

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE
HUMAN RECOMBINANT BB3 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DH1214

Lot Number: DH1214-122614

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 BRS3 cDNA (GenBank Accession Number NM_001727) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Stability: 1-2 days after thawing

Background: The bombesin-like peptides mediate a diverse spectrum of biological activities and have been implicated as autocrine growth factors in the pathogenesis and progression of cancer. The bombesin receptor subtype 3 (BB3 or BRS3) is expressed in the lung (normal and cancer), nasal mucosa, placenta, and uterus. Mice lacking BB3 receptor develop metabolic defects and obesity phenotype, suggesting that BB3 may be an important target for obesity research. In addition, BB3 may be involved in diabetes and hypertension.

Application: Functional assays

Figure 1

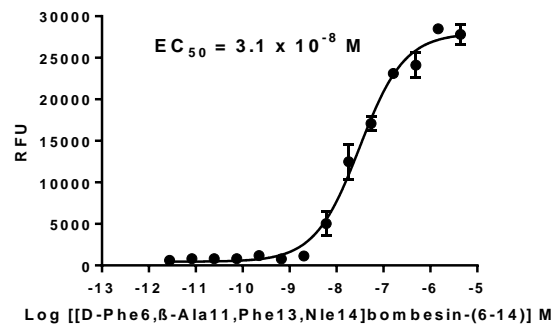


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Fathi *et al.* (1993) BRS-3: a novel bombesin receptor subtype selectively expressed in testis and lung carcinoma cells. *J Biol Chem* 268:5979-5984.

Mantey *et al.* (1997) Discovery of a high affinity radioligand for the human orphan receptor, bombesin receptor subtype 3, which demonstrates that it has a unique pharmacology compared with other mammalian bombesin receptors. *J Biol Chem* 272:26062-26071.

Maekawa *et al.* (2004) Leptin resistance and enhancement of feeding facilitation by melanin-concentrating hormone in mice lacking bombesin receptor subtype-3. *Diabetes* 53:570-576.

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.