

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT PK1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1128-1

Lot Number: C1128-1-031209

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Expression vector containing full-length human PK1 cDNA (GenBank Accession Number: NM_138964) with FLAG tag sequence at N-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 1 (PK1 or GPR73) is a receptor for the cysteine-rich secreted peptides prokineticin 1 and 2. RT-PCR detected GPR73 expression in the brain, testis, small intestine, ovary, thyroid, pituitary and salivary gland.

Application: Functional assays

Figure 1

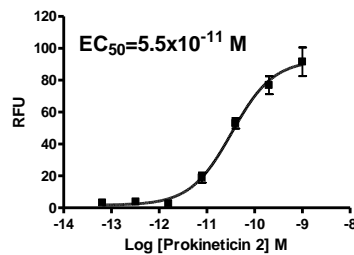


Figure 2

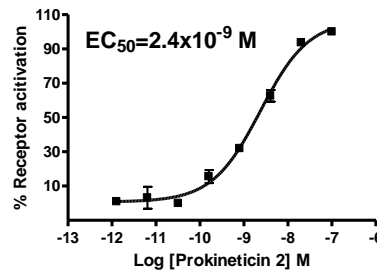


Figure 1. Dose-dependent stimulation of calcium flux measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Dose response of intracellular IP1 accumulation upon treatment with ligand, measured with IP-one Tb kit (Cisbio 62IPAPPEC).

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.