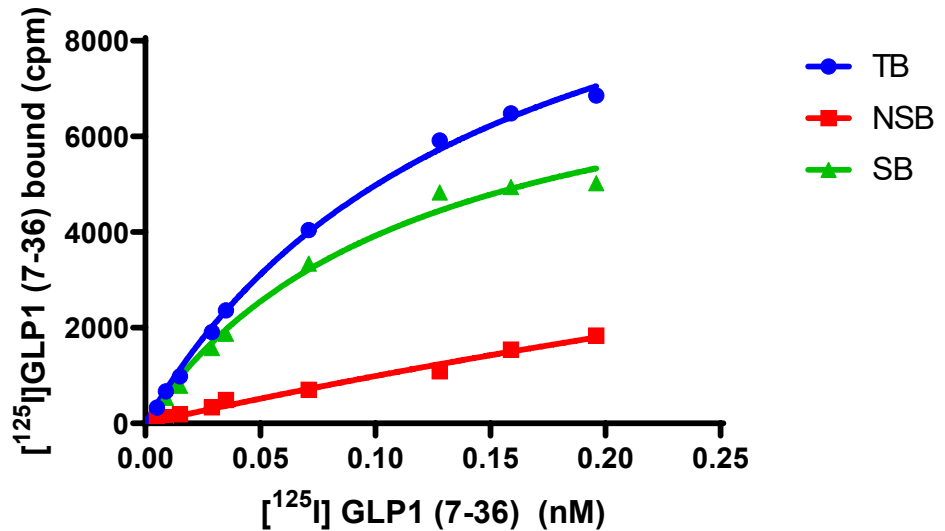


**PRODUCT DATASHEET**
**ChemiScreen™ GLP-1 Glucagon-Like Peptide Membrane Preparation**

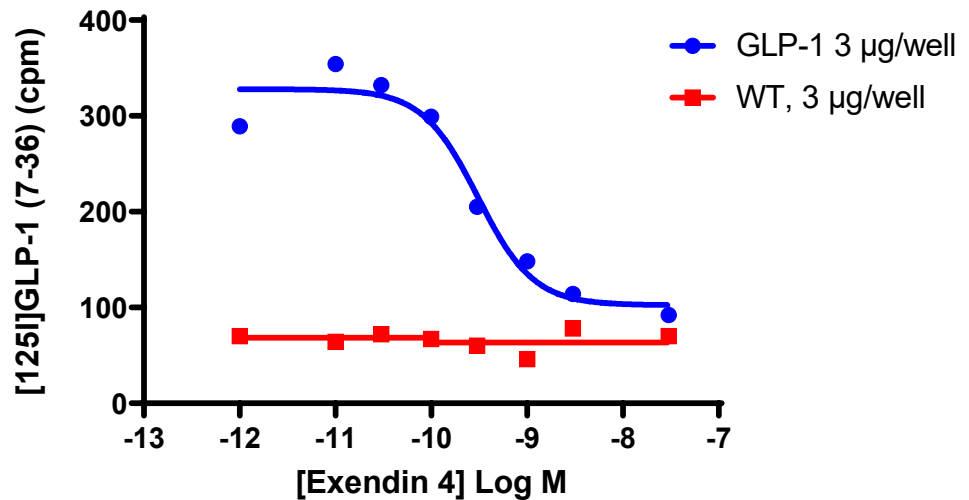
<b>CATALOG NUMBER:</b>	HTS163M	<b>QUANTITY:</b>	200 units
<b>LOT NUMBER:</b>	22M0715	<b>VOLUME/CONCENTRATION</b>	1 mL, 2 mg/mL

**BACKGROUND:** Glucagon-like peptide-I (GLP-1), part of the secretin peptide family, is a class B (class 2) receptor and is involved in glucose-dependent insulin secretion. Insulin secretion is predominantly controlled by glucose levels in the blood, however glucagon stimulates the secretion of insulin (Drucker *et al.* 1987). The GLP-1 receptor couples to G<sub>s</sub> to increase cAMP levels (Mayo *et al.* 2003). GLP-1 has been shown to delay gastric emptying and regulate appetite. Due to the involvement in insulin secretion GLP-1 has been identified as a therapeutic target in type II diabetes (Mayo *et al.* 2003, D'Alessio *et al.* 2004). GLP-1 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 antagonists of GLP-1 interactions with GLP-1.

**APPLICATIONS:** Radioligand binding assay and GTP<sub>γ</sub>S binding



**Figure 1. Saturation binding for GLP-1.** 3 µg/well GLP-1 Membrane Preparation was incubated with increasing amount of [<sup>125</sup>I]-GLP-1(7-36) in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled GLP-1(7-36). Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.



**Figure 2. Competition binding for GLP-1.** GLP-1 Membrane Preparation (3 µg/well) or Wild-Type membrane preparation (WT; Catalog # HTS000MC1) was incubated with 0.025 nM [<sup>125</sup>I]-GLP-1(7-36) and increasing concentrations of unlabeled Exendin-4, and more than 3-fold signal:background was obtained. Representative sample data.

**SPECIFICATIONS:** 1 unit = 10 µg membrane preparation  
 B<sub>max</sub>: 0.83 pmol/mg  
 K<sub>d</sub>: 0.117 nM  
 Signal:background: >3-fold

**TRANSFECTION:** Full-length human GLP-1 cDNA encoding GLP-1 (Accession Number: NM\_002062).

**HOST CELLS:** Chem-9, an adherent cell line expressing the promiscuous G-protein, G<sub>α</sub>15.

**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 1-2 h. Prior to filtration, an FC 96-well harvest plate (Millipore cat. # MAHF C1H) is coated with 0.33% polyethyleneimine for 30 min, then washed with 50mM HEPES, pH 7.4, 0.5% BSA. Binding reaction is transferred to the filter plate, and washed 3 times (1 mL per well per wash) with Wash Buffer. The plate is dried and counted.

**Binding buffer:** 50 mM Hepes, pH 7.4, 5 mM MgCl<sub>2</sub>, 1 mM CaCl<sub>2</sub>, 0.2% BSA, filtered and stored at 4°C

**Radioligand:** [<sup>125</sup>I] GLP-1(7-36) (Perkin Elmer # NEX308)

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

One package contains enough membranes for at least 200 assays (units), where an unit is the amount of membrane that will yield greater than 3-fold signal:background with <sup>125</sup>I-labeled GLP-1 at 0.025 nM.

**PRESENTATION:**

Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol and 1% BSA with no preservatives.

Packaging method: Membranes protein were adjusted to the indicated concentration in packaging buffer, rapidly frozen, and stored at -80°C.

**STORAGE/HANDLING:** Store at  $-70^{\circ}\text{C}$ . Product is stable for at least 6 months from the date of receipt when stored as directed. Do not freeze and thaw.

- REFERENCES:**
1. Drucker *et al.* (1987) Glucagon-like peptide I stimulates insulin gene expression and increases cyclic AMP levels in rat islet cell line. *Proc. Natl. Acad. Sci.*, 84: 3434-3438
  2. Mayo KE *et al.* (2003) International Union of Pharmacology. XXXV. The glucagon receptor family. *Pharmacol. Rev.* 55: 167-194.
  3. D'Alessio D *et al.* (2004) Glucagon-like peptide 1: evolution of an invretin into a treatment for diabetes. *Am. J. Physiol. Endocrinol. Metab.* 286: 882-890.

**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