



TAG-LITE® BETA2 STABLE CELL LINE

Cell line information - Version: rev 3 (Jul. 2020)

Part#: C1SU1BETA2

Storage temperature: Liquid nitrogen

Lot: 001

Passages number: 13

Total cells: 5 millions/vial

For in vitro research use only. Not for use in diagnostic procedures.

TARGET INFORMATION

Target: Adrenoceptor beta2 receptor. This receptor is fused in N-term position with SNAP-tag.

Species: Human

NCBI accession number: NM_000024.4

CELL CULTURE INFORMATION

Cellular background: HEK293 Adherent

Culture medium: D-MEM (Invitrogen #31966) ,10%FBS, 1% non essentials amino acids (Invitrogen #11140-035), 50U/ml penicillin, 50µg/ml streptomycin, 2mM Hepes, 0.6mg/mL G418

Growth conditions: 37°C, 5%CO₂

Freezing medium: 90% FCS+10%DMSO

Thawing procedure: Thaw the vials of cryopreserved cells rapidly in a water bath at 37°C. Carefully transfer the cells to a cell culture vessel containing the appropriate volume of pre-warmed culture medium without G418 so that cell density reaches 4.104 to 8.104 cells/cm². The day after, ensure that cells have properly attached and carefully replace with fresh culture medium supplemented with G418. Allow cells to grow until 80% confluency then transfer in a new cell culture vessel. Note that final DMSO concentration must not exceed 0.5% after the initial transfer. Use of T75 is recommended initially.

Cell adhesion may not happen overnight. If this is the case, waiting for 2-3 days may be necessary before being able to proceed with replacement by medium + G418 (you should wait until you have good cell adhesion before replacement for medium + G418). If the cells don't attach after 2-3 days, we recommend collecting the cells, centrifuging, resuspending in medium without G418, and seeding in a T25 flask. Supplementing with G418 will be possible once cells have attached to the flask. Then when 80% confluence is reached, it is possible to expand to larger flasks. G418 supplementation which is too early may considerably slow down the cells division.

Subculturing: Once cells have reached 80% confluency, carefully remove and discard culture medium. Rinse once with PBS then add the appropriate volume of cell dissociation solution. Observe under microscope until cell layer is detached. Add growth medium, then centrifuge 3min at 180g. Replace supernatant by fresh cell culture medium and pipet up and down to remove aggregates. Count cells, and split in a new culture vessel with a ratio 1:5 maximum. Replace medium once or twice a week.

CELL CULTURE INFORMATION

Biosafety level: 2

Warning: This product contains material of biological origin. Use for research purposes only. Do not use in humans or for diagnostic purposes. The purchaser assumes all risk and responsibility concerning reception, handling and storage. The use of the cell line will be done with appropriate safety and handling precautions to minimize health and environmental impact. The product is genetically modified and must be used according to biosafety level S2. The purchaser declares the authorization to manipulate GMO and agrees to apply all guidelines, laws and regulations.

TAG-LITE STABLE CELL LINE VALIDATION

Tag-lite cell line is a comprehensive cellular platform certified for:

- Ligand binding assay

FOR MORE INFORMATION

Europe and other countries +33(0)466-796-705 U.S. and Canada 1-888-963-4567 China +86-21-5018-9880

Japan +81-(0)43-306-8712 Visit www.cisbio.com to find a list of our regional distributors



cisbio