

PRODUCT DATASHEET
ChemiScreen™ CX₃CR1 Chemokine Membrane Preparation

CATALOG NUMBER:	HTS015M	QUANTITY:	200 units
LOT NUMBER:	SC260679	VOLUME/CONCENTRATION:	1 mL, 0.5 mg/mL

BACKGROUND: CX₃CR1 is a GPCR expressed on natural killer cells, cytotoxic T lymphocytes, and macrophages. The sole ligand for CX₃CR1, fractalkine, is an unusual chemokine that is expressed as a transmembrane molecule with a CX3C domain and a mucin domain (Imai *et al.*, 1997). Fractalkine is highly expressed on endothelial cells activated by TNF α and other proinflammatory cytokines, and fractalkine/CX₃CR1 interactions mediate recruitment of macrophages into the atherosclerotic plaque (Lesnick *et al.*, 2003; McDermott *et al.*, 2003). In addition, fractalkine and CX₃CR1 have been implicated in the pathogenesis of glomerulonephritis, HIV infection, and rheumatoid arthritis (Ito *et al.*, 2002; Faure *et al.*, 2003; Nanki *et al.*, 2002). CX₃CR1 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 CX₃CR1 interactions with fractalkine. The membrane preparations exhibit a K_d of 0.21 nM for [¹²⁵I]-Fractalkine. With 2.5 μ g/well of CX₃CR1 Membrane Prep and 0.05 nM [¹²⁵I]-Fractalkine, a greater than 5-fold signal-to-background ratio was obtained.

APPLICATIONS: Radioligand Binding Assay

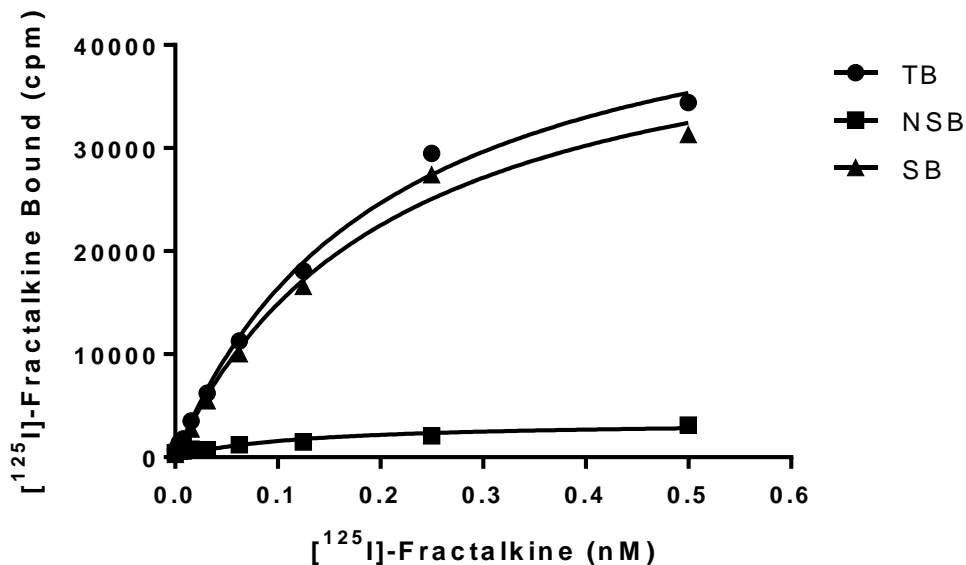


Figure 1. Saturation Binding for CX₃CR1. 2.5 μ g/well of CX₃CR1 Membrane Preparation were incubated with increasing amounts of [¹²⁵I]-Fractalkine in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 200-fold excess unlabeled human recombinant fractalkine. Specific binding (SB) was determined by subtracting NSB from TB. The data are from a representative sample of lot SC260679.

