

PRODUCT DATASHEET
ChemiScreen™ BB1 Bombesin Membrane Preparation

CATALOG NUMBER: HTS123M **QUANTITY:** 200 units
LOT NUMBER: **VOLUME/CONCENTRATION:** 1 mL, 1 mg/mL

BACKGROUND: Bombesin, a bioactive peptide first identified in amphibian skin, is related to two mammalian peptides, *gastrin*-releasing peptide (GRP) and neuromedin B (NMB). A family of 3 GPCRs, including NMB-R (BB₁), GRP-R (BB₂) and BRS-3 (BB₃), mediate the biological effects of the peptides. The receptors differ in their affinities for the peptides; BB₂ binds to GRP with 50-300-fold greater affinity than to NMB, whereas BB₁ binds to NMB with 10-800-fold greater affinity than to GRP (Tokita *et al.*, 2004). Binding of ligand to BB₁ activates G_q to increase intracellular calcium concentrations. The CNS is a major site of NMB and BB₁ expression, and BB₁ appears to be involved in thermoregulation (Ohki-Hamazaki *et al.*, 2005). BB1 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 BB1 interactions with neuromedin B. The membrane preparations exhibit a K_d of 0.35-0.43 nM for [¹²⁵I]-Bombesin. With 5 μg/well BB1 Membrane Prep and 0.3 nM [¹²⁵I]-bombesin, a greater than 40-fold signal-to-background ratio was obtained.

APPLICATIONS: Radioligand binding assay, and GTPγS binding

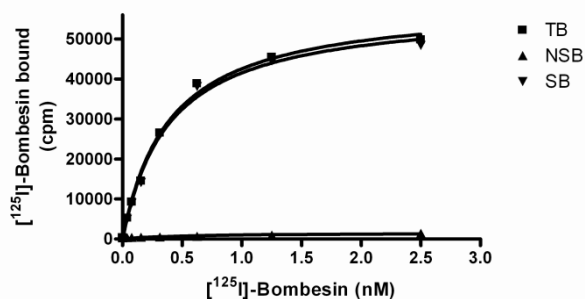


Figure 1. Saturation binding for BB1. 5 μg/well BB1 Membrane Preparation was incubated with increasing amount of [¹²⁵I]-bombesin in the absence (total binding, TB) or presence (nonspecific binding, NSB) of greater than 5000-fold excess unlabeled neuromedin B. Specific binding (SB) was determined by subtracting NSB from TB.

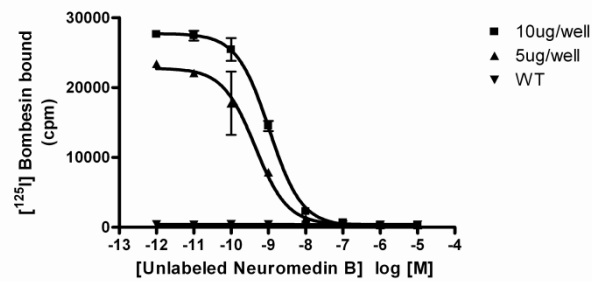


Figure 2. Competition binding for BB1. BB1 Membrane Preparation (5 or 10 μg /well) or Wild-Type Chem-1 membrane preparation (WT; Catalog # HTS000MC1) was incubated with 0.3 nM [^{125}I]-bombesin and increasing concentrations of unlabeled neuromedin B, and more than 40- fold signal:background was obtained.

Table 1. Signal:background and specific binding values obtained in a competition binding assay with varying amounts of BB1 membrane prep.

	10 μg /well	5 μg /well
Signal:background	82.5	74.8
Specific binding (cpm)	27377	22443

SPECIFICATIONS: 1 unit = 5 μg membrane preparation

B_{max} 3.1 pmol/mg

K_d 0.39 nM

Species: Full-length human NMBR cDNA encoding BB₁ (Accession Number: NM_002511)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous BB1 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 MgCl_2 , 1 mM CaCl_2 , 0.2% BSA, filtered and stored at 4°C

Radioligand: [^{125}I] bombesin (Perkin Elmer # NEX258)

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 40-fold signal:background with ^{125}I -labeled bombesin 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. Ohki-Hamazaki H *et al.* (2005) Development and function of bombesin-like peptides and their receptors. *Int. J. Dev. Biol.* 49: 293-300.
 2. Tokita K *et al.* (2004) Molecular basis of the selectivity of gastrin-releasing peptide receptor for gastrin-releasing peptide. *Mol. Pharmacol.* 61: 1435-1443.

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