

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT PK2 RECEPTOR, MUTANT

Data sheet

PRODUCT INFORMATION

Catalog Number: C1127CT-1

Lot Number: C1127CT-1-020409

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Expression vector containing human PROKR2 cDNA (GenBank Accession Number: NM_144773.2) with FLAG tag sequence at N-terminus and 15 bp truncated at the C-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM-F12,
10% FBS, 10 μ g/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.

The Multispan PK2 mutant cell line stably expresses the human PROKR2 with the last five amino acids truncated at the C-terminus.

Application: Functional assays

Figure 1

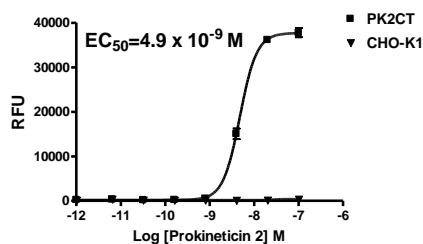


Figure 2

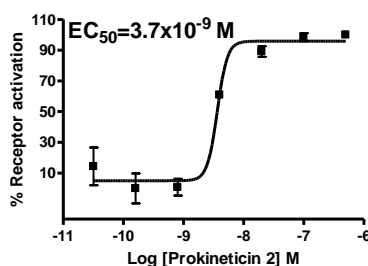


Figure 3

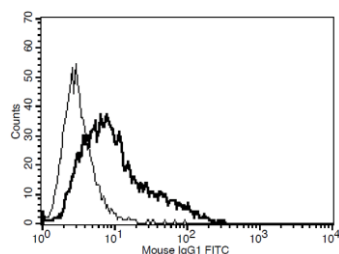


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Dose response of intracellular IP1 accumulation upon treatment with ligand, measured with IP-one Tb kit. **Figure 3.** Receptor expression 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.

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