

PRODUCT DATASHEET

ChemiScreen™ BLT₁ Bombesin Membrane Preparation

CATALOG NUMBER: HTS042M **QUANTITY:** 200 units
LOT NUMBER: 010714 **VOLUME/CONCENTRATION:** 1 mL, 2 mg/mL

BACKGROUND: BLT₁ receptor is a G_i-coupled GPCR that binds primarily to the eicosanoid leukotriene B₄. Macrophages, leukocytes and eosinophils express BLT₁, and in vivo, BLT₁ appears to mediate leukocyte migration and inflammation (reviewed in Brink *et al.*, 2003). EMD BLT₁ 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 BLT₁ interactions with Leukotriene B₄.

APPLICATIONS: Radioligand binding assay

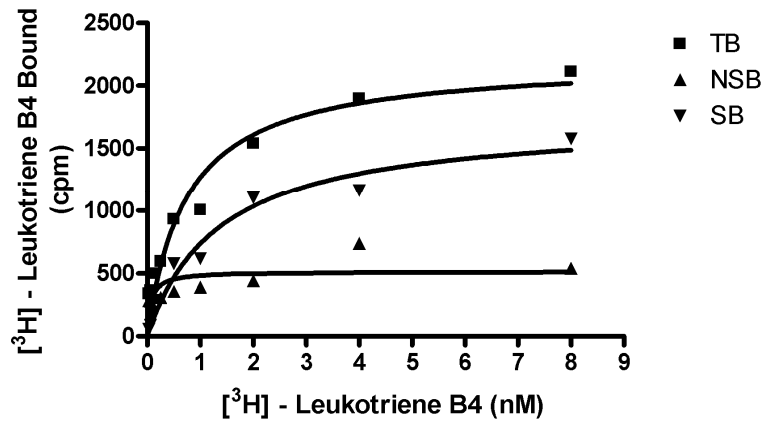


Figure 1. Saturation binding for BLT₁. 5 µg/well BLT₁ Membrane Preparation was incubated with increasing amount of ³H-labeled Leukotriene B₄ in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 200-fold excess unlabeled Leukotriene B₄. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.

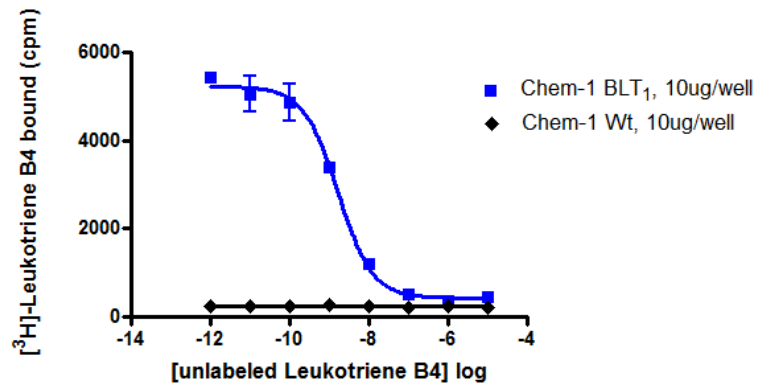


Figure 2. Competition binding for BLT₁. 10 mg/well BLT₁ Membrane Preparation (HTS042M) and Wild-Type Chem-1 Membrane Preparation (catalog # HTS000MC1) were incubated with 1 nM ³H-labeled Leukotriene B₄ and increasing concentrations of unlabeled Leukotriene B₄, and greater than 4-fold signal:background was obtained. Representative sample data.

SPECIFICATIONS: 1 unit = 10 µg membrane preparation
 B_{max} 2.53 pmol/mg
 K_d 1.3 nM
 Signal:background: ≥4-fold

Species: Full-length human BLT₁ (Accession number D89078)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous BLT₁ 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: [³H] Leukotriene (Perkin Elmer# NET852)

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 ³H-labeled Leukotriene B₄ at 1 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. Brink C *et al.* (2003) International Union of Pharmacology XXXVII. Nomenclature for leukotriene and lipoxin receptors. *Pharmacol. Rev.* 55: 195-227.

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