



# TAG-LITE® MC4 STABLE CELL LINE

Cell line information - Version: rev 0 (Sept. 2011)

**Part#:** C1SU1MC4

**Storage temperature:** Liquid nitrogen

**Lot:** 001

**Passages number:** 8

**Total cells:** 4 millions/vial

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

## TARGET INFORMATION

**Target:** Melanocortin MC4 receptor. This receptor is fused in N-term position with SNAP-tag.

**Species:** Human

**NCBI accession number:** NM\_005912.2

## 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%CO<sub>2</sub>

**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 4.10<sup>4</sup> to 8.10<sup>4</sup> cells/cm<sup>2</sup>. 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%.

**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
- cAMP 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](http://www.cisbio.com) to find a list of our regional distributors

Copyright 2015 Cisbio Bioassays. All rights reserved. HTRF, Tag-lite, EPIgeneous and the HTRF logo are trademarks or registered trademarks of Cisbio Bioassays. HaloTag is a registered trademark of Promega Corporation.

