

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT PK2 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1127-1a

**Lot Number:** C1127-1a-070711

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

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

**Host cell:** CHO-K1

**Transfection:** Expression vector containing full-length human PK2 cDNA (GenBank Accession Number NM\_144773.2) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

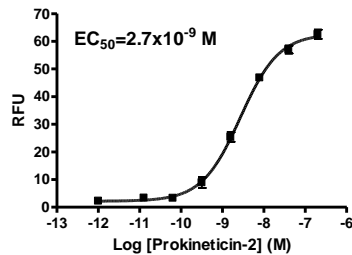
**Propagation Medium:** DMEM-F12, 10% FBS, 10 µg/ml puromycin

**Stability:** In progress

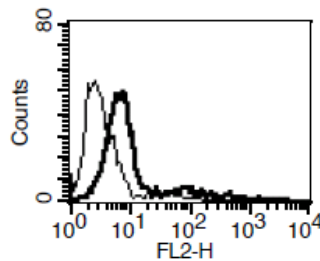
**Background:** Prokineticin receptor 2 (PK2 or GPR73L1) is a receptor for the cysteine-rich secreted peptides prokineticin 1 and 2. RT-PCR detected PK2 expression in the brain, testis, small intestine, ovary, thyroid, pituitary and salivary gland.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **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:

Lin *et al.* (2002) Identification and molecular characterization of two closely related G protein-coupled receptors activated by prokineticins/endocrine gland vascular endothelial growth factor. *J Biol Chem* 277:19276-19280.

**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.