

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE  
HUMAN RECOMBINANT BB2 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1215

**Lot Number:** DC1215-111623

**Quantity:** 1 vial (4 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** Cellbanker 2  
(Amsbio)

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human GRPR cDNA (GenBank accession number NM\_005314.2) with FLAG tag sequence at N-terminus

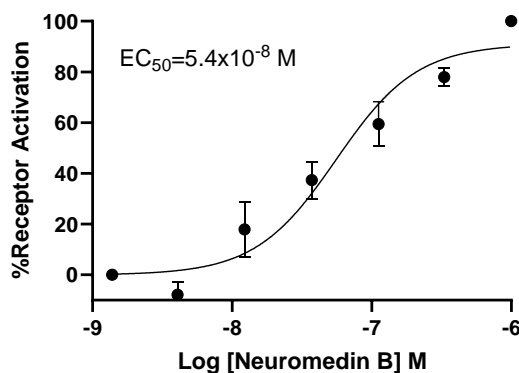
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS

**Background:** The bombesin receptor BB2 (or gastrin-releasing peptide receptor GRPR) is responsible for many physiological actions such as inhibition of feeding, smooth muscle contraction, exocrine and endocrine secretions, thermoregulation, blood pressure and sucrose regulations, and cell growth. BB2 is expressed in the brain, as well as in colon, lung, and prostate cancer cells. The development of potent receptor antagonists that block BB2 receptor responses has potential for new therapeutic treatments in cancer.

**Application:** Functional assays

**Figure 1**



**Figure 1.** Dose-dependent accumulation of intracellular IP1 upon treatment with ligand, measured with IP-one Tb kit.

**References:**

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

Ohki-Hamazaki *et al.* (2005) Development and function of bombesin-like peptides and their receptors. *Int J Dev Biol* 49:293-300.

Xiao *et al.* (2001) The human gastrin-releasing peptide receptor gene structure, its tissue expression and promoter. *Gene* 264:95-103.

**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.