

**MULTISCREEN™ MEMBRANE PREPARATION  
HUMAN RECOMBINANT GPR17 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** MC1076-3a

**Lot Number:** MC1076-3a-091115

**Quantity:** 19.2 mg/ml, 5 mg

**Packaging Buffer:** 20 mM Gly-Gly, 1 mM MgCl<sub>2</sub>, 25 mM Sucrose (pH 7.2)

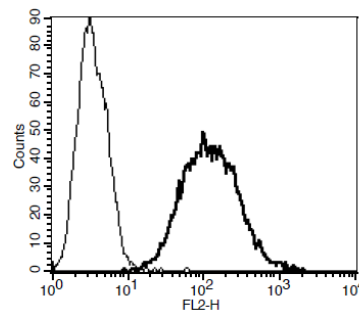
**Host cell:** 1321N1

**Transfection:** Expression vector containing full-length human GPR17cDNA (GenBank Accession Number NM\_005291) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

**Background:** G-protein coupled receptor 17 (GPR17) is closely related to the purinergic and cysteinyl-leukotriene receptor families. It is present on both neurons and a subset of oligodendrocyte precursor cells. Recent studies have shown that GPR17 plays an important role in both sensing and repairing brain damage. Thus, GPR17 represents a potential new target for the treatment of traumatic brain injuries as well as neurodegenerative diseases like Alzheimer's and multiple sclerosis.

**Figure 1**



**Figure 1.** 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:**

- Franke, Heike et al. "Changes of the GPR17 Receptor, a New Target for Neurorepair, in Neurons and Glial Cells in Patients with Traumatic Brain Injury." *Purinergic Signalling* 9.3 (2013): 451–462. *PMC*. Web. 18 Feb. 2015.
- Hennen, S et al. "Decoding Signaling and Function of the Orphan G Protein-Coupled Receptor GPR17 with a Small-Molecule Agonist." *Sci Signal* 2013 Oct 22;6(298)
- Lecca, Davide et al. "The Recently Identified P2Y-Like Receptor GPR17 Is a Sensor of Brain Damage and a New Target for Brain Repair." Ed. Kenji Hashimoto. *PLoS ONE* 3.10 (2008): e3579. *PMC*. Web. 18 Feb. 2015.

**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.