

HCT-116 VIM RFP (ATCC[®] CCL-247EMT[™]) epithelial-mesenchymal transition (EMT) induction instruction

Materials:

- 1. HCT-116 VIM RFP cells (ATCC[®] CCL-247EMTTM)
- 2. McCoy's 5A Medium (ATCC[®] 30-2007[™])
- 3. Fetal Bovine Serum (FBS) (ATCC[®] 30-2020TM)
- 4. Trypsin-EDTA Solution (1X) (ATCC[®] 30-2101[™])
- 5. DPBS (ATCC[®] 30-2200TM)
- 6. Anti-miR-200 LNA (Qiagen #339160 YFI 0450012)
- 7. Opti-MEM (Thermal Fisher #31985062)
- 8. RNAiMAX (Thermal Fisher #13778030)

Instructions:

A. The following directions are for a 12 well plate. They can be scaled up or down, accordingly.

Oligo stock should be re-suspended at a final concentration of 50 μ M.

- B. Culture cells for induction: Thaw one vial of HCT-116 VIM RFP cells and plate the cells in T-75 flask. Cell cultures generally reach 70-80% confluent in about 2-3 days.
- C. Induction protocol:
- 1. Prepare HCT-116 VIM RFP (ATCC[®] CCL-247EMTTM) cells for induction
 - 1.1 Warm culture medium (McCoy's 5A with 10% FBS) to 37 °C.
 - 1.2 Harvest HCT-116 VIM RFP cells using Trypsin-EDTA solution.
 - 1.3 Resuspend the cells in warmed culture medium. Centrifuge the cell suspension @ approximately 250 X g for 5 minutes. Aspirate the liquid.
 - 1.4 Gently resuspend the cell pellet in warmed culture media and count viable cells using Trypan blue.
 - 1.5 Reuspend cells at a density of 2.0 x 10⁵ cells/mL and put the cells in a 37°C + 5% CO₂ incubator.
- 2. Prepare siRNA transfection complexes:
 - 2.1 Bring RNAiMAX and Opti-MEM to room temperature.
 - 2.2 Label two sterile microcentrifuge tubes (–CONTROL, and +LNA) and pipette 100 μl Opti- MEM to each of them.
 - 2.3 Add 3 μ L of LNA to '+LNA' tube and 3 μ L of DPBS to '-CONTROL' tube. Gently mix.
 - 2.4 Add 3 µL of RNAiMAX to each tube. Mix thoroughly.
 - 2.5 Incubate transfection complexes at room temperature for 10 minutes.
- 3. Mix transfection complexes with cells
 - 3.1 Transfer the prepared '-CONTROL' and '+LNA' transfection complexes to wells of a 12-well plate.
 - 3.2 Mix cell suspension and add 500 μ l of the cell suspension to the appropriate wells.

American Type Culture Collection

P.O. Box 1549 Manassas, VA 20108 USA http://www.atcc.org

© 2018 ATCC. All rights reserved.



THE ESSENTIALS OF LIFE SCIENCE RESEARCH GLOBALLY DELIVERED **

- 3.3 Add an additional 500 μL of culture media to each well.
- 3.4 Gently rock plate back and forth at least 10 times.
- 3.5 Put the cells into a 37°C with 5% CO₂ incubator to complete the 1st induction.
- 4. When the cells grow to approximately 70-80% confluent (usually 3-4 days), repeat step 1 to 3 to complete the 2nd induction (Note: The number of live cells and the cell growth rate may decline due to toxicity of LNA to the cells).
- 5. When the cells grow to approximately 60-80% confluent (usually ~4 days), perform the 3rd induction. The 3rd induction uses half dose of LNA (100 μ L Opti-MEM + 1.5 μ L of LNA or DPBS + 1.5 μ L of RNAiMAX).
- 6. When the cells reach approximately 60-80% confluent, repeat step 5 to perform 4th induction cycle, and so on. Generally, RFP expression can be observed after the 6th induction.

Reference for preparing siRNA transfection complexes:

	-CONTROL	+LNA
Opti-MEM media	100 μL	100 μL
Oligos		3 μL or 1.5 μL
DPBS	3 μL or 1.5 μL	
RNAiMAX	3 μL or 1.5 μL	3 μL or 1.5 μL
Cell Suspension	500 μL	500 μL
Culture Media	500 μL	500 μL

Reference:

Park SM, *et al*. The miR-200 family determines the epithelial phenotype of cancer cells by targeting the E-cadherin repressors ZEB1 and ZEB2. *Genes Dev* 22: 894-907, 2008. PubMed: 18381893