

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

PRODUCT INFORMATION

Catalog Number: DC1283

Lot Number: DC1283-022316

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 GnRHR cDNA (GenBank Accession Number NM_000406) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM + 10% FBS

Stability: 1 – 2 days after thawing

Data sheet

Background: Gonadotropin-releasing hormone receptor (GnRHR) is a receptor for gonadotropin-releasing hormone (GnRH), also known as luteinizing hormone-releasing hormone (LHRH) and prolactin release-inhibiting factor (PIF). GnRHR is expressed on the surface of pituitary gonadotrope cells as well as lymphocytes, breast, ovary, and prostate. Gametes and preimplantation embryos also express GnRH and GnRHR. Binding of GnRH to GnRHR causes the release of follicle stimulating hormone (FSH) and luteinizing hormone (LH) from gonadotropic cells of the anterior pituitary. Mutations in GnRHR may cause hypogonadotropic hypogonadism and may be associated with anosmia.

Application: Functional assays

Figure 1

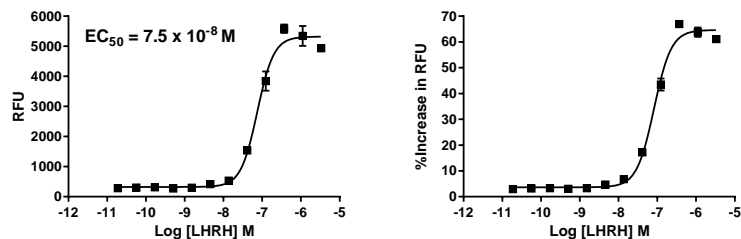


Figure legend: Figure 1. Dose-dependent calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Adelman *et al.* Two mammalian genes transcribed from opposite strands of the same DNA locus. *Science* 235:1514-1517, 1987.

Adelman *et al.* Isolation of the gene and hypothalamic cDNA for the common precursor of gonadotropin-releasing hormone and prolactin release-inhibiting factor in human and rat. *Proc Nat Acad Sci* 83:179-183, 1986.

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.