

PRODUCT DATASHEET
ChemiScreen™ CCK₂ Cholecystokinin Membrane Preparation

| | | | |
|------------------------|----------|------------------------------|---------------|
| CATALOG NUMBER: | HTS083M | QUANTITY: | 200 units |
| LOT NUMBER: | SC539801 | VOLUME/CONCENTRATION: | 1 mL, 1 mg/mL |

BACKGROUND: Cholecystokinins are a series of peptides of heterogeneous length (5 to 58 amino acids) that are derived from preprocholecystokinin and are found in gastrointestinal tissues and the central nervous system. Gastrin is a related peptide with 5 C-terminal amino acids identical to those of cholecystokinin. Two GPCRs, CCK₁ (CCK_A) and CCK₂ (CCK_B), bind to CCK and/or gastrin to mediate the biological effects of the peptides. CCK₂ binds to CCK and gastrin with similar affinity, whereas CCK₁ selectively binds sulfated CCK. Binding of ligands to CCK₂ stimulates mobilization of intracellular calcium by activation of G_{q/11}. In the periphery, CCK₂ is present in the stomach, where it mediates gastrin-stimulated gastric acid secretion. In the CNS, CCK₂ has been implicated in anxiety, depression, schizophrenia, depression and opioid analgesia (Noble *et al.*, 1999). CCK₂ 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 CCK₂ interactions and its ligands. The membrane preparations exhibit a K_d of 0.47 nM for [¹²⁵I]-cholecystokinin octapeptide (CCK8). With 0.3 nM [¹²⁵I]-CCK8, 5 µg/well of CCK₂ Membrane Prep yields greater than a 20-fold signal-to-background ratio.

APPLICATIONS: Radioligand Binding Assay

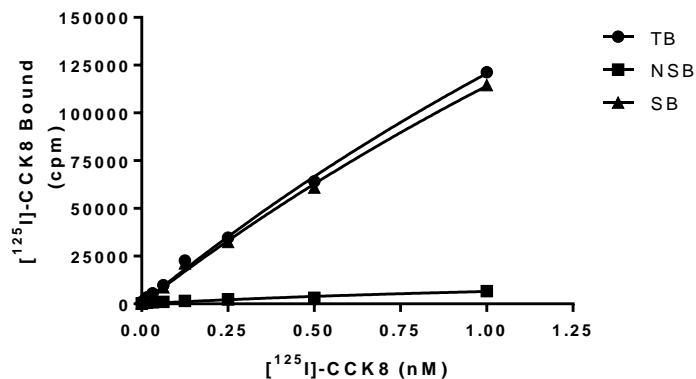


Figure 1. Saturation Binding for CCK₂. 5 µg/well CCK₂ Membrane Preparation was incubated with increasing amounts of [¹²⁵I]-CCK8 in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled gastrin. Specific binding (SB) was determined by subtracting NSB from TB. The data are from a representative sample from lot SC539801.

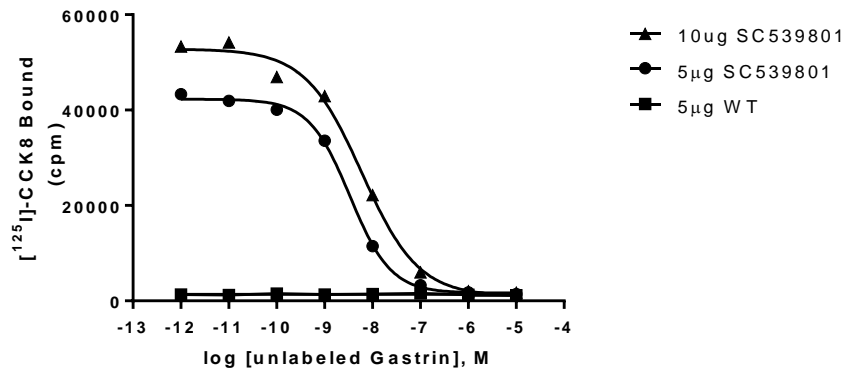


Figure 2. Competition Binding for CCK₂. CCK₂ Membrane Preparation (5 and 10 µg/well in a 96-well plate) were incubated with 0.3 nM [¹²⁵I]-CCK8 and increasing concentrations of unlabeled gastrin, and subjected to filtration binding. The data are from a representative sample from lot SC539801.

SPECIFICATIONS: 1 unit = 5 µg
 B_{max} for [¹²⁵I]-CCK8 binding: 51.5 pmol/mg protein
 K_d for [¹²⁵I]-CCK8 binding: 4.5 nM
 Signal:background: >20-fold

TRANSFECTION: Full-length human CCKB cDNA encoding CCK₂ (Accession Number: NM_176875).

HOST CELLS: Chem-1, an adherent mammalian cell line with no endogenous CCK₂ 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 2 h at room temperature. Prior to filtration, a GF/C 96-well filter plate is coated with 0.33% polyethyleneimine for 30 min, then washed with 50 mM HEPES, pH 7.4, 0.5% BSA. The binding reactions are transferred to the filter plate, and washed 3 times (1 mL per well per wash) with Wash Buffer. The wells are then 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]-CCK8 (PerkinElmer # NEX203).

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 20-fold signal:background ratio with [¹²⁵I]-CCK8 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: Membrane proteins were adjusted to 1.0 mg/mL 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. Avoid repeated freeze/thaw cycles.

REFERENCES:

1. Noble F *et al.* (1999). International Union of Pharmacology. XXI. Structure, distribution, and functions of cholecystinin receptors. *Pharmacol. Rev.* 51:745-781.

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.