

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT NPBW2 RECEPTOR**

PRODUCT INFORMATION

Catalog Number: C1133

Lot Number: C1133-061505

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T Gaqi5

Transfection: Full-length human NPBWR2 cDNA (GenBank Accession Number NM_005286)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 5 $\mu\text{g}/\text{mL}$ blasticidine, 0.625 $\mu\text{g}/\text{mL}$ puromycin

Stability: Stable after minimum of two months continuous growth

Data sheet

Background: Neuropeptide B/neuropeptide W receptor 2 (NPBW2 or GPR8) is one of the two receptors for neuropeptides B and W, which may be involved in neuroendocrine system regulation, food intake and the organization of other signals.

Application: Functional assays

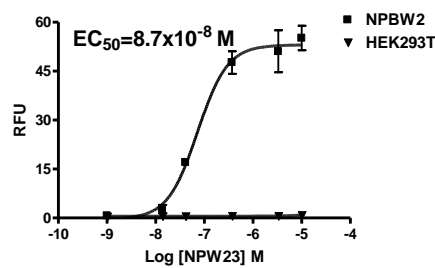


Figure legend: Dose-dependent calcium flux upon treatment with ligand, monitored with FlexStation.

References:

Tanaka *et al.* (2003) Characterization of a family of endogenous neuropeptide ligands for the G protein-coupled receptors GPR7 and GPR8. *Proc Natl Acad Sci USA* 100:6251-6256.

Levine *et al.* (2005) Injection of neuropeptide W into paraventricular nucleus of hypothalamus increases food intake. *Am J Physiol Regul Integr Comp Physiol* 288:R1727-R1732.

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.