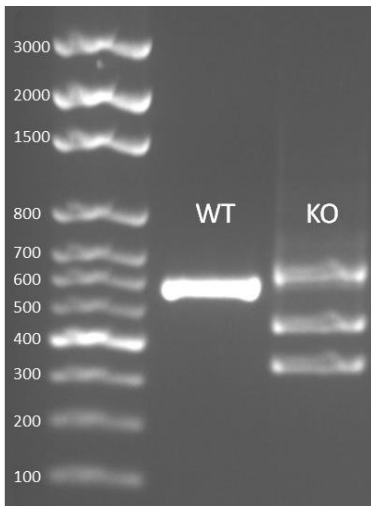


MULTISCREEN™ HETEROZYGOUS STABLE CELL LINE HUMAN RECOMBINANT GPR65 RECEPTOR CRISPR KNOCK OUT

PRODUCT INFORMATION

Catalog Number: C1121-CSP20
Lot Number: C1121-CSP20-051523
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



Data sheet

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.

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

```
ATGAACAGCACATGTATTGAAGAACAGCATGACCTGGATCACTATTGTTTCCCATTGT
TTACATCTTTGTGATATAGTCAGCATTCCAGCCAATATTGGATCTCTGTGTGTCTT
TCCTGCAAGCAAAGAAGGAAAGTGAAGTGAAGTGAAGTGAAGTGAAGTGAAGTGAAGT
GATTACTCTATGCATTAAGTCTCCCTTTATGGATTGATTATACCTGGAATAAAGACAA
CTGGACTTCTCTCCTGCCTTGTGCAAAGGGAGTGCCTTTCTCATGTACATGAATTTT
ACAGCAGCACAGCATTCCCTCACCTGCATTGCCGTTGATCGGTATTTGGCTGTTGTCTAC
CCTTTGAAGTTTTTTTTTCTAAGGACAAGAAGATTTGCACTCATGGTCAGCCTGTCCAT
CTGGATATTGGAACCATCTTCAATGCTGTCATGTTGGGAAGATGAAACAGTTGTTG
AATATATGCGATGCCGAAAAGTCTAATTTTACTTTATGCTATGACAAATACCCCTTAGAG
AAATGGCAAATCAACCTCAACTGTTTCAGGACGTGTACAGGCTATGCAATACCTTTGGT
CACCATCCTGATCTGCAACCGGAAAGTCTACCAAGCTGTGCCGCAATAAAGCCACGG
AAAACAAGGAAAAGAAGAAATCATAAACTACTTGTGTCAGCATCACAGTTACTTTTGTCT
TTATGCTTTACTCCCTTTTTCATGTGATGTTGCTGATTTCGCTGCATTTTAGAGCATGCTGT
GAACCTCGAAGACCACAGCAATTTGGGAAGCGAAGTACACAATGTATAGAATCACGG
TTGCATTAACAAGTTTAAATTTGTTGTTGCTGATCCAATTCTGTACTGTTTTGTAACCGAA
ACAGGAAGATATGATATGTGGAATATATTTAAATTTCTGCACTGGGAGGTGTAATACATC
ACAAAGACAAAGAAAACGCATACTTTCTGTGTCTACAAAAGATACTATGGAATTAGAGG
TCCTTGAGTAG
```

Figure 2: Verified GPR65 knock out protein sequence product

```
MNSTCIEEQHDLHDHYLFPIVYIFVIIVSIPANIGSLCVSFLQAKKESELGIYLFSLSL
DLLYALTLPLWIDYTWNKDNWTFSPALCKGSAFLMYMNFYSSTAFITCIQAMQYLLWSP
S*
```

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.