

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT PRRP RECEPTOR

PRODUCT INFORMATION

Catalog Number: C1037

Lot Number: C1037-080510

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO dhfr⁻

Transfection: Full-length human GPR10 cDNA (GenBank Accession Number BC095539)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: Alpha-MEM, 10% FBS, 500 µg/mL G418

Stability: Stable in culture for minimum of two months

Data sheet

Background: Prolactin-releasing peptide receptor PRRP (or GPR10) is detected in human brain, pituitaries, normal portions of adrenal glands and various tumor tissues. High level of expression is detected in pheochromocytomas. GPR10 was investigated regarding the effect of normal and high-fat diets on energy intake, body weight, and glucose homeostasis in wild-type and GPR10 knockout mice. The investigation provided direct evidence that GPR10 is involved in the regulation of energy balance.

Application: Functional assays

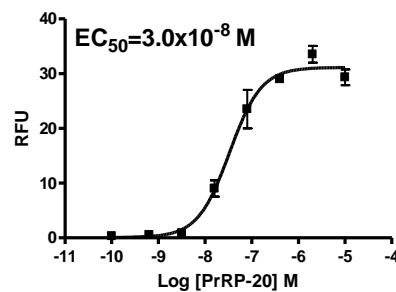


Figure legend: Dose-dependent calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Gu *et al.* (2004) The prolactin-releasing peptide receptor (GPR10) regulates body weight homeostasis in mice. *J Mol Neurosci* 22:93-103.

Takahashi *et al.* (2002) Expression of prolactin-releasing peptide and its receptor in the human adrenal glands and tumor tissues of adrenocortical tumors, pheochromocytomas and neuroblastomas. *Peptides* 23:1135-1140.

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.