

Datasheet for H1299/RFP Labeled A-Tubulin Cell Line

Catalog number: SL202

Product: H1299 cell line stably expressing RFP Labeled A-Tubulin.

Description: This product is a cell line stably expressing the A-Tubulin. A-Tubulins are

Microtubules of the eukaryotic cytoskeleton perform essential and diverse functions and are composed of a heterodimer of alpha and beta tubulins.. This

cell line also expresses RFP protein and the puromycin resistance gene.

Quantity: 1 vial of 2 x 10⁶ cells; frozen

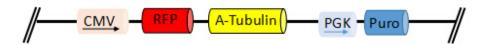
Shipping conditions: Dry ice

Storage conditions: Liquid nitrogen vapor phase. Remove the item from the dry ice packaging and

check all items for damage and leakage. Place immediately into storage at or

below -140 °C, preferably into the liquid nitrogen vapor phase, until use.

Transgene integration:



CMV-RFP-A-Tubulin-PGK-Puro

Source of parental line:

H1299

Organism: Homo sapiens, human

Tissue: lung

Cell type: Epithelial

Quality control: >95% viability before freezing. All cells were tested and found to be free of

mycoplasma, bacteria, viruses, and other toxins.

Safety instructions: To ensure safety, protective gloves, clothing, and a face mask should be worn

when handling frozen vials. Some leakage may occur into the vial during storage. The liquid nitrogen will be converted to gas upon thawing. Due to the nature of nitrogen gas, pressure may build within the vial and possibly result in the vial

exploding or losing its cap. This may cause flying debris.



Thawing procedure: The vial of cells should be thawed in a 37 °C water bath with gentle agitation. For optimal performance, the vial should be thawed in under two minutes. Ensure that the cap of the vial did not loosen upon thawing, and re-tighten as needed. Spray the vial with 70% EtOH and wipe off. Repeat. Using aseptic technique, add the contents of the vial to 9 ml of complete growth medium (without selection). Centrifuge for 5 min. at 125 x g. Aspirate the medium, being careful not to disturb the pellet. Resuspend in 10 mL of complete growth medium, and place into a culture vessel of your choice. Only add selection to the medium after 24 hours in culture.

Culture conditions:

Complete Growth Medium

The base medium for this cell line is DMEM. For optimal growth and maintenance of selection, add the following components to the base medium: dialyzed fetal bovine serum to a final concentration of 10%.

Selection

Puromycin to a final concentration of 2 µg/mL

Culture temperature:

37 °C with 5% CO₂

Subculture:

Replace culture medium with selection-free medium and incubate for up to 6 hours. Rinse the cells with PBS without cations, digest cells with 0.25% (w/v) Trypsin-EDTA (0.53 mM) solution and split at 1:3 to 1:10 ratio.



Cryopreservation: Freeze slowly in complete growth medium supplemented with 5% (v/v) DMSO.

Mycoplasma:

Negative

(MycoGuardTM Mycoplasma Bioluminescent Detection Kit from GeneCopoeia)

Product QC:

1. Transduction

(1) Plate cells

Typsinize and count the H1299 cells before one day. Adjust the cell concentration and seed H1299 cells in a 6-well plate ,The number of cells plated in each well should be determined so that are 50-60% confluent at the time of transduction.

(2) Add Lentiviral particles

Add LP506 Lentiviral particles in H1299 cells according to the MOI=1 (LP506, Lentiviral particles of RFP-A-Tubulin, GeneCopoeia: 9.25x10⁷ TU/ml).

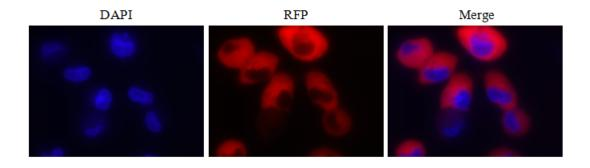
2. Drug selection

Typsinize the H1299 cells into 10cm^2 dish post-transduction 72 hours, add the puromycin in dish with final concentration 2 μ g/mL.observate the drug selection after 3 days,Change the final concentration of puromycin (1 μ g/mL) when the un-transduction cells were died.

3. Cells climbing to the carry sheet glass

Fluorescence:

Image of H1299/RFP Labeled A-Tubulin Cell Line. Observing the cells climbing to the carry sheet glass, which are stained with DAPI.



Citation of product: If use of this item results in a publication, please use this information: H1299/RFP Labeled A-Tubulin Cell Line (SL201, GeneCopoeia, Inc., Rockville, MD).



Limited Use License

A limited use license is granted to the Buyer of the Product. The Product shall be used by the Buyer for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products without prior written consent from GeneCopoeia. This Product should be used in accordance with NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

Copyright ©2016 GeneCopoeia, Inc.

C9SCL-DS-031116

GeneCopoeia, Inc. 9620 Medical Center Drive, #101 Rockville, MD 20850 USA

Tel: 301-762-0888; Fax: 301-762-3888 Email: support@genecopoeia.com Web: www.genecopoeia.com