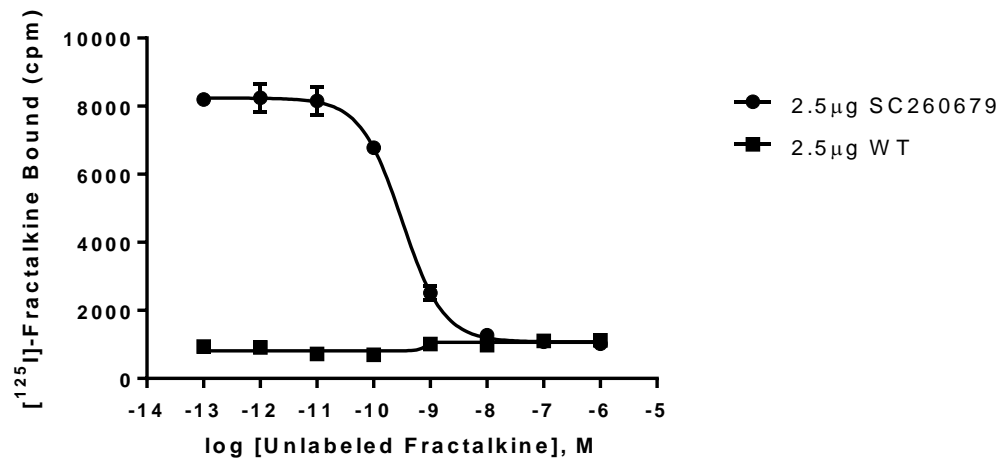


Figure 2. Competition Binding for CX₃CR1. 2.5 µg/well of CX₃CR1 Membrane Preparation or Wild-Type Chem-1 Membrane Preparation (WT; Catalog # HTS000MC1) were incubated with 0.05 nM [¹²⁵I]-Fractalkine and increasing concentrations of unlabeled fractalkine, and more than a 5-fold signal:background ratio was obtained. The data are from a representative sample of lot SC260679.

SPECIFICATIONS: 1 unit = 2.5 µg
 B_{max} for [¹²⁵I]-Fractalkine Binding: 7.5 pmol/mg protein
 K_d for [¹²⁵I]-Fractalkine Binding: 0.21 nM
 Signal:Background: ≥5-fold

SPECIES: Human CX₃CR1 (Accession number U28934)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous CX₃CR1 expression.

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

Binding Buffer: 50 mM HEPES, pH 7.4, 5 mM MgCl₂, 1 mM CaCl₂, 0.2% BSA, filtered and stored at 4°C

Radioligand: [¹²⁵I]-Fractalkine (PerkinElmer # NEX368)

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 5-fold signal:background ratio with [¹²⁵I]-Fractalkine at 0.05 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 0.5 mg/mL in packaging buffer, dispensed at 1 mL per vial, 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. Avoid repeated freeze/thaw cycles.

- REFERENCES:**
1. Faure S *et al.* (2003) Deleterious genetic influence of CX3CR1 genotypes on HIV-1 disease progression. *J Acquir. Immune Defic. Syndr.* 32:335-7.
 2. Imai T *et al.* (1997) Identification and molecular characterization of fractalkine receptor CX3CR1, which mediates both leukocyte migration and adhesion. *Cell* 91:521-30.
 3. Ito Y *et al.* (2002) Fractalkine expression and the recruitment of CX3CR1+ cells in the prolonged mesangial proliferative glomerulonephritis. *Kidney Int.* 61:2044-57.
 4. Lesnick P *et al.* (2003) Decreased atherosclerosis in CX3CR1-/- mice reveals a role for fractalkine in atherogenesis. *J. Clin. Invest.* 111:333-40.
 5. McDermott DH *et al.* (2003) Chemokine receptor mutant CX3CR1-M280 has impaired adhesive function and correlates with protection from cardiovascular disease in humans. *J. Clin. Invest.* 111:1241-50.
 6. Nanki T *et al.* (2002) Migration of CX3CR1-positive T cells producing type 1 cytokines and cytotoxic molecules into the synovium of patients with rheumatoid arthritis. *Arthritis Rheum.* 46:2878-83.

**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.