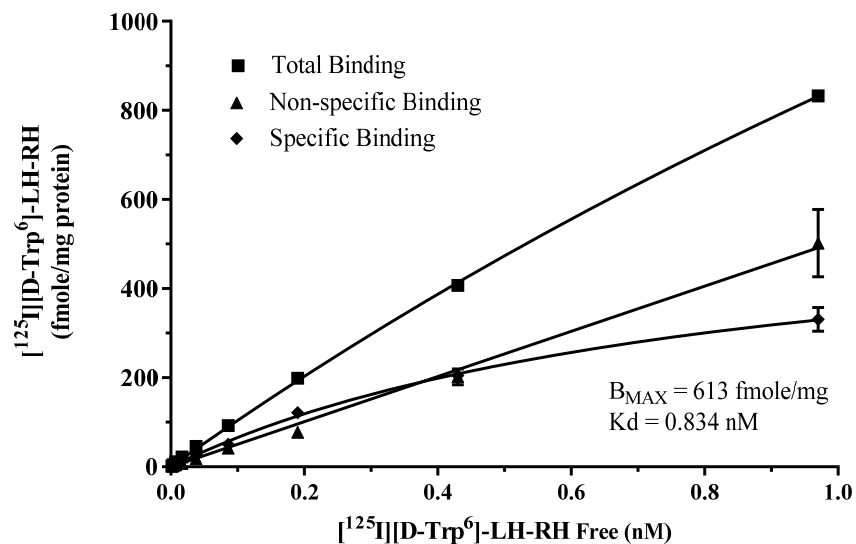


**PRODUCT DATASHEET**
**ChemiScreen™ Gonadotropin-Releasing Hormone Membrane Preparation**

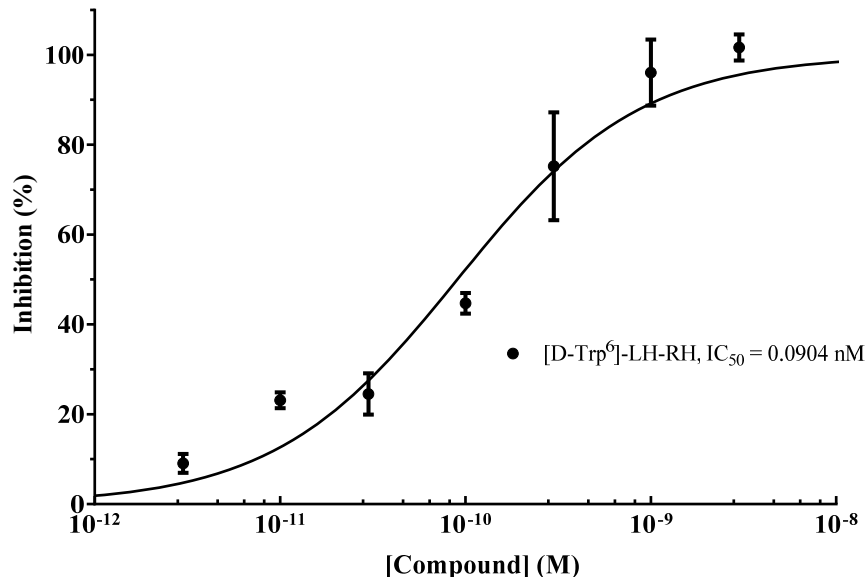
**CATALOG NUMBER:** HTS027M      **QUANTITY:** 200 units  
**LOT NUMBER:** 22C2403      **VOLUME/CONCENTRATION:** 1 mL, 2 mg/mL

