

PRODUCT DATASHEET
ChemiScreen™ PK₂ Prokineticins Membrane Preparation

CATALOG NUMBER:	HTS182M	QUANTITY:	200 units
LOT NUMBER:	JH1917577	VOLUME/CONCENTRATION:	1 mL, 2 mg/mL

BACKGROUND: Prokineticins, also known as endocrine gland vascular endothelial growth factors (EG-VEGF), are two ~10 kD secreted proteins originally described to mediate angiogenesis and gastrointestinal smooth muscle contraction (Li *et al.*, 2001; LeCouter *et al.*, 2003). Subsequently, prokineticins have been found to mediate central nervous system functions including circadian rhythms and olfactory bulb development (Cheng *et al.*, 2002; Ng *et al.*, 2005). Two G_q-coupled receptors, PK₁ and PK₂ (also known as GPR73a and GPR73b), mediate cellular responses to prokineticins (Lin *et al.*, 2002). PK₂ 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 agonists and antagonists at PK₂. The membrane preparations exhibit a K_d of 0.36 nM for [¹²⁵I]-MIT-1. With 0.3 nM [¹²⁵I]-MIT-1, 10 μg/well PK₂ Membrane Prep typically yields a 2-fold signal-to-background ratio.

APPLICATIONS: Radioligand binding assay

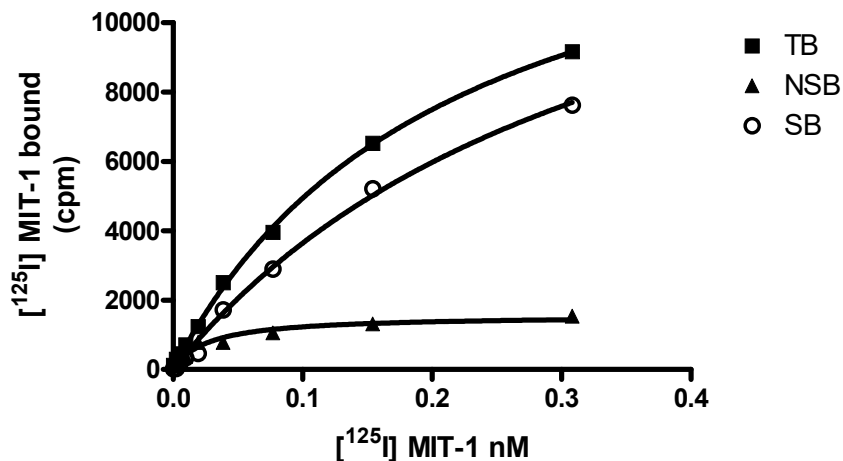


Figure 1. Saturation binding for PK₂. 10 μg/well PK₂ Membrane Preparation was incubated with increasing amount of [¹²⁵I]-labeled MIT-1 in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled EG-VEGF. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.

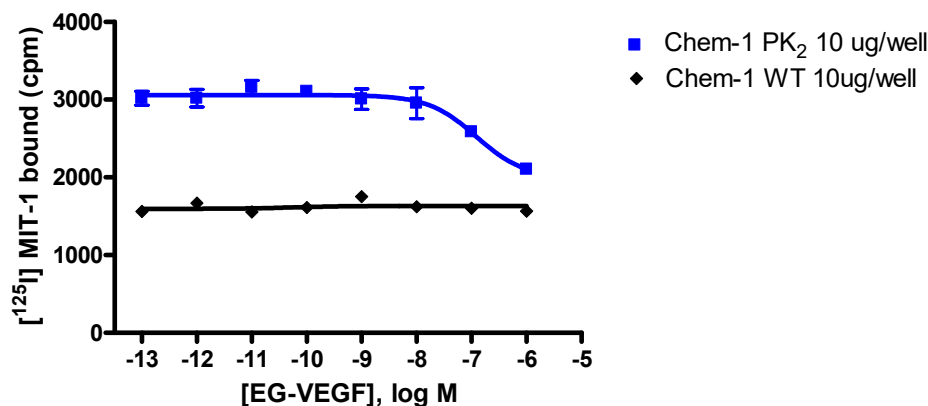


Figure 2. Competition binding for PK₂. 10 μ g/well PK₂ Membrane Preparation and wild-type Chem-1 Membrane Preparation (catalog # HTS000MC1) were incubated in a 96-well plate with 0.3 nM ¹²⁵I-labeled MIT-1 and increasing concentrations of unlabeled EG-VEGF. A 2-fold signal:background was obtained with unlabeled EG-VEGF. Representative sample data.

SPECIFICATIONS: 1 unit = 10 μ g
 B_{max} for [¹²⁵I]- MIT-1 binding: 0.52 pmol/mg protein
 K_d for [¹²⁵I]- MIT-1 binding: ~0.36 nM
Signal:background: >2-fold

TRANSFECTION: Full-length human GPR73L1 cDNA encoding PK₂ (Accession Number: NM_144773)

HOST CELLS: Chem-1, an adherent mammalian cell line with minimum amount of endogenous PK₂ 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 MgCl₂, 1 mM CaCl₂, 0.2% BSA, filtered and stored at 4°C

Radioligand: [¹²⁵I] MIT-1 (Perkin Elmer#: NEX-410)

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 a 2-fold signal:background with ¹²⁵I-labeled MIT-1 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 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. Cheng MY *et al.* (2002) Prokineticin 2 transmits the behavioural circadian rhythm of the suprachiasmatic nucleus. *Nature* 417: 405-10.
2. LeCouter J *et al.* (2003) Endocrine gland-derived VEGF and the emerging hypothesis of organ-specific regulation of angiogenesis. *Nat. Med.* 8: 913-7.
3. Li M *et al.* (2001) Identification of two prokineticin cDNAs: recombinant proteins potently contract gastrointestinal smooth muscle. *Mol. Pharmacol.* 59: 692-8.
4. Lin DC *et al.* (2002) Identification and molecular characterization of two closely related G protein-coupled receptors activated by prokineticins/endocrine gland vascular endothelial growth factor. *J. Biol. Chem.* 277: 19276-80.
5. Ng KL *et al.* (2005) Dependence of olfactory bulb neurogenesis on prokineticin 2 signaling. *Science* 308: 1923-7.

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