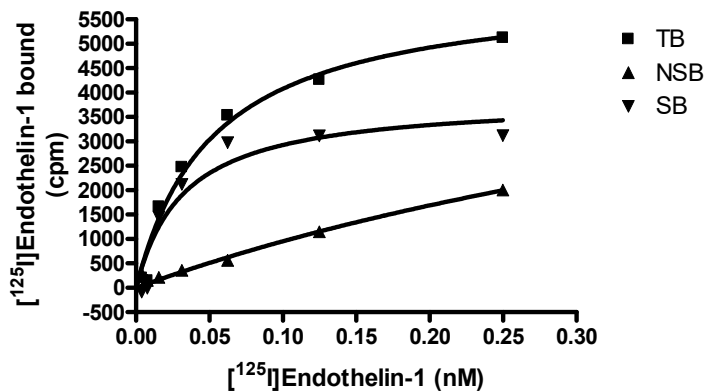


**PRODUCT DATASHEET**
**ChemiScreen™ ET<sub>B</sub> Endothelin Membrane Preparation**

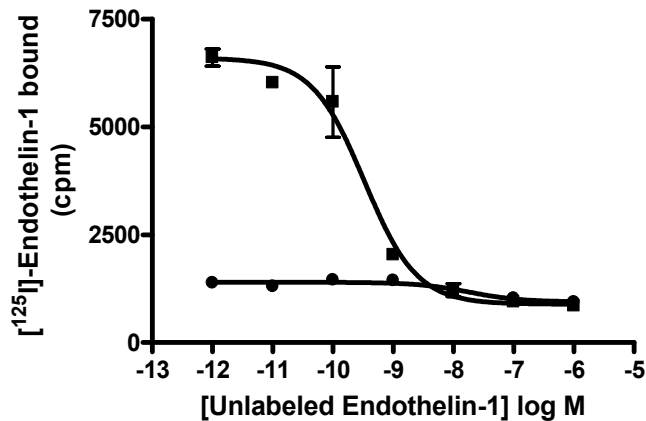
<b>CATALOG NUMBER:</b>	HTS046M	<b>QUANTITY:</b>	200 units
<b>LOT NUMBER:</b>	2071128	<b>VOLUME/CONCENTRATION:</b>	1 mL, 2 mg/mL

**BACKGROUND:** The endothelin family of peptides are potent vasoconstrictors synthesized by endothelial cells in both constitutive and inducible pathways. The biological actions of endothelins are mediated by two GPCRs, ET<sub>A</sub> and ET<sub>B</sub> (Davenport, 2002). Whereas ET<sub>A</sub> is the predominant receptor on smooth muscle cells and thus is the primary mediator of the vasoconstrictor activity, ET<sub>B</sub> is found on endothelial cells and mediates vasodilation (Nilsson *et al.*, 1997). In addition, ET<sub>B</sub> has also been linked to renal failure, congestive heart failure, atherosclerosis, and pulmonary hypertension (D'Orléans-Juste, *et al.*, 2002). A mutation in ET<sub>B</sub> has been found to cause Hirschsprung disease-2, a degenerative disease of the digestive tract (Puffenberger *et al.*, 1994). The cloned human ET<sub>B</sub>-expressing cell line is made in the CHO host, which supports high levels of recombinant ET<sub>B</sub> expression on the cell surface. Thus, the ET<sub>B</sub> Membrane Preparation is an ideal tool for screening for compounds binding to ET<sub>B</sub>. The membrane preparations exhibit a K<sub>d</sub> of 0.03-0.06 nM for [<sup>125</sup>I]-endothelin-1. With 0.2 nM [<sup>125</sup>I]-endothelin-1, 10 μg/well ET<sub>B</sub> Membrane Prep yields greater than 5 fold signal-to-background ratio.

**APPLICATIONS:** Radioligand binding assay



**Figure 1. Saturation Binding for ET<sub>B</sub>.** 10 μg/well ET<sub>B</sub> Membrane Preparation was incubated with increasing amount of [<sup>125</sup>I]Endothelin-1 in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled Endothelin-1. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.



**Figure 2. Competition binding for ET<sub>B</sub>.** ET<sub>B</sub> Membrane Preparation (10 µg/well) or Wild-Type Chem-1 membrane preparation (WT; Catalog # HTS000MC1) was incubated with 0.2 nM [<sup>125</sup>I]-Endothelin-1 and increasing concentrations of unlabeled Endothelin-1, and more than 5- fold signal:background was obtained. Representative sample data.

**Table 1.** Signal:background and specific binding values obtained in a competition binding assay with varying amounts of ET<sub>A</sub> membrane prep.

**SPECIFICATIONS:** 1 unit = 10 µg  
 B<sub>max</sub>: 0.49 pmol/mg  
 K<sub>d</sub>: 0.03-0.06 nM

**Species:** Human ET<sub>B</sub> isoform 1 (Accession Number: NM\_00015)

**HOST CELLS:** CHO-K1 (Chinese hamster ovary cells)

**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 (EMD 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]Endothelin-1 (Perkin Elmer # NEX259)

**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 5-fold signal:background with <sup>125</sup>I-labeled Endothelin-1 at 0.25 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°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. Davenport AP (2002) International Union of Pharmacology. XXIX. Update on Endothelin Receptor Nomenclature. *Pharmacol. Rev.* 54: 219-226.
2. Nilsson T, *et al.* (1997) Presence of contractile endothelin-A and dilatory endothelin-B receptors in human cerebral arteries. *Neurosurgery.* 40: 346-351;comment 351-353
3. D'Orléans-Juste P, *et al* (2002) Function of the endothelin<sub>B</sub> receptor in cardiovascular physiology and pathophysiology. *Pharmacol Ther.* 95: 221-238.
4. Puffenberger EG, *et al* (1994) A missense mutation of the endothelin-B receptor gene in multigenic Hirschsprung's disease. *Cell* 79: 1257-1266.

**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