort™ plates:
 Temperature: 37°C
 Time: 3-4 hours (**do not exceed 4 hours**).
- Pre-warm PreadyPort™ medium at 37°C before use.
- Avoid touching the cell monolayer when handling PreadyPort™ plates.
- Use a 24 or 96-well manifold (DRUMMOND ref. 3-000-097 for 24 wells and ref. 3-000-093 for 96 well) to remove cell culture medium.
- Use a multichannel pipette to dispense cell culture medium.
- Never change the shipping medium on more than **one** plate at a time.
- The Shipping medium from **both** the apical and the basal compartments must be changed as follows:



- While removing or replacing fluids, keep the basal and apical compartments separate.
- Plates should be returned to the incubator (37°C) immediately after medium has been replaced.

TEER MEASUREMENT

- Avoid contacting the cell monolayer with the electrode (WORLD PRECISION INSTRUMENTS ref. STX100C for 24 wells and ref. STX100M for 96 wells). If the recommended electrode is used, there will be no direct contact with cell monolayer.
- For optimal TEER results, cell culture medium should not be changed for 48 hours prior to measurement.
- Taking multiple TEER measurements in a short period of time can alter cell monolayer stability. We recommend leaving the plate in the incubator for a minimum of 20 minutes after carrying out TEER measurements.

For detailed instructions, please read carefully PreadyPort™ User's Manual
 For more information, please contact: reagents@advancell.net
support@solvo.com



SOLVO Biotechnology

2040-Hungary
 Gyár u. 2, Budaörs
 Fax: (36) 23 503-941
 Tel: (36) 23 503-940
 E-mail: sales@solvo.com
www.solvo.com



ADVANCELL

advanced in vitro cell technologies

PARC CIENTÍFIC DE BARCELONA
 C/ Baldri Reixac, 10 - 12
 08028 - BARCELONA - SPAIN
 TEL: (34) 93 403.45.45
 FAX: (34) 93 403.45.44
advancell@advancell.net

