

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT PK2 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1127-1

Lot Number: C1127-1-070711

Quantity: 1 vial (2 x 10⁶) 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: Stable in culture for minimum of two months

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

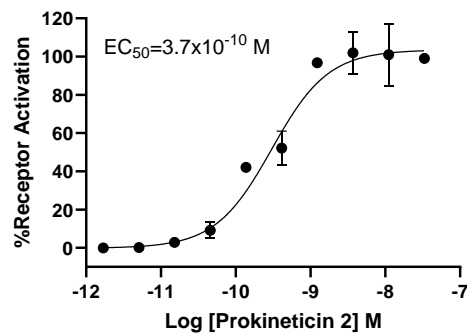


Figure 2

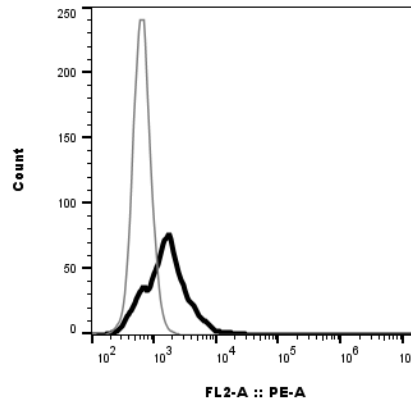


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Black line: parental cells; gray 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.

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