

## Tag-Lite® CCK1 stable cell line

[www.htrf.com](http://www.htrf.com)

### Cell line Information

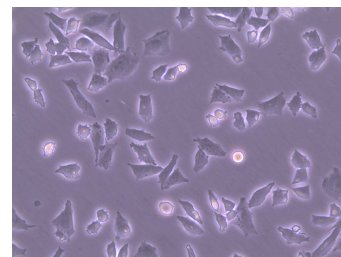
For in vitro research use only

Version Feb1A

<b>Cell line code</b>	C2SU1CCK1
<b>Storage</b>	Liquid nitrogen
<b>Batch</b>	21 dec 2009
<b>Passages number</b>	5
<b>Total cells</b>	9 millions/vial

### Target information

<b>Target:</b>	Cholecystokinin CCK1 receptor This receptor is fused in N-term position with SNAP-tag.
<b>Species:</b>	Human
<b>NCBI accession number:</b>	NM_000730.2



### Cell culture information

<b>Cellular background:</b>	CHO Adherent
<b>Culture medium:</b>	Ham F12+10%FCS+1mM Na-Pyruvate+2mM l-Glutamin+1mg/ml G418
<b>Growth conditions:</b>	37°C, 5%CO2
<b>Freezing medium:</b>	90% FCS+10%DMSO

#### Thawing procedure:

Thaw the vials of cryopreserved cells rapidly in a water bath at 37°C. Carefully transfer the cells in a cell culture vessel containing the appropriate volume of pre-warmed culture medium without G418 so that cell density reaches  $4 \cdot 10^4$  to  $8 \cdot 10^4$  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 for growth until 80% confluency then transfer in a new cell culture vessel.

#### 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.

### Biosafety and compliance

**Biosafety level:** 1

#### Warning:

This product contains material of biologic origin. Use for research purposes only. Do not use in humans or for diagnostic purposes. The purchaser assumes risk and responsibility concerning reception, handling and storage. The use of the cell line will be done in appropriate safety and handling precautions to minimize health and environmental impact. The product is genetically modified and must be used according to biosafety level S1. The purchaser declares the authorization to manipulate GMO and agrees to apply guidelines, laws and regulations.

### Tag-lite stable cell line validation

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

Ligand binding assay  
Receptor dimerization assays  
IP-One assay  
Phospho-ERK assay  
Internalization assay

For each application a protocol can be downloaded on [www.htrf.com](http://www.htrf.com)