

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

**PRODUCT INFORMATION**

**Catalog Number:** DC1283

**Lot Number:** DC1283-022316

**Quantity:** 1 vial ( $4 \times 10^6$ ) 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

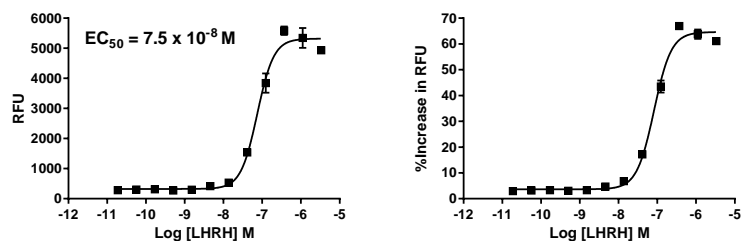
**Stability:** 1 – 2 days after thawing

**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.