

PRODUCT DATASHEET
ChemiScreen™ Wild-Type Chem-2 Membrane Preparation

CATALOG NUMBER: HTS000MC2 **QUANTITY:** 1 mL
LOT NUMBER: JH1813910 **VOLUME/CONCENTRATION:** 0.5 mg/mL

BACKGROUND: Chem-2 cells are a suspension cell line endogenously overexpressing G α 16 that is used as a host for the ChemiScreen™ GPCR Membrane Preparations. ChemiScreen™ Wild-Type Chem-2 Membrane Preparations are ideal controls for validation of radioligand binding assays employing ChemiScreen™ GPCR Membrane Preparations in the Chem-2 host cell line.

APPLICATIONS: Negative Control in Radioligand Binding Assay

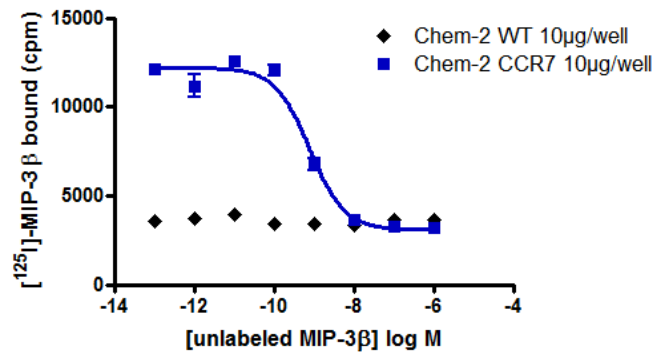


Figure 1. Competition binding of MIP-3 β to CCR7 Membrane Preparations. CCR7 Membrane Preparation (catalog # HTS012M) at 10 μ g/well or Wild-Type Chem-2 Membrane Preparation (catalog # HTS000MC2) at 2.5 μ g/well were incubated with 0.1 nM [125 I]-MIP-3 β and increasing concentrations of unlabeled MIP-3 β . 4-fold signal to background was obtained with the CCR7 Membrane Preparation, whereas negligible signal was obtained with the Wild-Type Chem-2 Membrane Preparation.

HOST CELLS: Chem-2.

PRESENTATION: Membranes were adjusted to 0.5 mg protein/ml in 50 mM Tris pH 7.4, 10% glycerol and 1% BSA, rapidly frozen, and stored at -80oC. Use at the same protein concentration as the ChemiScreen™ GPCR Membrane Preparations of interest.

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.

**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.

**Eurofins Pharma Bioanalytics
 Services US Inc.**

15 Research Park Drive
 St Charles MO 63304
 USA

T | +1 844 522 7787
 F | +1 636 362 7131
www.eurofins.com