

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT GPR97 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1396

**Lot Number:** C1396

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human GPR97 cDNA (GenBank accession Number BC064508) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

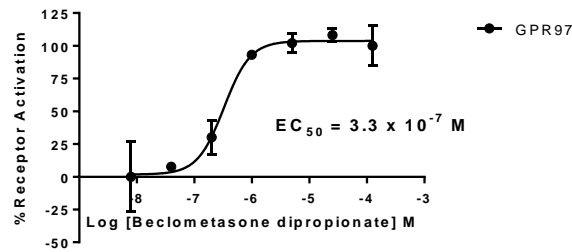
**Propagation Medium:** DME/F12, 10% FBS, 10  $\mu$ g/mL puromycin

**Stability:** In progress

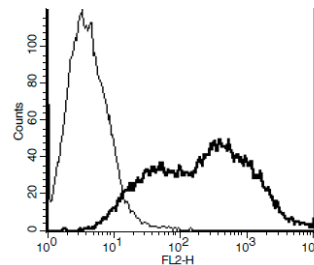
**Background:** Human G-protein coupled receptor 97 (ADGRG3), a member of the adhesion GPCR family, is a recently orphanized receptor for beclomethasone dipropionate. GPR97 is expressed in immune cells, specifically thymocytes and pre-B cells. Current research studies on knockout mice demonstrate the importance of GPR97 in B-cell development and its potential as a therapeutic target for the treatment of immunological disorders.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC).

**Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

#### References:

Fredriksson R, Lagerström MC, Höglund PJ, Schiöth HB (Nov 2002). "Novel human G protein-coupled receptors with long N-terminals containing GPS domains and Ser/Thr-rich regions". *FEBS Letters* **531** (3): 407–14.

Stacey M, Yona S (2011). *Adhesion-GPCRs: Structure to Function (Advances in Experimental Medicine and Biology)*. Berlin: Springer.

Wang, J., Zhang, L., Zhang, H., Shen, C., Lu, S., Kuang, Y., ... Wang, Z. (2013). Gpr97 is essential for the follicular versus marginal zone B-lymphocyte fate decision. *Cell Death & Disease*, **4**(10), e853–. doi:10.1038/cddis.2013.346.

**FOR RESEARCH USE ONLY.**

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com  
sales@multispaninc.com  
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817  
Fax: +1 (510) 887-0863  
26219 Eden Landing Road  
Hayward, CA 94545-3718  
U.S.A.