

PRODUCT DATASHEET
ChemiScreen™ CRF1 Neuropeptide Membrane Preparation

CATALOG NUMBER:	HTS023M	QUANTITY:	200 units
LOT NUMBER:	SC20190321	VOLUME/CONCENTRATION:	1 mL, 1 mg/mL

BACKGROUND: The CRF1 receptor is a G_s-coupled GPCR expressed in the brain and pituitary gland that binds to several neuropeptides, including corticotropin-releasing factor (CRF) and urocortin, and the amphibian peptide sauvagine (Chen *et al.*, 1993; Dautzenberg and Hauger, 2002; Bale and Vale, 2004). CRF plays a predominant role in stress response mediated by the hypothalamic-pituitary-adrenal axis, and alterations in CRF and its receptors CRF1 and CRF2 appear to be linked to depression and anxiety (Holsboer, 1999; Bale and Vale, 2004). A number of small molecule antagonists of the CRF1 receptor have been characterized, including R121919, SC241, NBI27914, antalarmin, DMP-696, and CP 154,526. When delivered in animal models of psychiatric disorders, these antagonists display effectiveness in reducing stress-related behaviors (Kehne and De Lombaert, 2002). CRF1 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 CRF1 interactions with its ligands. The membrane preparations exhibit a K_d of 1.76 nM for [¹²⁵I]-sauvagine.

APPLICATIONS: Radioligand binding assay

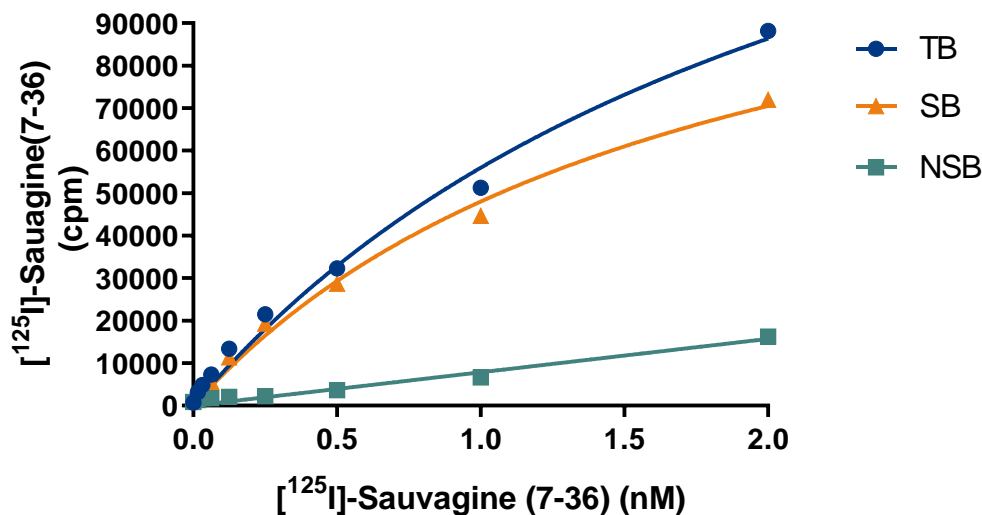


Figure 1. Saturation binding for CRF1. 5 µg/well CRF1 Membrane Preparation was incubated with increasing amount of ¹²⁵I-labeled Sauvagine in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 200-fold excess unlabeled sauvagine. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.

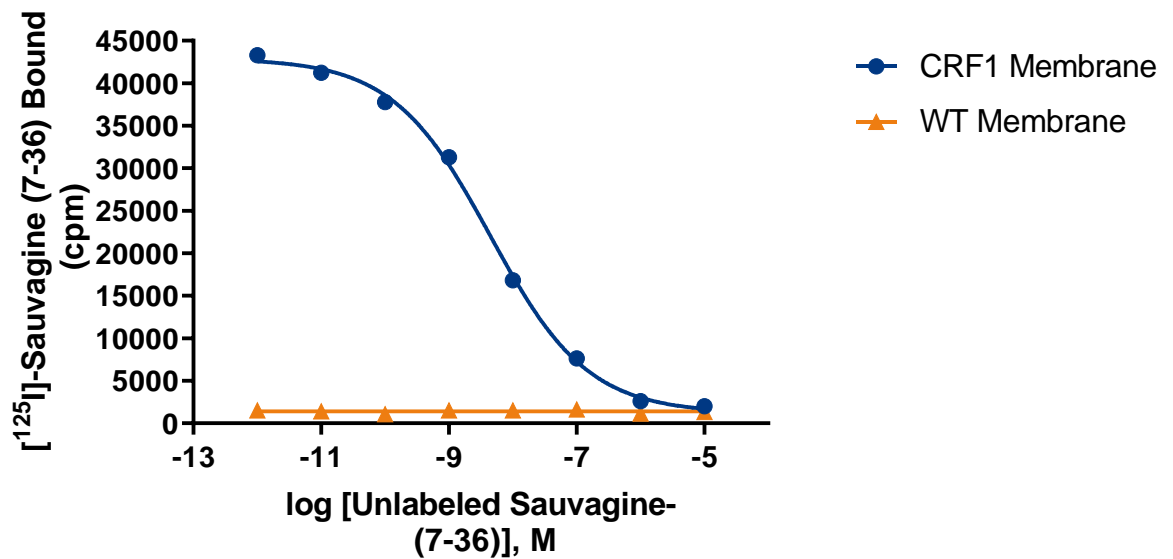


Figure 2. Competition binding for CRF1. 5 µg/well CRF1 Membrane Preparation (HTS023M) was incubated with 0.50 nM ¹²⁵I-labeled Savagine and increasing concentrations of unlabeled savagine, and more than 10- fold signal:background was obtained. Representative sample data.

SPECIFICATIONS: 1 unit = 5 µg
 B_{max} : 10.88 pmol/mg
 K_d : 1.76 nM
 Signal:background: ≥10-fold

Species: Human CRF1 (Accession number X72304)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous CRF1 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 (Costar 3605), 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: [¹²⁵I] savagine (Perkin Elmer#: NEX306)

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 10-fold signal:background with ¹²⁵I-labeled Savagine at 0.50 nM.

PRESENTATION:

Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol and 1% BSA no preservatives. Packaging method: Membranes protein were adjusted to the indicated concentration in 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. Bale TL and Vale WW (2004) CRF and CRF receptors: role in stress responsivity and other behaviors. *Annu. Rev. Pharmacol. Toxicol.* 44: 525-557.
2. Chen R. *et al.* (1993) Expression cloning of a human corticotropin-releasing factor receptor. *Proc. Natl. Acad. Sci. USA* 90: 8967-8971.
3. Dautzenberg FM and Hauger RL (2002) The CRF peptide family and their receptors: yet more partners discovered. *Trends Pharmacol. Sci.* 23: 71-77.
4. Holsboer F (1999) The rationale for corticotropin-releasing hormone receptor (CRH-R) antagonists to treat depression and anxiety. *J. Psychiatr. Res.* 33: 181-214.
5. Kehne J and De Lombaert S (2002) Non-peptidic CRF1 receptor antagonists for the treatment of anxiety, depression and stress disorders. *Curr. Drug Targets CNS Neurol. Disord.* 1: 467-493.

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