

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

**PRODUCT INFORMATION**

**Catalog Number:** C1133

**Lot Number:** C1133-061505

**Quantity:** 1 vial ( $2 \times 10^6$ ) 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

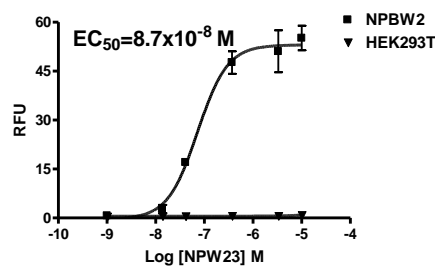
**Propagation Medium:** DMEM, 10% FBS, 5  $\mu\text{g}/\text{mL}$  blasticidine, 0.625  $\mu\text{g}/\text{mL}$  puromycin

**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, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

**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.**

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