

**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT KiSS1 RECEPTOR (GPR54)**

**PRODUCT INFORMATION**

**Catalog Number:** C1036

**Lot Number:** C1036-011411

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

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** CHO dhfr<sup>-</sup>

**Transfection:** Full-length human GPR54 cDNA (GenBank Accession Number NM\_032551)

**Recommended Storage:** Liquid nitrogen upon receiving

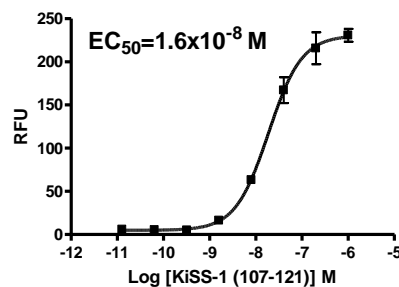
**Propagation Medium:** Alpha-MEM, 10% FBS, 5  $\mu$ g/mL blasticidine

**Stability:** Stable in culture for minimum of two months

**Data sheet**

**Background:** The Kisspeptin receptor KiSS1 is also known as metastin receptor or GPR54. Kisspeptin is a metastasis suppressor protein that suppresses metastasis in malignant melanomas and in some breast carcinomas without affecting tumorigenicity. The metastasis suppressor properties may be mediated in part by cell cycle arrest and induction of apoptosis in malignant cells. The KiSS1 receptor is involved in thyroid cancer, esophageal squamous cell carcinoma and hepatocellular carcinoma.

**Application:** Functional assays



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

**References:**

Shahab *et al.* (2005) Increased hypothalamic GPR54 signaling: a potential mechanism for initiation of puberty in primates. *Proc Natl Acad Sci USA* 102:2129-2134.

Becker *et al.* (2005) Activation of GPR54 promotes cell cycle arrest and apoptosis of human tumor cells through a specific transcriptional program not shared by other Gq-coupled receptors. *Biochem Biophys Res Commun* 326:677-686.

Ikeguchi *et al.* (2004) Clinical significance of the loss of KiSS-1 and orphan G-protein-coupled receptor (hOT7T175) gene expression in esophageal squamous cell carcinoma. *Clin Cancer Res* 10:1379-1383.

**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.

www.multispaninc.com  
sales@multispaninc.com  
support@multispaninc.com

Ver. June 2005

Phone: +1 (510) 887-0817  
Fax: +1 (510) 887-0863  
26219 Eden Landing Road  
Hayward, CA 94545-3718  
U.S.A.