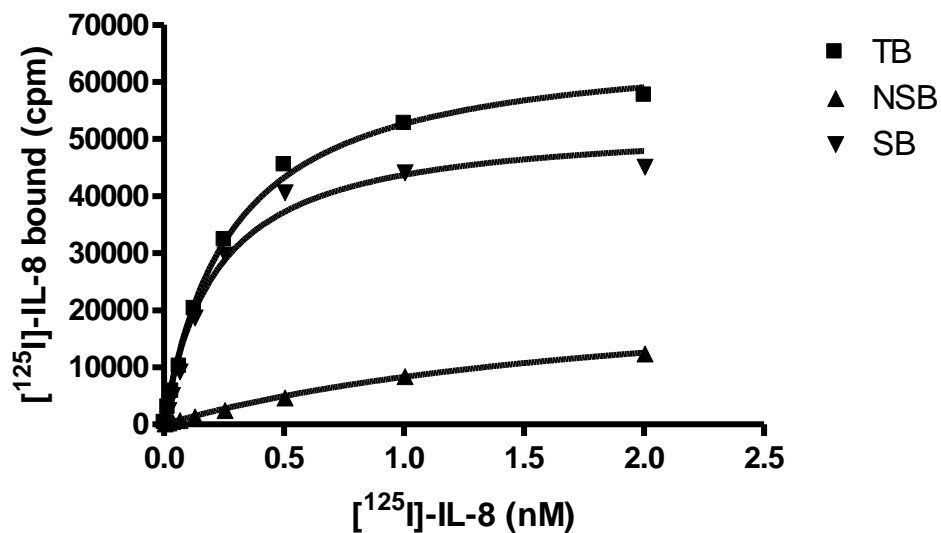


**PRODUCT DATASHEET**
**ChemiScreen™ CXCR1 Chemokine Membrane Preparation**

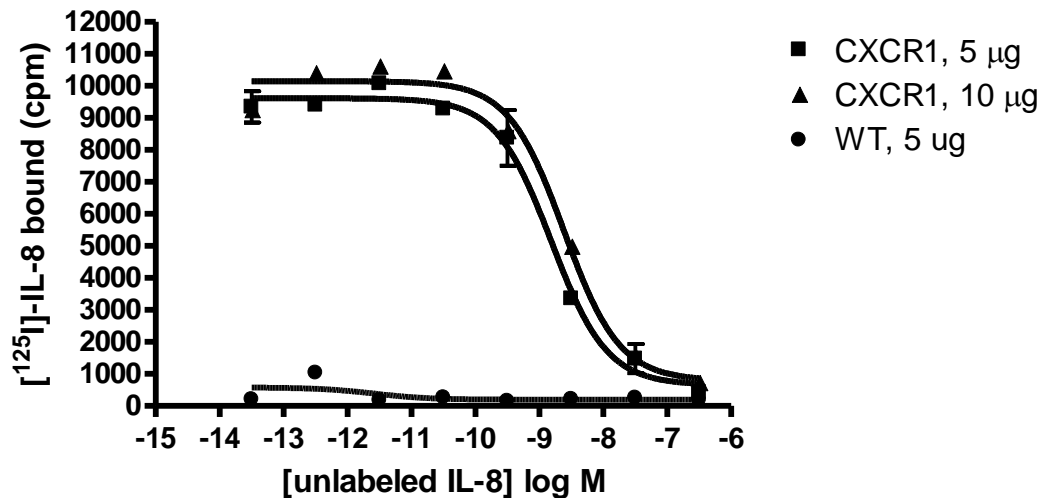
**CATALOG NUMBER:** HTS001M      **QUANTITY:** 200 units  
**LOT NUMBER:**      **VOLUME/CONCENTRATION:** 1 mL, 1 mg/mL

**BACKGROUND:** CXCR1 (also known as IL-8RA) is a Gi/o-coupled GPCR expressed on neutrophils, CD8(+) T cells, and intestinal epithelial cells (Murphy, 1997; Sturm *et al.*, 2005; Takata *et al.*, 2004). CXCR1 binds specifically to the chemokine interleukin-8 (IL-8), whereas a related receptor, CXCR2, binds promiscuously to IL-8 and several other chemokines. CXCR1 has been proposed to be the primary regulator of neutrophil response to IL-8 in sepsis (Cummings *et al.*, 1999). Antagonists of CXCR1 and CXCR2 have been developed and display efficacy in animal models of reperfusion injury (Bertini, *et al.*, 2004; Kaneider *et al.*, 2005). CXCR1 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 CXCR1 interaction with IL-8. The membrane preparations exhibit a Kd of 0.2 nM for [<sup>125</sup>I]-IL-8. With 5 µg/well CXCR1 Membrane Prep and 0.1 nM [<sup>125</sup>I]-IL-8, a greater than 5-fold signal-to-background ratio is obtained.

**APPLICATIONS:** Radioligand binding assay



**Figure 1. Saturation binding for CXCR1.** 10 µg/well CXCR1 Membrane Preparation was incubated with increasing amount of <sup>125</sup>I-labeled Interleukin-8 in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 200-fold excess unlabeled Interleukin-8. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.



**Figure 2. Competition binding for CXCR1.** CXCR1 Membrane Preparation at 5 and 10 µg/well and Wild-Type-Chem-1 Membrane Preparation (WT, cat. # HTS000MC1) were incubated with 0.1 nM <sup>125</sup>I-labeled Interleukin-8 and increasing concentrations of unlabeled Interleukin-8. With CXCR1 Membrane Preparation and unlabeled IL-8, more than 5-fold signal:background was obtained.

**Table 1.** Signal:background and specific binding values obtained in a competition binding assay with CXCR1 membrane preparation in a dose response curve with unlabeled IL-8.

	5 µg/well	10 µg/well
Signal:background	14.8	12.8
Specific binding (cpm)	8965	9624

**SPECIFICATIONS:** 1 unit = 5 µg  
 B<sub>max</sub>: 3.7 pmol/mg  
 K<sub>d</sub>: 0.2 nM

**Species:** Human CXCR1 Receptor (Accession number M68932)

**HOST CELLS:** Chem-1, an adherent mammalian cell line without any endogenous CXCR1 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 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] Interleukin-8 (Perkin Elmer #NEX277)

**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 a unit is the amount of membrane that will yield greater than 5-fold signal:background with <sup>125</sup>I-labeled IL-8 at 0.1 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 the indicated concentration in 1 ml 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. Bertini R *et al.* (2004) Noncompetitive allosteric inhibitors of the inflammatory chemokine receptors CXCR1 and CXCR2: prevention of reperfusion injury. *Proc. Natl. Acad. Sci. USA* 101(32):11791-6.
  2. Cummings CJ *et al.* (1999) Expression and function of the chemokine receptors CXCR1 and CXCR2 in sepsis. *J. Immunol.* 162(4):2341-6.
  3. Kaneider NC *et al.* (2005) Reversing systemic inflammatory response syndrome with chemokine receptor peptidicins. *Nat Med.* 11(6):661-5.
  4. Murphy PM (1997) Neutrophil receptors for interleukin-8 and related CXC chemokines. *Semin. Hematol.* 34: 311-8.
  5. Sturm A *et al.* (2005) CXCL8 modulates human intestinal epithelial cells through a CXCR1 dependent pathway. *Cytokine* 29(1):42-8.
  6. Takata H *et al.* (2004) Cutting edge: expression of chemokine receptor CXCR1 on human effector CD8+ T cells. *J. Immunol.* 173(4):2231-5.

**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