

MULTISCREEN™ HOMOZYGOUS STABLE CELL LINE HUMAN RECOMBINANT GPR65 RECEPTOR CRISPR KNOCK OUT

Data sheet

PRODUCT INFORMATION

Catalog Number: C1121-CSP75

Lot Number: C1121-CSP75-051523

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Cell Banker 2
(Amsbio 11891)

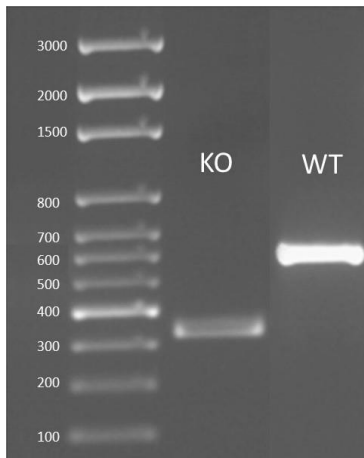
Host cell: Mia-Pa-Ca-2

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Stability: Stable for a minimum of 2 months in continuous culture

Figure 3: PCR confirmation of GPR65 knock-out clone by agarose gel electrophoresis



Background: GPR65 also known as T cell death-associated gene 8 (TDAG8) is a proton-sensing GPCR and plays a major role in pH homeostasis. This receptor is primarily expressed in lymphoid tissues (spleen, thymus, leukocytes and lymph nodes) and cancer tissues. The major function of this receptor is to reduce immune-mediated inflammation by regulating cytokine production from T cells and macrophages.

Application: Functional assays

Figure 1: Verified GPR65 coding sequence with deletion highlighted in grey.

```
ATGAACAGCACATGTATTGAAGAACAGCATGACCTGGATCACTATTTGTTTCCCATTGT
TTACATCTTTGTGATTATAGTCAGCATTCCAGCCAATATTGGATCTCTGTGTGTCTCT
TCCTGCAAGCAAAGAAGGAAAGTGAAGTGAAGTGAAGTGAAGTGAAGTGAAGTGAAGT
GATTTACTCTATGCATTAAGTCTCCCTTTATGGATTGATTATACCTGGAATAAAGACAA
CTGGACTTTCTCTCCTGCCTTGTGCAAAAGGGAGTGCTTTTCTCATGTACATGAATTTT
ACAGCAGCACAGCATTCCTCACCTGCATTGCCGTTGATCGGTATTTGGCTGTTGTCTAC
CCTTTGAAGTTTTTTTTTCTAAGGACAAAGAAGATTGCACTCATGGTCAGCCTGTCCAT
CTGGATATTGGAACCATCTTCAATGCTGTCTGTGTGGGAAGATGAAACAGTTGTG
AATATTGCGATGCCGAAAAGTCTAATTTTACTTTATGCTATGACAAATACCCTTTAGAG
AAATGGCAAATCAACCTCAACTTGTTCAGGACGTGTACAGGCTATGCAATACCCTTTGGT
CACCATCCTGATCTGCAACCGAAAGTCTACCAAGCTGTGCGGCACAATAAAGCCACGG
AAAACAAGGAAAAGAAGAGAATCATAAACTACTTGTGATCATCACAGTTACTTTTGTCT
TTATGCTTTACTCCCTTTCATGTGATGTTGCTGATTCGCTGCATTTTAGAGCATGCTGT
GAACCTCGAAGACCACAGCAATTCGGGAAGCGAACTTACACAATGTATAGAATCACGG
TTGCATTAACAAGTTAAATTTGTGTTGCTGATCCAATTCTGTACTGTTTTGTAAACCGAA
ACAGGAAGATATGATATGTGGAATATATTAATTTCTGCACTGGGAGGTGTAATACATC
ACAAAGACAAAGAAAACGCATACCTTTCTGTGTCTACAAAAGATACTATGGAATTAGAGG
TCCTTGAGTAG
```

Figure 2: Verified GPR65 knock out protein sequence product

```
MNSTCIEEQHDLHDHYLFPVIYIVFVIIVSIPANIGSLCVSFLQAKKESELGIYFLSLSL
DLLYALTLPLWIDYTNKDNWTFSPALCKGSAFLMYMNFYSSTAF LTCIQAMQYLWSPS
*
```

References:

Ludwig, M.-G., Vanek, M., Guerini, D., Gasser, J. A., Jones, C. E., Junker, U., Hofstetter, H., Wolf, R. M., Seuwen, K. Proton-sensing G-protein-coupled receptors. *Nature* 425: 93-98, 2003.

Saxena, H., Deshpande, D., Tiegs, B., Yan, H., Battafarano, R., Burrows, W., Penn, R. (2012). The GPCR OGR1 (GPR68) mediates diverse signalling and contraction of airway smooth muscle in response to small reductions in extracellular pH. *British Journal of Pharmacology*, 166(3), 981-990.

Satoshi Ishii, Yasuyuki Kihara and Takao Shimizu (2005) Identification of T Cell Death-associated Gene 8 (TDAG8) as a Novel Acid Sensing G-protein-coupled Receptor. *J Biol Chem* 280: 9083-9087

FOR RESEARCH USE ONLY.

All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.