

PRODUCT DATASHEET
ChemiScreen™ CCR6 Chemokine Membrane Preparation

CATALOG NUMBER: HTS011M **QUANTITY:** 200 units
LOT NUMBER: **VOLUME/CONCENTRATION:** 1 mL, 2 mg/mL

BACKGROUND: CCR6 is a GPCR that binds to a single chemokine, CCL20 (also known as MIP-3 α , LARC or Exodus-1), as well as several members of the beta-defensin family of antimicrobial peptides (Schutyser, 2003). Immature dendritic cells, B-cells, and effector/memory T-cells express CCR6, and the receptor promotes inflammation during the development of rheumatoid arthritis, inflammatory bowel disease, and graft-versus-host disease (Varona et al., 2003; Kleinewietfeld et al., 2005; Varona et al., 2005). CCR6 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 CCR6 interactions with MIP-3 α . The membrane preparations exhibit a K_d of 0.09 nM for [¹²⁵I]-MIP-3 α . With 0.3 nM [¹²⁵I]-MIP-3 α , 5 μ g/well and 10 μ g/well CCR6 Membrane Prep typically yield greater than 6-fold and 8-fold signal-to-background ratios, respectively.

APPLICATIONS: Radioligand binding assay

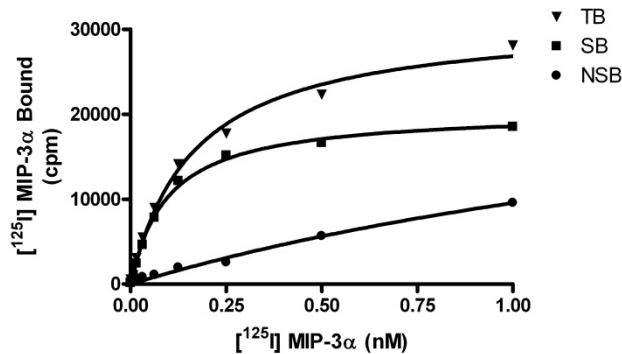


Figure 1. Saturation binding for CCR6. 5 μ g/well CCR6 Membrane Preparation was incubated with increasing amount of [¹²⁵I]-labeled MIP-3 α in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled MIP-3 α . Specific binding (SB) was determined by subtracting NSB from TB.

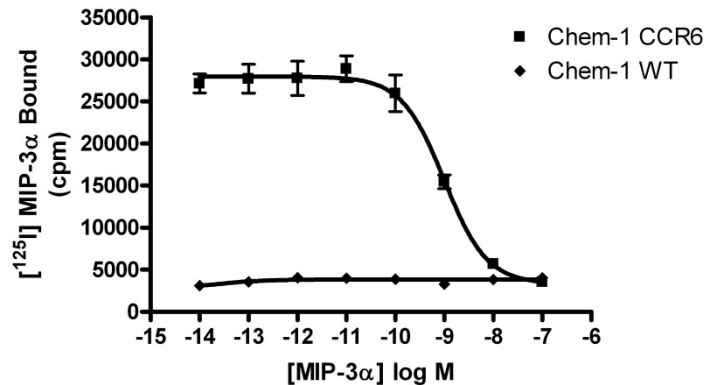


Figure 2. Competition binding for CCR6. 10 µg/well CCR6 Membrane Preparation and Wild-Type Chem-1 Membrane Preparation (catalog #HTS000MC1) were incubated with 0.3 nM ¹²⁵I-labeled MIP-3α and increasing concentrations of unlabeled MIP-3α, and more than 5- fold signal: background was obtained.

SPECIFICATIONS: B_{max}: 2.7 pmol/mg; K_d: 0.09 nM

Species: Human CCR6 (Accession number U45984)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous CCR6 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, a GF/C 96-well filter plate 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 MgCl₂, 1 mM CaCl₂, 0.2% BSA, filtered and stored at 4°C.

Radioligand: [¹²⁵I] MIP-3α. (Perkin Elmer# NEX371)

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 4-fold signal: background with ¹²⁵I-labeled MIP-3α at 0.3 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 2 mg/mL in packaging buffer, and dispensed at 1 mL/vial. Vials were 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. Kleinewietfeld M et al. (2005) CCR6 expression defines regulatory effector/memory-like cells within the CD25(+)CD4+ T-cell subset. *Blood* 105: 2877-86.
2. Schutyser E (2003) The CC chemokine CCL20 and its receptor CCR6. *Cytokine Growth Factor Rev.* 14: 409-26.
3. Varona R et al. (2003) CCR6 has a non-redundant role in the development of inflammatory bowel disease. *Eur. J. Immunol.* 33: 2937-46.
4. Varona R et al. (2005) CCR6 regulates CD4+ T-cell-mediated acute graft-versus-host disease responses. *Blood* 106: 18-26.

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