

Datasheet for HEK293/dCas9-VP64+MPH Cell Line

Catalog number: SL302

Product: HEK293 cell line stably expressing dCas9-VP64 and MS2-P65-HSF1

Description: This product is a cell line stably expressing the catalytically inactive, HA-tagged

dCas9 nuclease fused to the VP64 transcriptional activation domain and the activator helper complex MS2-P65-HSF1 (MPH). This cell line also contains blasticidin and hygromycin resistance genes. The target guide RNAs (with MS2-binding loops) can be transfected or transduced in the cell line to activate the

gene transcription.

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.

Source of parental line:

HEK293

Organism: Homo sapiens, human

Tissue: embryonic kidney Cell type: epithelial

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

when handling frozen vials. Some leakage of liquid nitrogen into the vial may occur 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. Spray the

vial with 70% EtOH and wipe off. Repeat. Using aseptic technique, add the contents of the vial to 9ml of complete growth medium (without selection). Centrifuge for 5min at 125xg. Aspirate the medium, being careful not to disturb the pellet. Re-suspend in 10mL 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: fetal bovine serum to a final concentration of 10%.

Selection

Blasticidin to a final concentration of 20 $\mu g/mL$, and hygromycin to a final concentration of 250 $\mu g/mL$.

Culture temperature:

37°C with 5% CO₂

Subculture:

Rinse the cells with PBS without cations, digest cells with 0.25% (w/v) Trypsin-EDTA (0.53mM) solution and split at 1:6 to 1:10 ratio.

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

Quality control: The genome integration, expression levels of dCas9-VP64 and MPH, and the

transcription activation of target genes using the cell line have been verified by qPCR. All other QC specifications have also been met. For detailed information,

please refer to the certificate of analysis.



Citation of product: If use of this item results in a publication, please use this information: CRISPR stable cell line HEK293/dCas9-VP64+MPH (SL302, 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 ©2018 GeneCopoeia, Inc.

C9SCL-DS-060618

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