

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE
HUMAN RECOMBINANT BB1 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1211

Lot Number: DC1211-122614

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human NMBR cDNA (GenBank Accession Number NM_002511.1) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Stability: 1-2 days after thaw

Background: The human BB1 receptor (or Neuromedin B receptor NMBR) is a receptor for neuromedin-B (NMB), which is a mammalian bombesin-like peptide distributed widely in the central nervous system. The BB1 pathway is involved in the regulation of a wide variety of behaviors, such as spontaneous activity, feeding and anxiety-related behavior. A study using BB1-deficient mice suggested that dysfunction in the BB1 pathway may constitute one of the risk factors of stress vulnerability.

Application: Functional assays

Figure 1

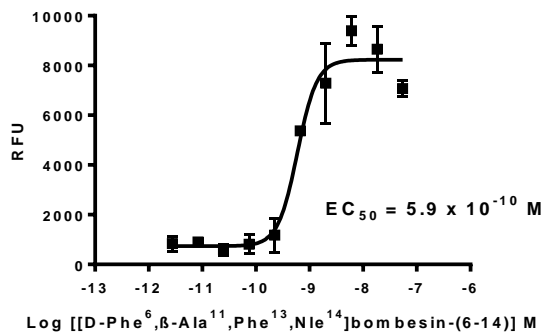


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Benya *et al.* (1995) Expression and characterization of cloned human bombesin receptors. *Mol Pharmacol* 47:10-20.

Moody *et al.* (2000) Nonpeptide neuromedin B receptor antagonists inhibit the proliferation of C6 cells. *Eur J Pharmacol* 409:133-142.

Yamada *et al.* (2002) Restraint stress impaired maternal behavior in female mice lacking the neuromedin B receptor (NMB-R) gene. *Neurosci Lett* 330:163-166.

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.