**BACKGROUND:** Gonadotropin-releasing hormone (GnRH), also known as luteinizing hormone-releasing hormone (LHRH), regulates the reproductive hormonal cascade in vertebrates. Upon release from the hypothalamus, GnRH stimulates secretion of luteinizing hormone and follicle-stimulating hormone from the pituitary. In humans, the type I GnRH receptor is a GPCR that is unusual in the lack of a cytoplasmic C-terminal tail. Although some species contain a type II GnRH receptor, the human version of this second receptor appears to contain mutations that render it inactive. GnRH analogs (agonists) are used in low doses in a pulsatile fashion in the treatment of infertility, delayed puberty and cryptorchidism. GnRH agonists at high doses desensitize the receptor, and are used along with antagonists in the treatment of hormone-dependent diseases (Millar *et al.*, 2004). GnRH membrane preparations are crude membrane preparations made from our proprietary stable recombinant cell lines to ensure high-level of GPCR surface expression; thus, they are ideal HTS tools for screening of agonists and antagonists at GnRH. The membrane preparations exhibit a  $K_d$  of 0.84 nM for  $[^{125}\text{I}]\text{-[D-Trp}^6\text{]-LHRH}$ . With 0.5 nM  $[^{125}\text{I}]\text{-[D-Trp}^6\text{]-LHRH}$ , 10  $\mu\text{g/well}$  of GnRH Receptor Membrane Prep yields greater than a 3-fold signal-to-background ratio.

**APPLICATIONS:** Radioligand Binding Assay



**Figure 1. Saturation Binding for GnRH.** 10  $\mu\text{g/well}$  of GnRH Receptor Membrane Preparation was incubated with increasing amounts of  $[^{125}\text{I}]\text{-LHRH}$  in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled LHRH. Specific binding (SB) was determined by subtracting NSB from TB. The data are from a representative sample lot.



**Figure 2. Competition Binding for GnRH.** 10  $\mu\text{g}/\text{well}$  GnRH Membrane Preparation was incubated with 0.5 nM [<sup>125</sup>I]-LHRH and increasing concentrations of unlabeled LHRH, and more than a 3-fold signal:background ratio was obtained. The data are from a representative sample lot.

**SPECIFICATIONS:** 1 unit = 10  $\mu\text{g}$   
 $B_{\text{max}}$  for [<sup>125</sup>I]-LHRH binding: 0.61 pmol/mg  
 $K_d$  for [<sup>125</sup>I]-LHRH binding: 0.83 nM  
 Signal:background: >3-fold

**Species:** Human Type I GnRH receptor (Accession number L03380)

**HOST CELLS:** Chem-1, an adherent mammalian cell line without any endogenous GnRH receptor expression.

**RECOMMENDED ASSAY CONDITIONS:** Membranes are mixed with radioactive ligand and unlabeled competitor (see Figures 1 and 2 for concentrations tested) in binding buffer in a nonbinding 96-well plate, and incubated for 2 h at room temperature. Prior to filtration, a GF/C 96-well filter plate is coated with 0.33% polyethyleneimine for 30 min, then washed with 50 mM HEPES, pH 7.4, 0.5% BSA. The binding reactions are transferred to the filter plate, and washed 3 times (1 mL per well per wash) with Wash Buffer. The wells are then dried and counted.

**Binding Buffer:** 50 mM HEPES, pH 7.4, 5 mM  $\text{MgCl}_2$ , 1 mM  $\text{CaCl}_2$ , 0.2% BSA, filtered and stored at 4°C.

**Radioligand:** [<sup>125</sup>I]-[D-Trp<sup>6</sup>]-LHRH (PerkinElmer # NEX365)

**Wash Buffer:** 50 mM HEPES, pH 7.4, 500 mM NaCl, 0.1% BSA, filtered and stored at 4°C.

One package contains enough membranes for at least 200 assays (units), where a unit is the amount of membrane that will yield greater than a 3-fold signal:background ratio with [<sup>125</sup>I]-LHRH at 0.5 nM.

- PRESENTATION:** Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol, and 1% BSA with no preservatives.
- Packaging method: Membrane proteins were adjusted to indicated concentration in 1 mL packaging buffer, rapidly frozen, and stored at -80°C.
- STORAGE/HANDLING:** Store at -80°C. Product is stable for at least 6 months from the date of receipt when stored as directed. Avoid repeated freeze/thaw cycles.
- REFERENCES:**
1. Millar RP *et al.* (2004). Gonadotropin-releasing hormone receptors. *Endocr. Rev.* 25:235-275.

**FOR RESEARCH USE ONLY; NOT FOR USE IN DIAGNOSTIC PROCEDURES. NOT FOR HUMAN OR ANIMAL CONSUMPTION.**

Unless otherwise stated in our catalog or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

No part of these works may be reproduced in any form without permission in writing.

**Eurofins Pharma Bioanalytics Services US Inc.** is an independent member of Eurofins Discovery Services.