

Tag-Lite® CXCR7 stable cell line

www.htrf.com

Cell line Information

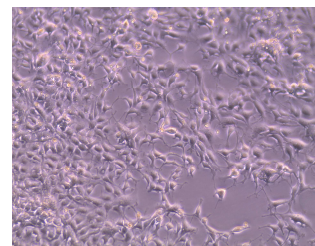
For in vitro research use only

Version Feb1A

Cell line code	C1SU1CXCR7
Storage	Liquid nitrogen
Batch	11 dec 2009
Passages number	12
Total cells	11 millions/vial

Target information

Target:	Chemokine CXCR7 receptor This receptor is fused in N-term position with SNAP-tag.
Species:	Human
NCBI accession number:	NM_020311.2



Cell culture information

Cellular background:	HEK293 Adherent
Culture medium:	DMEM (4.5g/l Glucose)10%FCS+4mM L-Glutamine+0.6µg/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 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/cm². 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:

Receptor dimerization assays

Phospho-ERK assay

Internalization assay

For each application a protocol can be downloaded on www.htrf.com